

# This product distributed by **Takara Bio USA, Inc.**

Ordering information 1.800.662.2566 ordersUS@takarabio.com Technical support 1.800.662.2566 techUS@takarabio.com

Clontech TakaRa cellartis

takarabio.com

use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Takara Bio Inc. @2018

# DNA isolation from FFPE samples

User manual

NucleoSpin<sup>®</sup> DNA FFPE XS

February 2018/Rev. 03





# **DNA Isolation from FFPE Samples**

# Protocol-at-a-glance (Rev. 03)

	NucleoSpin <sup>®</sup> DNA FFPE XS				
	DNA	isolati	Protocol 5.1: on with Paraffin Dissolver	D	<i>Protocol 5.2:</i> NA isolation with xylene
Sample preparation			For appropriate sample quantity see section 2.4.		For appropriate sample quantity see section 2.4
1 Deparaffinize			400 $\mu L$ Paraffin Dissolver		1 mL xylene
sample			60 °C, 3 min Mix hot sample		RT, 2 min Mix
				Ò	11,000 x <i>g</i> , 2 min Discard supernatant
			Let sample cool down	0	1 mL ~ 98 % ethanol Mix
				Ò	11,000 x g, 2 min Discard supernatant
					Dry at 60 °C, 3–10 min
2 Lyse sample	-		100 μL FL Mix vigorously		100 μL FL
		Ċ	11,000 x g, 1 min		-
	$ $ $\forall$	0	10 μL Proteinase K Mix lower phase		10 μL Proteinase K Mix
			RT, 3 hours or overnight		RT, 3 hours or overnight
3 Decrosslink		Ø	100 μL D-Link Mix gently		100 μL D-Link Mix gently
	V		11,000 x g, 30 s 90 °C, 30 min		90 °C, 30 min
4 Adjust binding conditions		Ċ	200 µL ~ 98 % ethanol Mix		200 μL ~ 98 % ethanol Mix
	V		1,000 x g, 1 s		-
5 Bind RNA	Ĵ		Load aqueous		Load lysate
		$\bigcirc$	2,000 x <i>g</i> , 30 s		2,000 x <i>g</i> , 30 s
6 Wash and dry silica membrane	8	1 <sup>st</sup>	400 μL B5 11 000 x α 30 s		400 μL B5 11.000 x α.30 s
		2 <sup>nd</sup>	400 uL B5		400 ul B5
	9	Ò	11,000 x g, 2 min		11,000 x <i>g</i> , 2 min
7 Elute DNA	-		20 uL BE		20 uL BE
		O	11,000 x <i>g</i> , 30 s		11,000 x <i>g</i> , 30 s
8 Optional: Remove residual ethanol			90 °C, 8 min		90 °C, 8 min

MACHEREY-NAGEL GmbH & Co. KG · Neumann-Neander-Str. 6–8 · 52355 Düren · Germany Tel.: +49 24 21 969-270 · Fax: +49 24 21 969-199 · tech-bio@mn-net.com · *www.mn-net.com* 



# **Table of contents**

1	Com	ponents	4
	1.1	Kit contents	4
	1.2	Reagents, consumables, and equipment to be supplied by user	5
	1.3	About this user manual	5
2	Prod	uct description	6
	2.1	The basic principle	6
	2.2	Kit specifications	7
	2.3	Handling, preparation, and storage of starting materials	8
	2.4	Quantities of FFPE sections	8
	2.5	Elution procedures	9
	2.6	Stability of isolated DNA	9
	2.7	Removal of residual traces of ethanol for highest sensitivity in downstream applications	10
3	Stora	age conditions and preparation of working solutions	12
4	Safe	ty instructions	13
5	Proto	pcols	15
	5.1	DNA purification from FFPE samples using Paraffin Dissolver	15
	5.2	DNA purification from FFPE samples with xylene deparaffinization	20
6	Appe	andix	24
	6.1	Comments on DNA quality and quantity	24
	6.2	Troubleshooting	25
	6.3	Ordering information	27
	6.4	Product use restriction / warranty	28

# 1 Components

## 1.1 Kit contents

	NucleoSpin <sup>®</sup> DNA FFPE XS		
REF	10 preps 740980.10	50 preps 740980.50	250 preps 740980.250
Paraffin Dissolver	5 mL	25 mL	125 mL
Lysis Buffer FL	8 mL	8 mL	8 mL
Decrosslink Buffer D-Link	8 mL	8 mL	30 mL
Wash Buffer B5 (Concentrate)*	6 mL	12 mL	50 mL
Proteinase K (lyophilized)*	6 mg	30 mg	75 mg
Proteinase Buffer PB	1.8 mL	1.8 mL	8 mL
Elution Buffer BE**	13 mL	13 mL	13 mL
NucleoSpin <sup>®</sup> DNA FFPE XS Columns (green rings plus Collection Tubes)	10	50	250
Collection Tubes (2 mL)	20	100	500
User manual	1	1	1

<sup>\*</sup> For preparation of working solutions and storage conditions see section 3.

<sup>\*\*</sup> Composition of Elution Buffer BE: 5 mM Tris/HCI, pH 8.5

# 1.2 Reagents, consumables, and equipment to be supplied by user

#### Reagents

- 96–100 % ethanol (undenaturated ethanol is preferable) to prepare Wash Buffer B5 and to adjust binding conditions.
- Optional for deparaffinisation without Paraffin Dissolver: Xylene, d-Limonene, mixtures of isoparaffinic hydrocarbons, or similar reagents for deparaffinization.

Consumables

- 1.5 mL microcentrifuge tubes (for sample lysis and DNA elution)
- Disposable pipette tips

#### Equipment

- Manual pipettors
- Centrifuge for microcentrifuge tubes
- Vortex mixer
- Thermal heating-block (adjustable to 60 °C and 90 °C)
- Personal protection equipment (e.g., lab coat, gloves, goggles)

### 1.3 About this user manual

It is strongly recommended that first time users of the **NucleoSpin® DNA FFPE XS** kit read the detailed protocol sections of this user manual. Experienced users, however, may refer to the Protocol at a glance instead. The Protocol at a glance is designed to be used only as a supplemental tool for quick referencing while performing the purification procedure.

