

RNase A

1 Contents

RNase A		
REF	740505	740505.50
RNaseA (lyophilized)	100 mg	50 mg

2 Product description

The RNase A is primarily intended to be used with NucleoSpin® and NucleoBond® kits. See individual protocols for detailed information. However, if the RNase A is used for another than the recommended purpose performance may differ considerably from its expected one.

3 RNase A specifications

Source: Bovine pancreas

Form: 100 mg / or 50 mg white freeze-dried material

RNase activity: 120 U/mg

4 Storage conditions and preparation of working solution

Lyophilized RNase A can be stored at room temperature (18–25 °C) for at least one year.

A: Dissolved in water: Store the RNase A solution at 4 °C for up to 3 months. For longer storage, the RNase A solution should be divided into small aliquots and stored at -20 °C.

B: Dissolved in Resuspension Buffer A1 / S1 / S1-EF / RES / RES-EF / RES-BAC: Buffer containing RNase A can be stored at 4 °C for at least 6 months.

RNase A		
REF	740505	740505.50
RNaseA (lyophilized)	100 mg	50 mg
	Dissolve in H2O or directly in Resuspension Buffer A1 / S1 / S1-EF / RES / RES-EF / RES-BAC*	

* Please refer to the individual protocols for detailed information.


5 Safety instructions

RNase A contains hazardous contents. Wear gloves and goggles!

GHS classification

Only harmful features must not be labeled with H and P phrases until 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
RNase A	RNase 90–100% RNase 90–100% CAS 9001-99-4	 DANGER GEFAHR	317, 334	261, 272, 280, 302+352, 304+340, 333+313, 342+311, 363

Hazard phrases

- H317 May cause an allergic skin reaction.
Kann allergische Hautreaktionen verursachen.
- H334 May cause allergy or asthma symptoms or breathing difficulties if inhaled.
Kann bei Einatmen Allergie, asthmaartige Symptome oder Atembeschwerden verursachen.

Precaution phrases

- P261 Avoid breathing vapours.
Dampf nicht einatmen.
- P272 Contaminated work clothing should not be allowed out of the workplace.
Kontaminierte Arbeitskleidung nicht außerhalb des Arbeitsplatzes tragen.
- P280 Wear protective gloves/protective clothing/eye protection/face protection.
Schutzhandschuhe/Schutzkleidung/Augenschutz/Gesichtsschutz tragen.
- P302+352 IF ON SKIN: Wash with plenty of water/...
BEI BERÜHRUNG MIT DER HAUT: Mit viel Wasser/... waschen.
- P304+340 IF INHALED: Remove person to fresh air and keep comfortable for breathing.
BEI EINATMEN: Die Person an die frische Luft bringen und für ungehinderte Atmung sorgen.
- P333+313 If skin irritation or rash occurs: Get medical advice/attention.
Bei Hautreizung oder -ausschlag: Ärztlichen Rat einholen/ärztliche Hilfe hinzuziehen.
- P337+313 If eye irritation persists: Get medical advice/attention.
Bei anhaltender Augenreizung: Ärztlichen Rat einholen/ärztliche Hilfe hinzuziehen.
- P342+311 If experiencing respiratory symptoms: Call a POISON CENTER/ doctor/...
Bei Symptomen der Atemwege: GIFTINFORMATIONSZENTRUM/Arzt/... anrufen.
- P363 Wash contaminated clothing before reuse.
Kontaminierte Kleidung vor erneutem Tragen waschen.

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



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.

Product use restriction / warranty

RNase A is developed, designed, and sold for research purpose only. All goods have to be checked immediately upon receipt. Damage claims are only acceptable in writing within 8 working days of receipt of all the goods. In case of legitimate claims the buyer can only require replacement of the goods. If replacement is not possible, the buyer has the right to choose between alternative products with same value or refund. The buyer cannot claim further compensation. All returns must first be authorized by us in writing. We offer a one year warranty from date of delivery that our products will conform to applicable specifications set forth in the product specifications if not sold to persons set forth in § 13 of German Civil Code. In such a case, the provisions of the German Civil Code shall be valid.

This warranty does not cover defects in or damage to the products which are due to improper installation or maintenance, misuse, neglect or any use other than ordinary commercial application.

Any discharge from liability will be void if a defect results from a heavily negligent or intentional breach of contract or if the buyer will be bodily injured as a result from a negligent or intentional breach.

The seller shall not be liable for the goods being fit for a particular purpose unless otherwise agreed upon, to which the buyer intends to put them.

This warranty is strictly exclusive. Any further damage compensation is impossible.

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; REPRODUCIBILITY; DURABILITY; FITNESS FOR A PARTICULAR PURPOSE OR USE; MERCHANTABILITY; CONDITION; OR ANY OTHER MATTER WITH RESPECT TO MACHEREY-NAGEL PRODUCTS OR THE SALE OF MACHEREY-NAGEL PRODUCTS.

If not superseded by these terms the current terms and conditions for the sale of goods (export version) are additionally valid.

**This product distributed by Clontech Laboratories, Inc.
A Takara Bio Company**

1290 Terra Bella Ave., Mountain View, CA 94043
www.clontech.com

For ordering information

1.800.662.2566
orders@clontech.com

For technical support

1.800.662.2566
tech@clontech.com



For research use only. Not for resale. Unless otherwise indicated, all trademarks are the property of Clontech Laboratories, Inc. ©2015