



# Protein Quantification Assay

## User Manual

*Visit our website  
for more details!  
click here...*



This product is 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

PT5032-1  
PROX3674



September 2010/Rev.02

**MACHERY-NAGEL**



# Protein Quantification Assay

## Protocol-at-a-glance (Rev.02)

		Microplate assay	Semi-micro cuvette assay	Micro cuvette assay	Low volume assay
1	<b>Prepare BSA reference protein dilution series</b>	Dispense <b>50 µL</b> PSB per tube #2 – #7	Dispense <b>250 µL</b> PSB per tube #2 – #7	Dispense <b>50 µL</b> PSB per tube #2 – #7	Dispense <b>20 µL</b> PSB per tube #2 – #6
		Pipette <b>50 µL</b> BSA stock solution into tube #2; then <b>50 µL</b> from #2 into #3 etc.*	Pipette <b>250 µL</b> BSA stock solution into tube #2; then <b>250 µL</b> from #2 into #3 etc.*	Pipette <b>50 µL</b> BSA stock solution into tube #2; then <b>50 µL</b> from #2 into #3 etc.*	Pipette <b>20 µL</b> BSA stock solution into tube #2; then <b>20 µL</b> from #2 into #3 etc.*
	<b>Dilution series sufficient for</b>	<i>Two calibration curves</i>	<i>One calibration curve</i>	<i>One calibration curve</i>	<i>Two calibration curves</i>
2	<b>Dispense dilution series</b>	20 µL	200 µL	40 µL	7.5 µL
3	<b>Dispense your protein sample</b>	20 µL (1 – 60 µL)	200 µL (10 – 600 µL)	40 µL (1 – 120 µL)	7.5 µL
4	<b>Fill up dilution series and sample with PSB</b>	40 µL (final vol. 60 µL)	400 µL (final vol. 600 µL)	80 µL (final vol. 120 µL)	–
5	<b>Add Quantification Reagent QR</b>	40 µL	400 µL	80 µL	5 µL
6	<b>Incubate</b>	30 min at room temperature			
7	<b>Measure light extinction</b>	At 570 nm (530 nm – 700 nm)			
8	<b>Calculate protein concentration</b>	Make sure that the signal of your sample lies within the range of the calibration curve.			

\* Keep tube #7 as BLANK – Do not add 50 µL from tube #6 into tube #7!

## Table of contents

1	Components	4
1.1	Kit contents	4
1.2	Consumables and equipment to be supplied by user	5
2	Product description	6
2.1	The basic principle	6
2.2	Kit specifications	6
2.3	Handling, preparation, and storage of starting materials	9
2.4	Calibration curves	10
2.5	Recommended sample volumes	10
2.6	Alternative wavelengths for extinction measurement	12
3	Storage conditions and preparation of working solutions	13
4	Safety instructions – risk and safety phrases	14
5	Protocols	15
5.1	Microplate assay procedure	15
5.2	Semi-micro cuvette assay procedure	18
5.3	Microcuvette assay procedure	21
5.4	Low volume assay procedure	24
6	Appendix	27
6.1	Troubleshooting	27
6.2	Ordering information	28
6.3	References	29
6.4	Product use restriction/warranty	29

# 1 Components

## 1.1 Kit contents

<b>Protein Quantification Assay</b>		
<b>REF</b>	<b>50 assays</b>	<b>250 assays</b>
	<b>740967.50</b>	<b>740967.250</b>
Protein Solving Buffer PSB	7.5 mL	30 mL
BSA (Bovine Serum Albumin; reference protein)*	1 mg	2 x 1 mg
Quantification Reagent QR	5 mL	20 mL
User Manual	1	1

\* For preparation of working solutions and storage conditions see section 3.

## 1.2 Consumables and equipment to be supplied by user

### Consumables

- Microplates, flat-bottom (e.g., UV-Star Microtiter plate, 96-well, F-bottom, Greiner bio-one REF 655801; similar non-UV transparent microtiter plates are also suitable) **or** semi-micro cuvettes (e.g., Plastibrand 1.5 mL semi-micro disposable cuvettes, Brand REF 759115) or micro-cuvettes (e.g., Plastibrand UV-Cuvette micro, Brand, REF 759220).
- 1.5 mL microcentrifuge tubes (to prepare dilution series for the calibration curve and to set up reactions when following the semi-micro cuvette assay procedure)
- Disposable pipette tips

### Equipment

- Manual pipettors
- Centrifuge for microcentrifuge tubes (to clean microcentrifuge lids if necessary)
- Vortex mixer
- Mixer or shaker for microplates
- Photometer set to 570 nm (570 nm is recommended, other wavelength settings in the range of 530 – 700 nm are also suitable), either for microplates (microplate assay procedure), for semi-micro/microcuvettes (semi-micro cuvette and/micro cuvette assay procedure) or for low volume analysis (e.g., NanoDrop (Thermo Scientific), NanoVue (GE Healthcare), or NanoPhotometer™ (Implen)).
- Personal protection equipment (e.g., lab coat, gloves, goggles)