All technical literature is available on the internet at www.mn-net.com.

Please contact Technical Service regarding information about changes of the current user manual compared to previous revisions.

# 2 Product description

Formalin-fixed, paraffin-embedded (FFPE) tissue samples are routinely prepared from human surgical tissue samples by fixation with formalin and embedding in paraffin. Thin sections of FFPE samples are commonly subjected to histophathological analysis and remaining paraffin-tissue blocks are usually archived. Existing extensive archives of FFPE tissue samples represent a valuable source for retrospective studies of gene expression patterns and mutation analysis. However, the use of such samples for DNA analysis is limited due to chemical modification by formaldehyde and fragmentation of the DNA during tissue processing (sampling, fixing, embedding) and storage (humidity, time, temperature) of the samples. Standard DNA isolation procedures often result in low DNA yield or poor performance in downstream applications (e.g., PCR). A special purification system taking the unique requirements of FFPE tissue into account is inevitably necessary for successful analysis of nucleic acids from FFPE samples.

## 2.1 The basic principle

The **NucleoSpin® DNA FFPE XS** kit provides a convenient, reliable, and fast method to isolate DNA from formalin-fixed, paraffin-embedded (FFPE) tissue specimen. The procedure omits the use of flammable and malodorous xylene or d-limonene commonly used for deparaffinization. Further, the procedure omits the difficult removal of organic solvent from often barely visible tissue pellets. **NucleoSpin® DNA FFPE XS** employs the odorless Paraffin Dissolver (patent pending) and allows efficient lysis in a convenient two-phase system.

First, the paraffin of FFPE sections is dissolved in the Paraffin Dissolver. Tissue is then digested by proteinase to solubilize the fixed tissue and release DNA into solution. Subsequently, heat incubation with specially designed buffer effectively eliminates crosslinks from the previously released DNA. After addition of ethanol, the lysate is applied to the **NucleoSpin® DNA FFPE XS Column**. DNA is bound to the silica membrane. Two washing steps help remove salts, metabolites, and macromolecular cellular components. Pure DNA is finally eluted under low ionic strength conditions in a small volume (20  $\mu$ L) of Elution Buffer BE, yielding highly concentrated DNA.

DNA preparation using **NucleoSpin® DNA FFPE XS** kits can be performed at room temperature. The eluate, however, should be treated with care, because the Elution Buffer BE does not contain DNase inhibitors like EDTA. To ensure DNA stability store frozen DNA at -20 °C.

## 2.2 Kit specifications

- The NucleoSpin<sup>®</sup> DNA FFPE XS kit is recommended for the isolation of DNA from formalin-fixed, paraffin-embedded (FFPE) tissue samples. Samples are typically thin sections (approx. 3–20 µm thickness) of human or animal origin usually obtained by tissue resection or biopsy.
- Sample amount: The maximum sample size is determined by a) the amount of tissue and b) by the amount of paraffin.
   NucleoSpin<sup>®</sup> DNA FFPE XS is suitable for up to 5 mg tissue.
   The amount of paraffin is limited to 15 mg, when using the standard protocol with Paraffin Dissolver (ca. 7 sections of 10 µm x 250 mm<sup>2</sup>). However, larger amounts of paraffin samples may be processed by using either additional Paraffin Dissolver or by deparaffinization using xylene (see also section 2.4).
- DNA yield strongly depends on the sample type, quality, quantity, and time of storage. Further, measured DNA yield may vary considerably between different quantification methods. Yield determined by absorption measurement at 260 nm or by a fluorescent dye (e.g., PicoGreen<sup>®</sup>) may deviate from values obtained by quantification with PCR. Even quantification values obtained via PCR with a short (e.g., 80 bp) and a long (e.g., 300 bp) amplicon may also differ considerably. The deviation of quantification also depends on DNA size distribution as well as on efficiency of decrosslinking (or extent of remaining crosslinks). Please also see section 6.1 for considerations on determining DNA quality and quantity.
- The innovative column design with a funnel shaped thrust ring and a small silica membrane area allows elution of DNA in as little as 5–30 μL. Thus, eluted DNA is highly concentrated and ready to use in all common downstream applications (e.g., PCR).
- **DNA size distribution**: DNA isolated from formalin-fixed, paraffin-embedded tissue shows size distribution from 50 to 5,000 bases. Often short sized DNA from ca. 100–300 bases predominate, especially when the sample material is old. However, samples which were subjected to good tissue fixation, embedding, and storage conditions can yield DNA even larger than 5,000 bases.
- **DNA preparation time** strongly depends on the sample and the required lysis time. For best results lysis is performed at room temperature for at least three hours. For some kinds of sample a longer lysis (e.g., overnight) will even result in remarkably higher DNA yield.

Table 1: Kit specifications at a glance			
Parameter	NucleoSpin <sup>®</sup> DNA FFPE XS		
Sample material*	Up to 7 sections, 10 $\mu m,$ surface of 250 $mm^2$		
Typical yield	Strongly depends on sample quality and amount		
Elution volume	5–30 μL		
Maximum loading volume	600 μL		
Format	Mini spin column – XS design		

## 2.3 Handling, preparation, and storage of starting materials

Many factors influence the yield and usability of DNA obtained from FFPE samples. The procedure of tissue sampling, post sampling delay before fixation, fixation time, embedding, and storage conditions have a high impact on DNA quality and yield.

Starting from a paraffin embedded tissue block, samples should be sectioned under clean conditions. Paraffin sections may be stored at +4 °C or lower for at least several weeks without observable effects on DNA yield or usability. Long term storage of paraffin sections may have a negative effect on the DNA due to air oxidation.

Wear gloves at all times during the preparation. Change gloves frequently.

### 2.4 Quantities of FFPE sections

The standard protocol (section 5.1.) allows the preparation of FFPE samples with approximately 15 mg (ca. 17  $\mu$ L) paraffin. This corresponds to:

 $\sim 17~sections$  of 10  $\mu m$  thickness and 100  $mm^2$  area

 $\sim 7$  sections of 10  $\mu m$  thickness and 250  $mm^2$  area

 $\sim\!5$  sections of 10  $\mu m$  thickness and 325  $mm^2$  area

~4 sections of 10  $\mu m$  thickness and 400  $mm^2$  area

 $\sim\!3$  sections of 10  $\mu m$  thickness and 575  $mm^2$  area

 $\sim\!2$  sections of 10  $\mu m$  thickness and 840  $mm^2$  area

 $\sim$  1 section of 10  $\mu m$  thickness and 1680  $mm^2$  area

Larger amounts of paraffin can be dissolved by adding a higher volume of Paraffin Dissolver (REF 740968.25) to the sample initially (30  $\mu$ L Paraffin Dissolver per mg paraffin), or by using xylene for deparaffinization as described in section 5.2. When using more than 400  $\mu$ L Paraffin Dissolver per preparation, it is necessary to use a collection tube larger than 1.5 mL to enable removal of the lower, aqueous phase after the decrosslink step without spillage.

<u>Note:</u> The NucleoSpin<sup>®</sup> DNA FFPE XS standard procedure is recommended for samples containing **up to 15 mg** paraffin (to ensure efficient dissolving of the paraffin with the

<sup>\*</sup> When using the standard procedure with Paraffin Dissolver. Processing larger quantities is possible with protocol modifications, see section 2.4.

indicated volume of Paraffin Dissolver) and **up to 5 mg tissue** (to avoid an overloading of the membrane).

- Three sections of 20 mm x 25 mm area and 10 μm thickness can contain up to 15 mg paraffin (especially, if only minor parts of the section contain tissue).
- One section of 20 mm x 25 mm area and 10 µm thickness does contain approximately 5 mg tissue, if the section contains area-wide tissue.

### 2.5 Elution procedures

High DNA concentration in the elution fraction is desirable for all typical downstream applications. With regard to limited volumes of reaction mixtures, high template concentration can be a crucial criterion. Due to a large default elution volume, standard kits often result in low concentrated DNA, when small samples are processed. Such DNA samples may even require a subsequent concentration to be suitable for the desired application.

**NucleoSpin<sup>®</sup> DNA FFPE XS** kits allow efficient elution in very small volumes resulting in highly concentrated DNA. Elution volumes in the range of 5–30  $\mu$ L are recommended, the default volume is 20  $\mu$ L.



Figure 1 Correlation between elution volume and DNA concentration

(NucleoSpin<sup>®</sup> DNA FFPE XS Columns)

## 2.6 Stability of isolated DNA

Due to its composition, the Elution Buffer does not inhibit DNases, i.e., it does not contain substances (e.g., EDTA) to complex divalent cations. Therefore, be aware not to contaminate Elution Buffer with DNase!

For short term DNA solution may be stored at 0–4  $^\circ C$  and for long term storage at -20  $^\circ C$  is recommended.

<sup>\*</sup> The maximum percentage of template volume in a PCR reaction may vary depending on the robustness of the PCR system; 40 % template volume were tested using LightCycler<sup>®</sup> PCR (Roche) with the DyNAmo™ Capillary SYBR<sup>®</sup> Green qPCR Kit (Finnzymes).

# 2.7 Removal of residual traces of ethanol for highest sensitivity in downstream applications

The default elution volume of **NucleoSpin® DNA FFPE XS** is 20 µL. The kit allows even lower elution volumes down to 5 µL to increase the DNA concentration (see section 2.5). Be aware that a reduction of the 20 µL default elution volume will also increase the concentration of residual ethanol in the eluate.

For the default elution volumes a heat incubation of the eluate is recommended if the eluate comprises more than 20% of the final PCR volume (incubate eluate with open lid for 8 min at 90 °C). Inhibition of sensitive downstream reactions can be avoided by this precautional measure.

In this context, please mind the remarks below:

a) An incubation of the elution fraction at higher temperatures will increase signal output in PCR. This is especially of importance if the template represents more than 20% of the total PCR reaction volume (e.g., more than 4  $\mu$ L eluate used as template in a PCR reaction with a total volume of 20  $\mu$ L).

The template may represent up to 40  $\%^*$  of the total PCR reaction volume, if the eluate is incubated at 90 °C for 8 min as described above.

- b) Typically 20  $\mu$ L eluate will evaporate to 12–14  $\mu$ L during heat incubation for 8 min at 90 °C. If higher final volumes are required, please increase the volume of elution buffer (e.g., from 20  $\mu$ L to 30  $\mu$ L).
- c) An incubation of the elution fraction for 8 min at 90 °C will denature DNA. If non denatured DNA is required (for downstream applications other than PCR; e.g., ligation or cloning), we recommend incubating at a temperature below 80 °C for a longer time as most DNA has a melting point above 80 °C. Suggestion: incubate for 17 min at 75 °C.
- d) The incubation of the eluate at higher temperatures may be adjusted according to Figure 2. The incubation time- and temperature conditions shown will reduce an elution volume of 20  $\mu$ L to about 12–14  $\mu$ L and will effectively remove traces of ethanol as described above.
- e) If the initial volume of elution buffer applied to the column is less than 20  $\mu$ L, heat incubation time should be reduced in order to avoid complete dryness. If the elution volume is for example 5  $\mu$ L, a heat incubation of the eluate for 2 min at 80 °C will adequately remove residual ethanol.



#### Figure 2 Removal of residual ethanol from the elution fraction by heat treatment.

In order to obtain highest PCR sensitivity, heat incubation of the eluate is recommended. Heat incubation may be performed at temperatures of 70–90 °C in a heat block, with or without shaking. Effective conditions (temperature, time, and shaking rate) for ethanol removal can be read from the diagram; an initial volume of 20  $\mu$ L will evaporate to 12–14  $\mu$ L during the described incubation.

# 3 Storage conditions and preparation of working solutions

#### Attention:

Buffers FL contains chaotropic salts. Wear gloves and goggles!

- All kit components should be stored at room temperature (18–25 °C) and are stable up to one year. Storage at lower temperatures may cause precipitation of salts.
- Check that 96–100 % ethanol is available (undenaturated ethanol is preferable) to adjust the binding conditions in the lysate and to prepare Wash Buffer B5 (see below).