## 2 Product description

### 2.1 The basic principle

The **Protein Quantification Assay** is a convenient and reliable kit for the determination of protein concentration in samples typically used for SDS-PAGE (sodium dodecylsulfate polyacrylamide gel electrophoresis). It is mainly designed for proteins solved in Protein Solving Buffer PSB or Protein Loading Buffer PLB (components of NucleoSpin® RNA/Protein and NucleoSpin® TriPrep), but will also work with proteins solved in buffer as described by Laemmli (1970), or similar. These protein sample buffers usually contain SDS, a reducing agent, dye, and a component to increase the buffer density. The majority of protein quantification assays\* are either influenced by or incompatible with SDS, reducing agents, or dyes commonly present in protein sample buffers. The **Protein Quantification Assay** however, is well suited for such buffer systems. It is a fast and sensitive assay, based on a modification of a protocol described by Karlsson *et al.* (1994). The samples are mixed with Protein Solving Buffer PSB and subsequently incubated for 30 minutes with Quantification Reagent QR. After incubation light extinction is measured photometrically. Light extinction is caused by turbidity appearing after addition of Quantification Reagent QR. The protein concentration is determined in reference to a BSA (Bovine Serum Albumin) calibration curve (BSA is provided with the **Protein Quantification Assay**).

### 2.2 Kit specifications

- **Protein Quantification Assay** allows the determination of protein concentration in samples containing up to 10% SDS and comprising reducing agent (e.g.,  $\beta$ -mercaptoethanol (BME), dithiothreitol (DTT), dithioerythritol (DTE) or tris-(2-carboxyethyl) phosphine hydrochloride (TCEP)), buffering salts (e.g., TRIS or BIS-TRIS), dye (bromphenol blue), and a component to create a high density of the solution (e.g., glycerol or sucrose).
- **Protein Quantification Assay** is designed for the determination of protein concentration in samples with low nucleic acid concentration, as obtained with NucleoSpin® RNA/Protein or NucleoSpin® TriPrep. For samples rich in nucleic acids the quantification is less accurate.
- 

---

\* For example: Coomassie Brilliant Blue G-250, Bradford 1979; copper tartrate solution and Folin reagent, Lowry *et al.* 1951;  $\text{Cu}^{2+}/\text{Cu}^{1+}$  - BCA interaction, Smith *et al.* 1985.

- **Protein Quantification Assay** is suited for samples comprising protein solved in buffers, commonly used for SDS-PAGE (e.g., Laemmli buffer). Accuracy depends on nucleic acid content of the sample. For typical cultured cells (e.g., HeLa) accuracy is affected by approximately 5 – 20% due to nucleic acid content\*.
- The kit REF **740967.50** is sufficient for 50 protein determinations plus six calibration curves with seven calibration points each (approx. 100 reactions in total), according to the microplate assay. Alternatively the kit is sufficient for approx. 10 reactions according to the semi-micro cuvette assay (three protein determinations plus seven calibration points).
- The kit REF **740967.250** is sufficient for 250 protein determinations plus 25 calibration curves with seven calibration points each (approx. 450 reactions in total), according to the microplate assay. Alternatively, the kit is sufficient for approx. 50 reactions according to the semi-micro cuvette assay (26 protein determinations plus three calibration curves with seven calibration points each).
- Following the **microplate assay procedure** the kit allows the determination of protein amount (exemplary BSA) in the range of 0.6 – 20 µg per assay provided in a standard volume of 20 µL Protein Solving Buffer PSB (alternatively 1 – 60 µL). This corresponds to a protein concentration of 30 – 1000 ng/µL. This concentration range can be expanded to 10 – 20,000 ng/µL if alternative sample volumes (1 – 60 µL) are applied.
- Following the **semi-micro cuvette assay procedure** the kit allows the determination of protein (exemplary BSA) amount in the range of 6 – 200 µg per assay provided in a standard volume of 200 µL Protein Solving Buffer PSB. This corresponds to a protein concentration of 30 – 1000 ng/µL.

\* One microgram DNA causes ca. 50 – 70% of the extinction signal caused by one microgram protein (BSA).  
One microgram RNA causes ca. 10 – 40% of the extinction signal caused by one microgram protein (BSA).

DNA, RNA, and protein content of a typical cell and influence on the protein quantification:				
Molecule	Content per cell	Content per one million cells	Extinction signal obtained with the Protein Quantification Assay relative to the reference protein BSA	Extinction signal obtained relative to total protein
DNA	6 pg	6 µg	~ 50 – 70%	3 – 5%
RNA	10 – 30 pg	10 – 30 µg	~ 10 – 40%	1 – 12%
Protein	100 – 200 pg	100 – 200 µg	~ 100%	100%
				Total: 104 – 117%

Signal obtained from total cell extract containing RNA and DNA, relative to nucleic acid free total protein: 104% – 117%.