Before starting protocol prepare the following:

- Proteinase K: Add the indicated volume of Proteinase Buffer PB (see following table or on the vial) to dissolve lyophilized Proteinase K. Proteinase K solution is stable at -20 °C for 6 months.
- Wash Buffer B5: Add the indicated volume of 96–100 % ethanol (see following table or on the bottle) to Buffer B5 Concentrate. Store Wash Buffer B5 at room temperature (18–25 °C) for up to one year.

	NucleoSpin <sup>®</sup> DNA FFPE XS			
REF	10 preps 740980.10	50 preps 740980.50	250 preps 740980.250	
Wash Buffer B5 (Concentrate)	6 mL Add 24 mL 96–100 % ethanol	12 mL Add 48 mL 96–100 % ethanol	50 mL Add 200 mL 96–100 % ethanol	
Proteinase K (lyophilized)	6 mg Add 260 μL Proteinase Buffer PB	30 mg Add 1.35 mL Proteinase Buffer PB	75 mg Add 3.35 mL Proteinase Buffer PB	

# 4 Safety instructions

The following components of the NucleoSpin  $^{\otimes}$  DNA FFPE XS kits contain hazardous contents.

Wear gloves and goggles and follow the safety instructions given in this section.

## **GHS classification**

Only harmful features do not need to be labeled with H and P phrases up to 125 mL or 125 g. *Mindergefährliche Eigenschaften müssen bis 125 mL oder 125 g nicht mit H- und P-Sätzen gekennzeichnet werden.* 

Component	Hazard contents	GHS symbol	Hazard phrases	Precaution phrases
Inhalt	Gefahrstoff	GHS-Symbol	H-Sätze	P-Sätze
Paraffin Dissolver	paraffin dissolver 90–100 % Paraffin Dissolver 90–100 %		304	301+310, 311
	CAS 9003-98-9	DANGER <i>GEFAHR</i>		
Proteinase K	Proteinase K 90–100 % Proteinase K 90–100 %	♦ ♦	317, 334	261sh, 280sh, 342+311
	CAS 39450-01-6	WARNING ACHTUNG		

The symbol shown on labels refers to further safety information in this section. Das auf Etiketten dargestellte Symbol weist auf weitere Sicherheitsinformationen dieses Kapitels hin.

#### Hazard phrases

H 304	May be fatal if swallowed and enters airways. Kann bei Verschlucken und Eindringen in die Aternwege tödlich sein.				
H 317	May cause an allergic skin reaction. Kann allergische Hautreaktionen verursachen.				
H 334	May cause allergy or asthma symptoms or breathing difficulties if inhaled. Kann bei Einatmen Allergie, asthmaartige Symptome oder Atembeschwerden verursachen				
Precaution ph	Precaution phrases				
P 261sh	Avoid breathing dust/vapors. Einatmen von Staub/Dampf vermeiden.				
P 280sh	Wear protective gloves/eye protection/face protection. Schutzhandschuhe/Augenschutz tragen.				

- P 301+310 IF SWALLOWED: Immediately call a POISON CENTER/doctor/.... BEI VERSCHLUCKEN: Sofort GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.
- P 311 Call a POISON CENTER/doctor/.... GIFTINFORMATIONSZENTRUM

P 342+311 If experiencing respiratory symptoms: Call a POISON CENTER/doctor/... Bei Symptomen der Atemwege: GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.

For further information please see Material Safety Data Sheets (*www.mn-net.com*). Weiterführende Informationen finden Sie in den Sicherheitsdatenblättern (www.mn-net.com).

# 5 Protocols

NucleoSpin<sup>®</sup> DNA FFPE XS kits offer two different methods for sample deparaffinization. One utilizes the Paraffin Dissolver (included in the kit) and one utilizes xylene or comparable organic solvents (not supplied with the kit). Both methods show same results and effiency.

Deparaffinization with Paraffin Dissolver:	Section 5.1
Deparaffinization with xylene:	Section 5.2

### 5.1 DNA purification from FFPE samples using Paraffin Dissolver

Before starting the preparation:

- Check if Proteinase K and Buffer B5 were prepared according to section 3.
- Check if 96–100 % ethanol is available.
- Set incubator(s) at 60 °C (for paraffin melting) and 90 °C (for decrosslink step).

#### Sample preparation

Insert FFPE section(s) in microcentrifuge tube (not supplied).

For appropriate sample amounts see section 2.4.

#### 1 Deparaffinize sample

Add 400 µL Paraffin Dissolver to the sample.

Incubate **3 min** at **60** °C (to melt the paraffin).

Vortex the sample immediately (at 60 °C) at a vigorous speed to dissolve the paraffin.

<u>Note:</u> Ideally, use a thermoshaker for incubation!

Cool down sample to room temperature.

Make sure that paraffin completely melts during the heat incubation step and mix well after melting to completely dissolve the paraffin.

Insufficient mixing of the heated sample may cause recurrence of solid paraffin particles. Make sure the sample does not comprise more than 15 mg paraffin or adjust the volume of Paraffin Dissolver (see section 2.4).

For samples comprising more than 15 mg paraffin, use  $30 \ \mu L$  Paraffin Dissolver per 1 mg paraffin. If more than  $400 \ \mu L$  Paraffin Dissolver is necessary, place sample in a 2 mL tube (not provided).



Vortex hot sample

#### 2 Lyse sample + 100 µL FL Add 100 µL Buffer FL. Mix Vortex vigorously. Centrifuge at 11,000 x g for 1 min 11,000 x q. Two phases will be formed: a lower (aqueous) phase 1 min and an upper (organic) phase. Tissue material will be transferred to the lower (aqueous) phase. Optional: The upper organic phase can be removed and discarded after centrifugation. Pipette 10 µL Proteinase K solution directly into the lower (aqueous) phase. + 10 µL Proteinase K Mix the aqueous phase by pipetting up and down several times. (Pipette only the lower, aqueous phase up Mix by and down. Avoid mixing lower phase and upper phase pipetting up excessively.) and down (lower phase) Make sure that the Proteinase K is mixed well with the lvsis buffer. If multiple samples are processed, preparation of a Buffer FL/Proteinase K premix is recommended. Add 110 µL of the premix to the reaction tube, mix, and centrifuge to achieve phase formation and to transfer the tissue into the aqueous (lower) phase. Pipette aqueous phase up and down several times in order to disperse the tissue in the lysis buffer. Incubate at room temperature for 3 hours to lyse sample tissue. RT. 3 hours If residual unlysed tissue particles are visible after 3 hours, add additional 10 µL Proteinase K solution and continue digestion for further 3 hours or overnight Note: Yield of amplifiable DNA typically increases with elongated lysis time. During this incubation step protein is digested and DNA is released into solution. Vortex 5 s. Vortex 5 s Set heating block to 90 °C. Convenient stopping point: At this point, the procedure can temporarily be stopped. If pausing, we recommend to store the samples at -20 °C.