**Table 1: Kit specifications at a glance\***

<b>Protein Quantification Assay</b>	
Sample size	1 – 600 µL containing 0.6 – 200 µg protein (BSA equivalents)
Microplate assay	0.6 – 20 µg protein (BSA equivalents) in 20 µL, corresponding to 30 – 1000 ng/µL
Semi-micro cuvette assay	6 – 200 µg protein (BSA equivalents) in 200 µL, corresponding to 30 – 1000 ng/µL
Sample type	Protein solved in Protein Solving Buffer PSB, Laemmli buffer or equivalent, preferable free of nucleic acids
Protein concentration	Approx. 30 – 1,000 ng/µL (standard range) or Approx. 10 – 20,000 ng/µL (extended range)
Correlation coefficient	0.97 – 1.00
Wavelength for light extinction measurement	570 nm (530 – 700 nm)
Time	Approx. 40 min

\* Kit specifications vary depending on the type of assay. Please find more detailed information in the tables below:

Type of assay	Required sample volume	Protein amount per assay	Determinable protein concentration
Microplate	20 µL (1 – 60 µL)	0.6 – 20 µg	30 – 1000 ng/µL
Semi-micro cuvette	200 µL (10 – 600 µL)	6 – 200 µg	30 – 1000 ng/µL
Micro cuvette	40 µL (1 – 120 µL)	1.2 – 40 µg	30 – 1000 ng/µL
Low volume	7.5 µL	0.47 – 7.5 µg	60 – 1000 ng/µL

Type of assay	740967.50			740967.250		
	Protein determination	Calibration curves (7 points per curve)	Total number of reactions	Protein determination	Calibration curves (7 points per curve)	Total number of reactions
Microplate	50	6	Approx. 100	250	25	Approx. 450
Semi-micro cuvette	3	1	Approx. 10	26	3	Approx. 50
Micro cuvette	35	3	Approx. 55	130	15	Approx. 235
Low volume	800	25	Approx. 1000	3000	70	Approx. 3500



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

After dissolving protein in Protein Solving Buffer PSB or Protein Loading Buffer PLB (with or without Reducing Agent TCEP), Laemmli buffer, or analogs, freeze your protein samples for long term storage or keep samples at 4°C for short term storage. Before use, make sure that the samples are free of precipitates. If necessary heat to approximately 30°C in order to dissolve any possible SDS precipitate. Subsequently, spin sample briefly to remove any further insoluble matter.

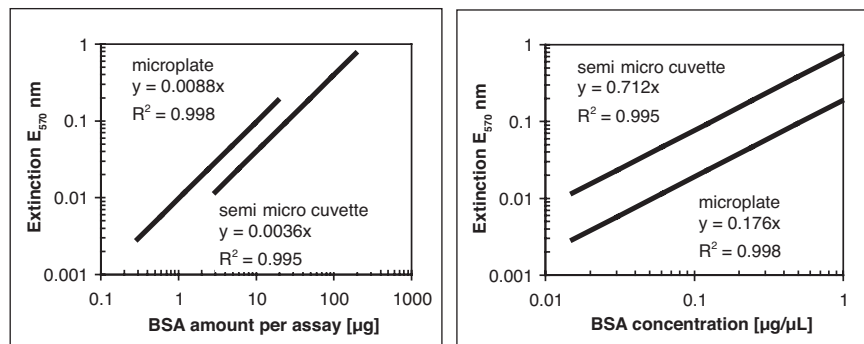
Protein samples obtained with NucleoSpin® RNA/Protein or NucleoSpin® TriPrep and dissolved in Protein Solving Buffer PSB or Protein Loading Buffer PLB (with or without Reducing Agent TCEP), are optimal for determination of protein concentration with the **Protein Quantification Assay**.

Quantification of protein samples obtained by boiling cells or tissue directly in either PSB or PLB (with or without Reducing Agent TCEP), Laemmli buffer, or analogs is possible, but the measurement may be less accurate due to the presence of nucleic acids, which interfere with the assay. The extent of interference depends on the content of protein and nucleic acid in the sample. Many samples, like, for example, cultured HeLa cells or liver tissue, contain much more protein than nucleic acid and thus nucleic acids cause only small interference (see footnote page 7).

*Wear gloves at all times during the handling to reduce risk of sample contamination with skin keratins.*

## 2.4 Calibration curves

Reference protein (BSA) dilution series give good correlations with measured light extinction. Typical correlation coefficients of 0.97 – 1.00 are obtained in the range of approx. 0.03 – 1  $\mu\text{g}/\mu\text{L}$  BSA concentration. BSA concentration versus extinction and BSA amount versus extinction are shown in Figure 1.



**Figure 1: Correlation between BSA amount and extinction signal as well as between BSA concentration and extinction signal.**

For the microplate assay BSA was supplied in 20  $\mu\text{L}$ ; path length for extinction measurement was 3 mm. For the semi-micro cuvette assay BSA was supplied in 200  $\mu\text{L}$ ; path length for extinction measurement was 10 mm.

## 2.5 Recommended sample volumes

As guidance, follow the recommendations of Table 2 – Table 5 to choose an appropriate volume of your sample for measuring. For the initial determination of protein concentration in samples containing hard-to-estimate protein amounts, measurement of multiple sample volumes (e.g., 2  $\mu\text{L}$ , 5  $\mu\text{L}$ , 50  $\mu\text{L}$ ) is recommended. This will increase the probability that one of the measured protein amounts lies within the range of the calibration curve.