3	Decrosslink		
	Add 100 $\mu$ L Decrosslink Buffer D-Link to the tube and vortex gently to mix Buffer D-Link into the aqueous (lower) phase.	Ţ	+ 100 µL D-Link Vortex
	Centrifuge at <b>11,000 x</b> <i>g</i> for <b>30 s</b> to obtain phase formation.	Ò	11,000 x <i>g</i> , 30 s
	Incubate at 90 °C for exactly 30 min.		
	Vortex 5 s and let cool down to room temperature (approx. 2 min).		90 °C, 30 min
	If necessary, spin down briefly to clear the lid (approx. 1 s at 1,000 x g).		Vortex
	<u>Note:</u> This decrosslink step is necessary to remove the crosslinks (chemical modification caused by formalin) from the DNA, which was released into solution by the previous lysis step. Decrosslinked DNA generally shows better performance in downstream applications.		
4	Adjust binding conditions		
	Add <b>200 <math>\mu</math>L ethanol (96–100 %)</b> to the tube and mix by vortexing (2 x 5 s).		+ 200 μL ethanol
	Spin down briefly (approx. 1 s at $1,000 \times g$ ) to achieve complete phase separation.	Ċ	Vortex 1,000 x <i>g</i> ,
	<u>Note:</u> Avoid to centrifuge at much higher g-force, because nucleic acid might precipitate.		1 s
	The ethanol will merge with the aqueous (lower) phase		

only.

#### 5 Bind DNA

For each preparation, take one NucleoSpin<sup>®</sup> DNA FFPE XS Column (green ring) placed in a CollectionTube (2 mL).

Pipette aqueous (lower) phase completely into the NucleoSpin<sup>®</sup> DNA FFPE XS Column.

It is recommended to pipette a volume of 450  $\mu$ L on the column, to ensure that the complete aqueous (lower) phase is transferred (the volume of the aqueous phase is approx. 410  $\mu$ L). Small carry-over of the organic (upper) phase has no negative effect on the binding procedure.

Centrifuge for **30 s** at **2,000 x** *g*. If the solution does not flow through completely, centrifuge for **30 s** at **11,000 x** *g* until the complete solution passed the column.

The recommended centrifugation at  $2,000 \times g$  is more efficient than centrifugation at  $11,000 \times g$ .

Discard Collection Tube with flowthrough and place the column in a new Collection Tube (2 mL).

	· · · · ·	
	Wash and dry silica membrane	6
μ	1 <sup>st</sup> wash	
	Add 400 µL Buffer B5 to the NucleoSpin <sup>®</sup> DNA FFPE XS Column.	
11,000 x <i>g</i> ,	Centrifuge for 30 s at 11,000 x g.	
30 S	Discard Collection Tube with flowthrough and place the column into a new Collection Tube (2 mL).	
	2 <sup>nd</sup> wash	
+ 400 μL B5	Add 400 µL Buffer B5 to the NucleoSpin <sup>®</sup> DNA FFPE XS Column.	
⊘_ 11,000 x <i>g</i> ,	Centrifuge for 2 min at 11,000 x $g$ to dry the membrane.	
2 min	Discard the Collection Tube with flowthrough and place the column into a new nuclease-free microcentrifuge tube (1.5 mL, not provided).	



Load

aqueous

lower) phase

vided).

7	Elute DNA	
	Pipette <b>20 <math>\mu</math>L Buffer BE</b> directly to the center of the silica membrane of the column.	C + 20 μL BE
	Elution volume may be varied from 5–30 μL. For the correlation of elution volume, DNA concentration, and DNA yield see section 2.5.	[₹]
	Centrifuge for 30 s at 11,000 x g.	
8	Optional: Remove residual ethanol	
	Incubate the eluate (20 $\mu L$ ) with open lid for 8 min at 90 °C.	90 °C, 8 min
	See section 2.7 for detailed information and recommendations for removal of residual ethanol.	

# 5.2 DNA purification from FFPE samples with xylene deparaffinization

Before starting the preparation:

- Check if Proteinase K and Buffer B5 were prepared according to section 3.
- Check if 96–100 % ethanol is available.
- Check if xylene (or a similar reagent\*) is available for deparaffinisation.
- Set incubator(s) at 60 °C (for ethanol evaporation) and 90 °C (for decrosslink step).

#### Sample preparation

Insert FFPE section(s) in a microcentrifuge tube (not supplied).

For appropriate sample amounts see section 2.4.

1 Deparaffinize sample

Add <b>1 ml_xviene</b> (or alternative reagent*) to the sample	🌱 1 mL xylene
Incubate at <b>room temperature</b> until the paraffin is completely dissolved (usually approx. 2 min) and vortex vigorously (10 s).	RT 2 min Vortex
Make sure that the paraffin is completely dissolved.	11,000 x <i>g</i> ,
Centrifuge for <b>2 min</b> at <b>11,000 x</b> <i>g</i> .	2 min
<b>Discard the supernatant</b> by pipetting. Do not remove any of the pellet.	Discard supernatant
Add <b>1 mL ethanol (96–100%)</b> to the pellet and vortex (5 s).	°── 1 mL ethanol
Centrifuge for <b>2 min</b> at <b>11,000 x</b> <i>g</i> .	Vortex
<b>Discard the supernatant</b> by pipetting. Do not remove any of the pellet.	
	Discard supernatant
Incubate the open tube at <b>60</b> ° <b>C</b> for <b>3–10 min</b> to dry the pellet.	60 °C,
It is important to evaporate all residual ethanol. Residual ethanol may reduce DNA yield.	3–10 min

<sup>\*</sup> Examples of alternatives to xylene are: d-Limonene (e.g., Roti®-Histol, Hemo-De) or mixtures of isoparafinnic hydrocarbons (e.g., Roticlear<sup>®</sup>, Micro-ClearTM, Neo-Clear<sup>®</sup>).

2	Lyse sample		
	Add <b>100 µL Buffer FL</b> and <b>10 µL Proteinase K</b> to the pellet. Vortex vigorously (5 s).	+ 100 μL FL + 10 μL	
	If multiple samples are processed, preparation of a Buffer FL/Proteinase K premix is recommended. Add 110 $\mu$ L of the premix to the pellet.	Proteinase K Vortex	
	Centrifuge briefly (approx. 1 s at 1,000 x g).		
	Solid section residuals at the tube wall should be flushed back into the solution by pipetting. Pipette solution up and down in order to homogenize sections.		
	Incubate at <b>room temperature</b> for <b>3 hours</b> to lyse sample tissue.	RT,	
	If residual unlysed tissue particles are visible after 3 hours incubation, add additional 10 $\mu L$ Proteinase K solution and continue digestion for further 3 hours or overnight.	3 hours	
	<u>Note:</u> Yield of amplifiable DNA typically increases with elongated lysis time. During this incubation step protein is digested and DNA is released into solution.		
	Vortex tube 5 s.		
	Convenient stopping point: At this point, the procedure can temporarily be stopped. If pausing, we recommend to store the samples at -20 °C.		
	Set heating block to 90 $^\circ\mathrm{C}$ (for subsequent decrosslink step).		
3	Decrosslink		
	Add <b>100 µL Decrosslink Buffer D-Link</b> to the lysate and <b>vortex</b> vigorously (5 s).	+ 100 μL D-Link Vortex	
	Incubate at 90 °C for exactly 30 min.		
	Vortex 5 s let cool down to room temperature (approx. 2 min.	90 °C, 30 min	
	If necessary, spin down briefly to clear the lid (approx. 1 s at $1,000 \times g$ )		
	<u>Note:</u> This decrosslink step is necessary to remove crosslinks (chemical modification caused by formalin) from the DNA which is released in solution by the previous lysis step. Decrosslinked DNA generally shows better performance in downstream applications.		

4	Adjust binding conditions		
	Add <b>200 <math>\mu</math>L ethanol (96–100 %)</b> to the lysate and <b>mix</b> by vortexing (2 x 5 s).		+ 200 μL ethanol
	Spin down briefly to clear the lid (approx. 1 s at 1,000 x $g$ ).	U	Vortex
5	Bind DNA		
	For each preparation, take one NucleoSpin <sup>®</sup> DNA FFPE XS Column (green ring) placed in a Collection Tube (2 mL).		Load lysate
	Pipette lysate up and down two times before loading the lysate.		
	Load the lysate into the column.		
	Centrifuge for <b>30 s</b> at <b>2,000 x</b> <i>g</i> . If the solution does not flow through completely, centrifuge for <b>30 s</b> at <b>11,000 x</b> <i>g</i> until the complete solution passed the column.	Ö	2,000 x <i>g</i> , 30 s
	The recommended centrifugation at 2,000 x g is more efficient than centrifugation at 11,000 x g.		
	Discard Collection Tube with flowthrough and place the column in a new Collection Tube (2 mL).		
6	Wash and dry silica membrane		
	1 <sup>st</sup> wash		+ 400 μL B5
	Add <b>400 µL Buffer B5</b> to the NucleoSpin <sup>®</sup> DNA FFPE XS Column.		
	Centrifuge for <b>30 s</b> at <b>11,000 x</b> <i>g</i> .	Ò	11,000 x <i>g</i> , 30 s
	Discard Collection Tube with flowthrough and place the column into a new Collection Tube (2 mL).		
	2 <sup>nd</sup> wash		
	Add <b>400 µl Buffer B5</b> to the NucleoSpin <sup>®</sup> DNA FFPE XS		+ 400 µl B5
	Column.		11,000 x <i>g</i> ,
	Centrifuge for <b>2 min</b> at <b>11,000 x</b> <i>g</i> to dry the membrane.		2 min
	Discard the Collection Tube with flowthrough and place the column into a nuclease-free Collection Tube (1.5 mL; not provided).		

7	Elute DNA	
	Pipette <b>20 µL Buffer BE</b> directly to the center of the silica membrane of the column.	κοι το με BE
	Elution volume may be varied from 5–30 µL. For the correlation of elution volume, DNA concentration, and DNA yield see section 2.5.	A
	Centrifuge for <b>30 s</b> at <b>11,000 x</b> <i>g</i> .	
8	Optional: Remove residual ethanol	
	Incubate the eluate (20 μL) with open lid for <b>8 min</b> at <b>90 °C</b> .	90 °C, 8 min
	See section 2.7 for detailed information and recommendations for removal of residual ethanol.	

# 6 Appendix

## 6.1 Comments on DNA quality and quantity

Due to tissue fixation, nucleic acids in FFPE samples are commonly fragmented and chemically modified by formaldehyde. Formaldehyde modifications of DNA cannot be detected by standard quality control assays, such as gel electrophoresis, spectrophotometry, fluorometry, or microfluidics analysis. However, efficiency of enzymatic reactions (e.g., PCR) with chemically modified DNA is significantly decreased.

Affected DNA analysis methods and applications are for example:

- Spectrophotometry (e.g., absorption measurement A<sub>230</sub>, A<sub>260</sub>, A<sub>280</sub>)
- Fluorometry (e.g., RiboGreen<sup>®</sup>)
- Denaturing agarose gel electrophoresis
- Mirofluidics analysis (e.g., Agilent 2100 Bioanalyzer, BioRad´s Experion Automated Electrophoresis System)
- PCR
- Array analysis (e.g., DNA microarrays)

The following aspects should be considered when applying one of the listed methods, especially when comparing efficiency of different DNA isolation and decrosslink procedures or the usability of the isolated DNA:

- A high DNA yield, as determined by A<sub>260</sub> readings or by fluorescent dye (e.g., PicoGreen<sup>®</sup>) analysis does not necessarily result in good performance of the DNA in a PCR. DNA may be highly degraded (i.e., smaller fragments than the PCR target) or insufficiently decrosslinked.
- Low or no DNA yield as determined by A<sub>260</sub> readings will most likely result in poor PCR results, but it is still possible to achieve a good performance. There may be a small amount DNA which is decrosslinked sufficiently and shows good reactivity.
- DNA of high molecular weight does not guarantee a good amplifiability in PCR or reactivity in other enzymatic reactions. DNA may be insufficiently decrosslinked although it has high molecular weight.
- DNA of low molecular weight, i.e. highly degraded DNA with fragment sizes exclusively below 200 nucleotides will certainly not enable amplification of fragments exceeding this size. However, it is still likely that small sized target sequences (e.g., 80–150 bp) can be amplified successfully, especially if the DNA is well decrosslinked.