For protein samples obtained with NucleoSpin® RNA/Protein or NucleoSpin® TriPrep, see the respective user manual for a first estimation of the protein yield.

**Table 2: Microplate assay –  
Recommended sample volumes for protein quantification**

Expected protein concentration	Recommended sample volume	Protein amount per well
0.01 – 0.33 µg/µL	60 µL	0.6 – 20 µg
0.03 – 1.0 µg/µL	20 µL	0.6 – 20 µg
0.6 – 20 µg/µL	1 µL	0.6 – 20 µg

**Table 3: Semi-micro cuvette assay –  
Recommended sample volumes for protein quantification**

Expected protein concentration	Recommended sample volume	Protein amount
0.01 – 0.33 µg/µL	600 µL	6 – 200 µg
0.03 – 1.0 µg/µL	200 µL	6 – 200 µg
0.6 – 20 µg/µL	10 µL	6 – 200 µg

**Table 4: Microcuvette assay –  
Recommended sample volumes for protein quantification**

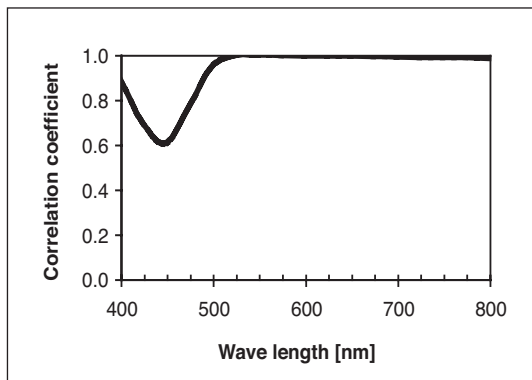
Expected protein concentration	Recommended sample volume	Protein amount per microcuvette
0.01 – 0.33 µg/µL	120 µL	1.2 – 40 µg
0.03 – 1.0 µg/µL	40 µL	1.2 – 40 µg
1.2 – 40 µg/µL	1 µL	1.2 – 40 µg

**Table 5: Low volume assay –  
Recommended sample volumes for protein quantification**

Expected protein concentration	Recommended sample volume	Protein amount per microcuvette
0.06 – 1 µg/µL	7.5 µL	0.47 – 7.5 µg

## 2.6 Alternative wavelengths for extinction measurement

A wavelength in the range of 530 – 700 nm is recommended for light extinction measurements. Figure 2 shows the dependency of correlation coefficient on the wavelength, used for light extinction measurement.



**Figure 2: Dependency of correlation coefficient on the wavelength, used for extinction measurement.**

Light extinction of BSA samples in the range of 0.3 – 20  $\mu\text{g}$  was measured for wavelength between 400 nm and 800 nm. The correlation coefficient was calculated from the BSA amount per assay (0.3 – 20  $\mu\text{g}$  per assay) and corresponding extinction signal.

### 3 Storage conditions and preparation of working solutions

**Attention:**

*Quantification Reagent QR contains hydrochloric acid. Wear gloves and goggles!*

- All kit components should be stored at room temperature (18 – 25°C). Storage at lower temperatures may cause precipitation in the Protein Solving Buffer PSB. Kit components are stable up to one year.

Before starting the **Protein Quantification Assay** prepare the following:


- Dissolve the reference protein (BSA, 1 mg) in 1 mL Protein Solving Buffer PSB to obtain a 1 mg/mL BSA stock solution. Freeze BSA stock solution for long term storage; for short term storage keep solution at 4°C. If necessary, dissolve any precipitate by heating the reference solution (approx. 30°C) before use. BSA stock solution (1 mg/mL BSA in PSB) is stable at -20°C to +20°C for six months.

Protein Quantification Assay		
	50 assays	250 assays
REF	740967.50	740967.250
BSA (reference protein)	1 mg add 1 mL PSB	2 x 1 mg add 1 mL PSB to each vial

## 4 Safety instructions – risk and safety phrases

Quantification Reagent QR contains hydrochloric acid.

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

Component	Hazard contents	Hazard symbol		Risk phrases	Safety phrases
QR	Hydrochloric acid <20%	 Xi*	Irritating to eyes, respiratory system and skin	R 36/37/38	S 26-45

### Risk phrases

R 36/37/38 Irritating to eyes, respiratory system and skin

### Safety phrases

- S 26 In case of contact with eyes, rinse immediately with plenty of water and seek medical advice
- S 45 In case of accident or if you feel unwell, seek medical advice immediately (show the label where possible)

\* Hazard labeling not necessary if quantity per bottle below 25 g or mL (certificate of exemption according to 67/548/EEC Art. 25, 1999/45/EC Art. 12 and German GefStoffV § 20 (3) and TRGS 200 7.1). For further information see Material Safety Data Sheet.

## 5 Protocols

### 5.1 Microplate assay procedure

#### Before starting the preparation:

- Check if the BSA reference protein stock solution was prepared according to section 3.
- Make sure that there are no precipitates in Protein Solving Buffer PSB and in the reference protein (BSA) solution (if necessary, heat to approx. 30°C).