Neither DNA yield, molecular weight, absorbance ratios, nor size distribution can reliably predict the performance in downstream PCR applications, especially if different purification and decrosslinking systems are compared.

The major quality indicator for DNA isolated from FFPE samples is its performance in the intended downstream application.

Problem	Possible cause and suggestions		
DNA is	Poor sample quality		
degraded/no DNA obtained	<ul> <li>Sample quality has a high impact on quality and amount of the DNA.</li> </ul>		
	Reagents not applied or restored properly		
	<ul> <li>Reagents not properly restored. Add the indicated volume of ethanol to Buffer B5 Concentrate and mix. Reconstitute and store Proteinase K according to instructions given in section 3.</li> </ul>		
	<ul> <li>Sample and reagents have not been mixed completely. Always vortex vigorously after each reagent has been added.</li> </ul>		
	<ul> <li>No ethanol has been added after lysis. Binding of DNA to the silica membrane is only effective in the presence of ethanol.</li> </ul>		
	Kit storage		
	• Reconstitute and store Proteinase K according to instructions given in section 3.		
	• Store kit components as described in section 3.		
Poor DNA quality or yield	<ul> <li>Keep bottles tightly closed in order to prevent evaporation or contamination.</li> </ul>		
	lonic strength and pH influence $A_{\rm 260}$ absorption as well as ratio $A_{\rm 260}/A_{\rm 280}$		
	<ul> <li>For absorption measurement, use 5 mM Tris pH 8.5 as diluent. Please also see:         <ul> <li>Manchester, K. L. 1995. Value of A<sub>260</sub>/A<sub>280</sub> ratios for measurement of purity of nucleic acids. Biotechniques 19, 208-209.</li> <li>Wilfinger, W. W., Mackey, K. and Chomczyski, P. 1997. Effect of pH and ionic strength on the spectrophotometric assessment of nucleic acid purity. Biotechniques 22, 474-481.</li> </ul> </li> </ul>		
	Proteinase K digestion time		
	<ul> <li>Depending of the nature of the sample, an optimal digestion time from 3 to 16 hours has to be determined empirically. If residual unlysed tissue is still visible after 3 h continue the incubation for up to 16 hours. After the first 3 h incubation, additional Proteinase K</li> </ul>		

# 6.2 Troubleshooting

may be added to the sample.

	Sample material			
Clogged NucleoSpin <sup>®</sup> DNA FFPE XS Column/	<ul> <li>Too much starting material was used. Overloading may lead to a decrease of DNA yield. Reduce the quantity of sample material or use larger volumes of Paraffin Dissolver and/or Lysis Buffer FL.</li> </ul>			
Poor DNA quality or yield	<ul> <li>Insufficient disruption and / or homogenization of starting material. Perform only an overnight incubation, if the tissue was not completely digested after 3 hours.</li> </ul>			
	Carry-over of ethanol or salt			
	• Do not let the flowthrough touch the column outlet after the second wash with Buffer B5. Be sure to centrifuge at the corresponding speed for the respective time in order to remove ethanolic Buffer B5 completely.			
Suboptimal performance of DNA in	<ul> <li>Check if Buffer B5 has been equilibrated to room temperature before use. Washing at lower temperatures decreases efficiency of salt removal by Buffer B5.</li> </ul>			
experiments	<ul> <li>Depending on the robustness of the used PCR system, PCR might be inhibited if too much eluate is applied. Use less eluate as template.</li> </ul>			
	Store isolated DNA properly			
	• Eluted DNA should always be kept on ice for optimal stability since possible traces of DNases will degrade the isolated DNA.			
	Silica abrasion from the membrane			
Discrepancy between A <sub>260</sub> quantification values and PCR quantification values	• Due to the typically low DNA content in small FFPE samples and the resulting low total amount of isolated DNA, a DNA quantification via $A_{260}$ absorption measurement is often hampered due to the low sensitivity of the absorption measurement. When performing absorption measurements close to the detection limit of the photometer, the measurement may be influenced by minor amounts of silica abrasion. In order to prevent incorrect $A_{260}$ -quantification of small DNA amounts centrifuge the eluate for 30 s at >11.000 x g and take an aliquot for measurement without disturbing any sediment. Alternatively, use a silica abrasion insensitive RNA quantification method (e.g., PicoGreen <sup>®</sup> fluorescent dye).			
	Measurement not in the range of photometer detection limit			
Unexpected A <sub>260</sub> /A <sub>280</sub> ratio	• In order to obtain a significant A <sub>260</sub> /A <sub>280</sub> ratio it is necessary that the initially measured A <sub>260</sub> and A <sub>280</sub> values are significantly above the detection limit of the photometer used. An A <sub>280</sub> value close to the background noise of the photometer will cause unexpected A <sub>260</sub> /A <sub>280</sub> ratios.			

# 6.3 Ordering information

Product	REF	Pack of
NucleoSpin <sup>®</sup> DNA FFPE XS	740980.10/.50/.250	10/50/250
NucleoSpin <sup>®</sup> totalRNA FFPE	740982.10/.50/.250	10/50/250
NucleoSpin <sup>®</sup> totalRNA FFPE XS	740969.10/.50/.250	10/50/250
NucleoSpin <sup>®</sup> Tissue XS	740901.10/.50/.250	10/50/250
Paraffin Dissolver	740968.25	25 mL
Paraffin Dissolver blue	740343.60	60 mL
Collection Tubes (2 mL)	740600	1000
Decrosslink Buffer D-Link	740979.30	30 mL

Visit *www.mn-net.com* for more detailed product information.