#### 1 Prepare a BSA (reference protein) dilution series

Number seven reaction tubes according to column A (see table below; #1: BSA stock solution).

Add **50 µL Protein Solving Buffer PSB** to tubes #2 – #7 (column B).

Add **BSA solution** to tubes #2 – #6 according to column C.

The resulting protein concentration and amount are shown in columns D and E.

PSB contains detergent! When pipetting BSA and PSB solutions avoid bubble formation and foaming as far as possible.

A	B	C	D	E
Tube	Add PSB to tube	Add BSA solution to tube	Resulting BSA concentration	Resulting BSA in 20 µL
#1	BSA stock solution		1 µg/µL	20 µg
#2	50 µL	50 µL from tube #1	0.5 µg/µL	10 µg
#3	50 µL	50 µL from tube #2	0.25 µg/µL	5 µg
#4	50 µL	50 µL from tube #3	0.125 µg/µL	2.5 µg
#5	50 µL	50 µL from tube #4	0.063 µg/µL	1.25 µg
#6	50 µL	50 µL from tube #5	0.031 µg/µL	0.625 µg
#7	50 µL	–	0 µg/µL	0 µg

The prepared BSA dilutions series is sufficient for the determination of **two** calibration curves. Freeze BSA stock solution for storage. Keep dilutions series at room temperature during use and dispose all dilutions at the end of a working day.

**2 Dispense dilution series into microplate**

Add **20 µL** of **each dilution series solution** (#1 – #7) into microplate wells.

**20 µL of dilution series**

(#1: BSA stock solution; #2 – #6: BSA dilutions; #7: BSA-free PSB)

---

**3 Dispense your protein samples**

Pipette **20 µL** of your **samples** to empty wells.

**20 µL of samples**

*Alternatively, 1 – 60 µL of sample can be applied.*

---

**4 Fill up dilution series and protein samples**

Add **40 µL PSB** to each well (dilution series and protein samples). Final volume is 60 µL.

**+ 40 µL PSB**

*Alternatively, when applying other sample volumes than 20 µL in step 3, fill up with PSB to a final volume of 60 µL (e.g., 10 µL sample + 50 µL PSB).*

---

**5 Add Quantification Reagent QR**

Add **40 µL Quantification Reagent QR** to each well (dilution series and protein samples).

**+ 40 µL QR**

**Shake microplate** until a complete color change from blue to yellow occurs.

**Shake microplate**

*Caution: Quantification Reagent QR contains hydrochloric acid. Wear protective clothing and goggles.*

---

**6 Incubate**

Incubate microplate for **30 min** at room temperature.

Gently shake microplate after incubation, but avoid bubble formation and foaming. For optimal measurement the solution surface in the microplate well should be free of bubbles and foam. Light scattering caused by foam has impact on the measurement.

**Incubate 30 min**

*A variation in incubation time may result in reduced signal and loss of sensitivity. An incubation time of 30 ± 5 min is recommended.*

---



---

**7 Measure light extinction**

Measure light extinction photometrically at **570 nm**.

*Light extinction can be measured in the range of 530 – 700 nm. Typical correlation coefficients (concentration of BSA versus extinction value) of 0.97 – 1.00 are obtained within this wavelength range.*

**Measure  
extinction  
at 570 nm**

---

**8 Calculate protein concentration**

Calculate protein concentration of samples in relation to the BSA dilution series.

*Make sure that the protein concentration of your sample lies within the range of the largest (#1) and the smallest (#6) concentration of the calibration curve in order to obtain valid measurements. Do not extrapolate beyond this range.*

**Calculate  
protein  
concentration**

## 5.2 Semi-micro cuvette assay procedure

### Before starting the preparation:

- Check if the BSA reference protein stock solution was prepared according to section 3.
- Make sure that there are no precipitates in Protein Solving Buffer PSB and in the reference protein (BSA) solution (if necessary, heat to approx. 30°C).

### 1 Prepare a BSA (reference protein) dilution series

Number seven reaction tubes according to column A (see table below; #1: BSA stock solution).

Add **250 µL Protein Solving Buffer PSB** to tubes #2 – #7 (column B).

Add **BSA solution** to tubes #2 – #6 according to column C.

The resulting protein concentration and amount are shown in columns D and E.

A	B	C	D	E
Tube	Add PSB to tube	Add BSA solution to tube	Resulting BSA concentration	Resulting BSA in 20 µL
#1	BSA stock solution		1 µg/µL	200 µg
#2	250 µL	250 µL from tube #1	0.5 µg/µL	100 µg
#3	250 µL	250 µL from tube #2	0.25 µg/µL	50 µg
#4	250 µL	250 µL from tube #3	0.125 µg/µL	25 µg
#5	250 µL	250 µL from tube #4	0.063 µg/µL	12.5 µg
#6	250 µL	250 µL from tube #5	0.031 µg/µL	6.25 µg
#7	250 µL	–	0 µg/µL	0 µg

The prepared BSA dilutions series is sufficient for the determination of **one** calibration curve. Freeze BSA stock solution for storage. Keep dilutions series at room temperature during use and dispose all dilutions at the end of a working day.

---

## 2 Dispense dilution series into microcentrifuge tubes

Pipette **200 µL** of each **dilution series solution** (#1 – #7) into 1.5 mL microcentrifuge tubes (not supplied).

**200 µL of  
dilution series**

(#1: *BSA stock solution*; #2 – #6: *BSA dilutions*; #7: *BSA-free PSB*)

---

## 3 Dispense your protein samples

Pipette **200 µL** of your **samples** to (new) microcentrifuge tubes.

**200 µL of  
samples**

*Alternatively, 10 – 600 µL of sample can be applied.*

---

## 4 Fill up dilution series and protein samples

Add **400 µL PSB** to each microcentrifuge tube (dilution series and protein samples). Final volume is 600 µL.

**+ 400 µL PSB**

*Alternatively, when applying other sample volumes than 200 µL in step 3, fill up with PSB to a final volume of 600 µL (e.g., 100 µL sample + 500 µL PSB).*

---

## 5 Add Quantification Reagent QR

Add **400 µL Quantification Reagent QR** to each microcentrifuge tube (dilution series and protein samples).

**+ 400 µL QR**

**Shake tubes** until a complete color change from blue to yellow occurs.

**Shake tubes**

*Caution: Quantification Reagent QR contains hydrochloric acid. Wear protective clothing and goggles.*

---

## 6 Incubate

Incubate microcentrifuge tubes for **30 min** at room temperature.

Shake tubes after incubation. **Do not centrifuge** tubes at this point.

**Incubate  
30 min**

*A variation in incubation time may result in reduced signal and loss of sensitivity. An incubation time of 30 ± 5 min is recommended.*

---

---

## 7 Measure light extinction

Transfer the solution of each tube to a suitable semi-micro cuvette. Measure light extinction photometrically at **570 nm**.

*Light extinction can be measured in the range of 530 – 700 nm. Typical correlation coefficients (concentration of BSA versus extinction value) of 0.97 – 1.00 are obtained within this wavelength range.*

**Measure  
extinction  
at 570 nm**

---

## 8 Calculate protein concentration

Calculate protein concentration of samples in relation to the BSA dilution series.

*Make sure that the protein concentration of your sample lies within the range of the largest (#1) and the smallest (#6) concentration of the calibration curve in order to obtain valid measurements. Do not extrapolate beyond this range.*

**Calculate  
protein  
concentration**

### 5.3 Microcuvette assay procedure

#### Before starting the preparation:

- Check if the BSA reference protein stock solution was prepared according to section 3.
- Make sure that there are no precipitates in Protein Solving Buffer PSB and in the reference protein (BSA) solution (if necessary, heat to approx. 30°C).

#### 1 Prepare a BSA (reference protein) dilution series

Number seven reaction tubes according to column A (see table below; #1: BSA stock solution).

Add **50 µL Protein Solving Buffer PSB** to tubes #2 – #7 (column B).

Add **BSA solution** to tubes #2 – #6 according to column C.

The resulting protein concentration and amount are shown in columns D and E.

A	B	C	D	E
Tube	Add PSB to tube	Add BSA solution to tube	Resulting BSA concentration	Resulting BSA in 20 µL
#1	BSA stock solution		1 µg/µL	20 µg
#2	50 µL	50 µL from tube #1	0.5 µg/µL	10 µg
#3	50 µL	50 µL from tube #2	0.25 µg/µL	5 µg
#4	50 µL	50 µL from tube #3	0.125 µg/µL	2.5 µg
#5	50 µL	50 µL from tube #4	0.063 µg/µL	1.25 µg
#6	50 µL	50 µL from tube #5	0.031 µg/µL	0.625 µg
#7	50 µL	–	0 µg/µL	0 µg

The prepared BSA dilutions series is sufficient for the determination of **one** calibration curve. Freeze BSA stock solution for storage. Keep dilutions series at room temperature during use and dispose all dilutions at the end of a working day.

**2 Dispense dilution series into microcentrifuge tubes**

Pipette **40  $\mu$ L** of **each dilution series solution** (#1 – #7) into 1.5 mL microcentrifuge tubes (not supplied).

**40  $\mu$ L of  
dilution series**

*(#1: BSA stock solution; #2 – #6: BSA dilutions; #7: BSA-free PSB)*

---

**3 Dispense your protein samples**

Pipette **40  $\mu$ L** of your **samples** to (new) microcentrifuge tubes.

**40  $\mu$ L of  
samples**

*Alternatively, 1 – 120  $\mu$ L of sample can be applied.*

---

**4 Fill up dilution series and protein samples**

Add **80  $\mu$ L PSB** to each well (dilution series and protein samples). Final volume is 120  $\mu$ L.

**+ 80  $\mu$ L PSB**

*Alternatively, when applying other sample volumes than 40  $\mu$ L in step 3, fill up with PSB to a final volume of 120  $\mu$ L (e.g., 10  $\mu$ L sample + 110  $\mu$ L PSB).*

---

**5 Add Quantification Reagent QR**

Pipette **80  $\mu$ L Quantification Reagent QR** to each tube (dilution series and protein samples).