## 6.4 Product use restriction / warranty

**NucleoSpin® DNA FFPE XS** kit components are intended, developed, designed, and sold FOR RESEARCH PURPOSES ONLY, except, however, any other function of the product being expressly described in original MACHEREY-NAGEL product leaflets.

MACHEREY-NAGEL products are intended for GENERAL LABORATORY USE ONLY! MACHEREY-NAGEL products are suited for QUALIFIED PERSONNEL ONLY! MACHEREY-NAGEL products shall in any event only be used wearing adequate PROTECTIVE CLOTHING. For detailed information please refer to the respective Material Safety Data Sheet of the product! MACHEREY-NAGEL products shall exclusively be used in an ADEQUATE TEST ENVIRONMENT. MACHEREY-NAGEL does not assume any responsibility for damages due to improper application of our products in other fields of application. Application on the human body is STRICTLY FORBIDDEN. The respective user is liable for any and all damages resulting from such application.

DNA/RNA/PROTEIN purification products of MACHEREY-NAGEL are suitable for *IN VITRO*-USES ONLY!

ONLY MACHEREY-NAGEL products specially labeled as IVD are also suitable for *IN VITRO*diagnostic use. Please pay attention to the package of the product. *IN VITRO*-diagnostic products are expressly marked as IVD on the packaging.

IF THERE IS NO IVD SIGN, THE PRODUCT SHALL NOT BE SUITABLE FOR *IN VITRO*-DIAGNOSTIC USE!

ALL OTHER PRODUCTS NOT LABELED AS IVD ARE NOT SUITED FOR ANY CLINICAL USE (INCLUDING, BUT NOT LIMITED TO DIAGNOSTIC, THERAPEUTIC AND/OR PROGNOSTIC USE).

No claim or representations is intended for its use to identify any specific organism or for clinical use (included, but not limited to diagnostic, prognostic, therapeutic, or blood banking). It is rather in the responsibility of the user or - in any case of resale of the products - in the responsibility of the reseller to inspect and assure the use of the DNA/RNA/protein purification products of MACHEREY-NAGEL for a well-defined and specific application.

MACHEREY-NAGEL shall only be responsible for the product specifications and the performance range of MN products according to the specifications of in-house quality control, product documentation and marketing material.

This MACHEREY-NAGEL product is shipped with documentation stating specifications and other technical information. MACHEREY-NAGEL warrants to meet the stated specifications. MACHEREY-NAGEL's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Supplementary reference is made to the general business terms and conditions of MACHEREY-NAGEL, which are printed on the price list. Please contact us if you wish to get an extra copy.

There is no warranty for and MACHEREY-NAGEL is not liable for damages or defects arising in shipping and handling (transport insurance for customers excluded), or out of accident or improper or abnormal use of this product; defects in products or components not manufactured by MACHEREY-NAGEL, or damages resulting from such non-MACHEREY-NAGEL components or products.

MACHEREY-NAGEL makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE

WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, REPRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO MACHEREY-NAGEL PRODUCTS.

In no event shall MACHEREY-NAGEL be liable for claims for any other damages, whether direct, indirect, incidental, compensatory, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of MACHEREY-NAGEL products to perform in accordance with the stated specifications. This warranty is exclusive and MACHEREY-NAGEL makes no other warranty expressed or implied.

The warranty provided herein and the data, specifications and descriptions of this MACHEREY-NAGEL product appearing in MACHEREY-NAGEL published catalogues and product literature are MACHEREY-NAGEL's sole representations concerning the product and warranty. No other statements or representations, written or oral, by MACHEREY-NAGEL's employees, agent or representatives, except written statements signed by a duly authorized officer of MACHEREY-NAGEL are authorized; they should not be relied upon by the customer and are not a part of the contract of sale or of this warranty.

Product claims are subject to change. Therefore please contact our Technical Service Team for the most up-to-date information on MACHEREY-NAGEL products. You may also contact your local distributor for general scientific information. Applications mentioned in MACHEREY-NAGEL literature are provided for informational purposes only. MACHEREY-NAGEL does not warrant that all applications have been tested in MACHEREY-NAGEL laboratories using MACHEREY-NAGEL products. MACHEREY-NAGEL does not warrant the correctness of any of those applications.

Last updated: 07/2010, Rev. 03

Please contact: MACHEREY-NAGEL GmbH & Co. KG Tel.: +49 24 21 969-270 tech-bio@mn-net.com

DyNAmo is a trademark of Finnzymes Oy

Trademarks:

LighCycler is a registeres trademark of a member of the Roche Group

Micro-Clear is a registered trademark of Micron Environmental Industries

Neo-Clear is a registered trademark of Merck KGaA

NucleoSpin is a registered trademark of MACHEREY-NAGEL GmbH & Co KG

PicoGreen is a registered trademark of Molecular Probes Inc.

Roticlear is a registered trademark of CARL ROTH GmbH & Co KG

Roti is a registered trademark of CARL ROTH GmbH & Co KG

SYBR is a registered trademark of Molecular Probes, Inc.

All used names and denotations can be brands, trademarks, or registered labels of their respective owner – also if they are not special denotation. To mention products and brands is only a kind of information (i.e., it does not offend against trademarks and brands and can not be seen as a kind of recommendation or assessment). Regarding these products or services we can not grant any guarantees regarding selection, efficiency, or operation.



#### www.mn-net.com

# **MACHEREY-NAGEL**

(MN) EN BO 1001 EN BO 13485 CEPTIFIED

DE/International: 
 Tel.:
 +49 24 21 969-0
 Tel.:
 +41 62 388 55 00

 Fax:
 +49 24 21 969-199
 Fax:
 +41 62 388 55 05
 E-mail: info@mn-net.com

CH: E-mail: sales-ch@mn-net.com E-mail: sales-fr@mn-net.com

MACHEREY-NAGEL GmbH & Co. KG · Neumann-Neander-Str. 6-8 · 52355 Düren · Germany FR: Tel.: +33 388 68 22 68 Fax: +33 388 51 76 88

US:

Tel.: +1 484 821 0984 Fax: +1 484 821 1272 E-mail: sales-us@mn-net.com

A039236/8020.5