**+ 80  $\mu$ L QR**

**Shake tube** until a complete color change from blue to yellow occurs.

**Shake tube**

*Caution: Quantification Reagent QR contains hydrochloric acid. Wear protective clothing and goggles.*

---

**6 Incubate**

Incubate tubes for **30 min** at room temperature.

**Incubate  
30 min**

Shake tubes after incubation. **Do not centrifuge** tubes at this point!

*A variation in incubation time may result in reduced signal and loss of sensitivity. An incubation time of 30  $\pm$  5 min is recommended.*

---

---

## 7 Measure light extinction

Transfer the solution of each tube to a suitable microcuvette. Measure light extinction photometrically at **570 nm**.

*Light extinction can be measured in the range of 530 – 700 nm. Typical correlation coefficients (concentration of BSA versus extinction value) of 0.97 – 1.00 are obtained within this wavelength range.*

**Measure  
extinction  
at 570 nm**

---

## 8 Calculate protein concentration

Calculate protein concentration of samples in relation to the BSA dilution series.

*Make sure that the protein concentration of your sample lies within the range of the largest (#1) and the smallest (#6) concentration of the calibration curve in order to obtain valid measurements. Do not extrapolate beyond this range.*

**Calculate  
protein  
concentration**

## 5.4 Low volume assay procedure

### Before starting the preparation:

- Check if the BSA reference protein stock solution was prepared according to section 3.
- Make sure that there are no precipitates in Protein Solving Buffer PSB and in the reference protein (BSA) solution (if necessary, heat to approx. 30°C).

### 1 Prepare a BSA (reference protein) dilution series

Number six reaction tubes according to column A (see table below; #1: BSA stock solution).

Add **20 µL Protein Solving Buffer PSB** to tubes #2 – #6 (column B).

Add **BSA solution** to tubes #2 – #5 according to column C.

The resulting protein concentration and amount are shown in columns D and E.

A	B	C	D	E
Tube	Add PSB to tube	Add BSA solution to tube	Resulting BSA concentration	Resulting BSA in 20 µL
#1	BSA stock solution		1 µg/µL	7.5 µg
#2	20 µL	20 µL from tube #1	0.5 µg/µL	3.75 µg
#3	20 µL	20 µL from tube #2	0.25 µg/µL	1.88 µg
#4	20 µL	20 µL from tube #3	0.125 µg/µL	0.94 µg
#5	20 µL	20 µL from tube #4	0.063 µg/µL	0.47 µg
#6	20 µL	–	0 µg/µL	0 µg

The prepared BSA dilutions series is sufficient for the determination of **two** calibration curves. Freeze BSA stock solution for storage. Keep dilutions series at room temperature during use and dispose all dilutions at the end of a working day.



---

**2 Dispense dilution series into microcentrifuge tubes**

Pipette **7.5 µL** of **each dilution series solution** (#1 – #6) into 1.5 mL microcentrifuge tubes (not supplied).

**7.5 µL of dilution series**

(#1: BSA stock solution; #2 – #5: BSA dilutions; #6: BSA-free PSB)

---

**3 Dispense your protein samples**

Pipette **7.5 µL** of your **samples** to (new) microcentrifuge tubes.

---

**7.5 µL of samples**

**4 Fill up dilution series and protein samples**

*Not necessary! Proceed directly with step 5.*

---

**5 Add Quantification Reagent QR**

Add **5 µL Quantification Reagent QR** to each tube (dilution series and protein samples).

**+ 5 µL QR**

**Mix (e.g., by pipetting up and down)** until a complete color change from blue to yellow occurs.

**Mix**

*Caution: Quantification Reagent QR contains hydrochloric acid. Wear protective clothing and goggles.*

---

**6 Incubate**

Incubate tubes for **30 min** at room temperature.

**Incubate 30 min**

Shake tubes after incubation. **Do not centrifuge** at this point!

*A variation in incubation time may result in reduced signal and loss of sensitivity. An incubation time of  $30 \pm 5$  min is recommended.*

---

## 7 Measure light extinction

Transfer 10  $\mu\text{L}$  of the solution of each tube to a suitable **low volume photometer with 1 mm path length**. Measure light extinction photometrically at **570 nm**. Avoid bubbles in the solution because they severely disturb the measurement.

*Caution: The solution to be measured contains HCl; check the compatibility of your instrument with HCl. Do not spill. Immediately remove solution from the photometer after measurement.*

*Light extinction can be measured in the range of 530 – 700 nm. Typical correlation coefficients (concentration of BSA versus extinction value) of 0.97 – 1.00 are obtained within this wavelength range.*

**Measure  
extinction  
at 570 nm**

## 8 Calculate protein concentration

Calculate protein concentration of samples in relation to the BSA dilution series.

*Make sure that the protein concentration of your sample lies within the range of the largest (#1) and the smallest (#5) concentration of the calibration curve in order to obtain valid measurements. Do not extrapolate beyond this range.*

**Calculate  
protein  
concentration**

## 6 Appendix

### 6.1 Troubleshooting

Problem	Possible cause and suggestions
Lowest value of calibration curve cannot be measured	<p data-bbox="346 344 598 367"><i>Storage of dilution series</i></p> <ul data-bbox="346 379 978 427" style="list-style-type: none"> <li>• Do not store dilution series of the BSA reference protein. Prepare fresh dilution series.</li> </ul> <p data-bbox="346 472 710 494"><i>Photometer, microplates or cuvettes</i></p> <ul data-bbox="346 507 978 608" style="list-style-type: none"> <li>• Sensitivity of the assay may be influenced by the type of photometer, microplates, or cuvettes used. If the lowest calibration point is not discriminated against background, prepare a calibration series with higher BSA amounts.</li> </ul>
Samples appear turbid after addition of Quantification Reagent QR	<p data-bbox="346 678 611 700"><i>High protein concentration</i></p> <ul data-bbox="346 713 978 762" style="list-style-type: none"> <li>• As long as the measured extinction of your sample falls within the range of the calibration curve, this is acceptable.</li> </ul>
Varying results upon multiple measurements	<p data-bbox="346 831 978 879"><i>Samples not mixed immediately before extinction measurement</i></p> <ul data-bbox="346 892 978 1241" style="list-style-type: none"> <li>• Shake microplate immediately before extinction measurement.</li> <li>• Shake reaction tubes after incubation and before transfer to semi-micro cuvettes. After transfer of samples to semi-micro cuvettes, measure extinction immediately.</li> <li>• Strictly keep to the recommended incubation time.</li> <li>• Do not centrifuge at any time after addition of Quantification Reagent QR.</li> <li>• Avoid bubble formation and foaming, especially for protocol section 5.1. (microplate assay procedure). Light scattering caused by foam has impact on turbidity measurements.</li> </ul>
Protein Solving Buffer PSB appears turbid	<p data-bbox="346 1295 598 1318"><i>Low storage temperature</i></p> <ul data-bbox="346 1331 665 1353" style="list-style-type: none"> <li>• Warm PSB to approx. 30°C.</li> </ul>

Similar extinction for all dilution series samples

*Fill-level of semi-micro or microcuvette not compatible with photometer*

- Make sure that the sample volume in the semi-micro cuvette is high enough to let the light beam pass through the solution. Consult your photometer user manual. Check the compatibility of disposable cuvettes used with your photometer – consider light beam center height and cuvette fill volume.
- 

## 6.2 Ordering information

Product	REF	Number of assays or preparations
Protein Quantification Assay	740967.50/.250	50/250
NucleoSpin® RNA/Protein	740933.10/50/.250	10/50/250
NucleoSpin® TriPrep*	740966.10/50/.250	10/50/250
Porablot transfer membranes	see <a href="http://www.mn-net.com/bioanalysis">www.mn-net.com/bioanalysis</a>	
Blotting paper	see <a href="http://www.mn-net.com/bioanalysis">www.mn-net.com/bioanalysis</a>	

---

\* DISTRIBUTION AND USE OF NUCLEOSPIN® TRIPREP IN THE USA IS PROHIBITED FOR PATENT REASONS.

## 6.3 References

**Bradford MM** (1976): A rapid and sensitive method for the quantitation of microgram quantities of protein utilizing the principle of protein-dye binding. *Anal. Biochem.* 72, 248-254.

**Karlsson JO *et al.*** (1994): A method for protein assay in Laemmli buffer. *Analytical Biochemistry* 219, 144-146.

**Laemmli UK** (1970): Cleavage of structural proteins during the assembly of the head of bacteriophage T4. *Nature* 227, 680-685 (1970).

**Lowry OH *et al.*** (1951): Protein measurement with the Folin phenol reagent. *J. Biol. Chem.* 193, 265-275.

**Smith PK *et al.*** (1985): Measurement of protein using bicinchoninic acid. *Anal. Biochem.* 150(1), 76-85.

## 6.4 Product use restriction/warranty

**Protein Quantification Assay** kit components were developed, designed, distributed, and sold **FOR RESEARCH PURPOSES ONLY** They are suitable **FOR IN-VITRO USES ONLY**. No claim or representation is intended for its use to identify any specific organism or for clinical use (diagnostic, prognostic, therapeutic, or blood banking).

It is rather the responsibility of the user to verify the use of the **Protein Quantification Assay** kit for a specific application range as the performance characteristic of this kit has not been verified to a specific organism.

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 an extra copy.

MACHEREY-NAGEL does not warrant against 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; against defects in products or components not manufactured by MACHEREY-NAGEL, or against 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.

Please contact:  
MACHEREY-NAGEL Germany  
Tel.: +49 (0) 24 21 969 270  
e-mail: TECH-BIO@mn-net.com

Last updated: 12/2006, Rev.02

---

**Trademarks:**

NanoPhotometer is a trademark of IMPLEN Ltd.  
NucleoSpin is a registered trademark of MACHEREY-NAGEL GmbH & Co KG

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.