

# Overall Exosome Capture and Quantification Assay Kit ab287851

4 Images

Overview

Product name	Overall Exosome Capture and Quantification Assay Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 10%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			CV%
				< 13%
Sample type	Cell culture supernatant			
Assay type	Sandwich			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	<p>The Overall Exosome Capture and Quantification Assay Kit ab287851 (Previously known as ExoQuant Overall Exosome Capture and Quantification AssayKit (Cell Media, Colorimetric) Kit), is a double sandwich enzyme-linked immunoassay for quantitative and qualitative analysis of exosomes. It is a successful platform for exosomes quantification and characterization from small amount of human biological fluids (plasma, serum, urine, saliva) or cell media. It may be exploited to identify exosomes released by cancer cells in the plasma and urine of tumor patients in various disease conditions. The kit consists of ELISA plates pre-coated with proprietary pan-exosome antibodies enabling specific capture of exosomes from cell culture supernatants. Quantification and characterization of exosomal proteins is subsequently performed using appropriate detection antibodies against exosome associated antigens that can be for either generic or cell/tissue-specific exosomes. The kit contains Lyophilized Exosome Standards for assay calibration. Lyophilized Exosome Standards, characterized for protein content and particle number allow the quantification of unknown sample by a standard calibration curve. Transparent and white plates are available depending on the downstream detection approach (colorimetric and luminometric respectively)</p>			

**Applications:** Overall exosome immunocapture and quantification from cell culture supernatant

**Features & Benefits:** Double sandwich enzyme-linked immunoassay for quantitative and qualitative analysis of exosomes

**Platform**

Microplate (12 x 8 well strips)

**Properties**

**Storage instructions**

Store at +4°C. Please refer to protocols.

Components	96 tests
1X Sample buffer	2 x 10ml
25X Washing buffer	1 x 25ml
Exosome Standards	2 x 100µg
HRP-Conjugate	1 x 5µl
Immunoplate (Transparent)	1 unit
Primary Antibody	1 x 20µl
Stop Solution	1 x 10ml
Substrate Chromogenic Solutions	1 x 10ml

**Function**

Involved in platelet activation and aggregation. Regulates paranodal junction formation. Involved in cell adhesion, cell motility and tumor metastasis. Required for sperm-egg fusion.

**Tissue specificity**

Expressed by a variety of hematopoietic and epithelial cells.

**Sequence similarities**

Belongs to the tetraspanin (TM4SF) family.

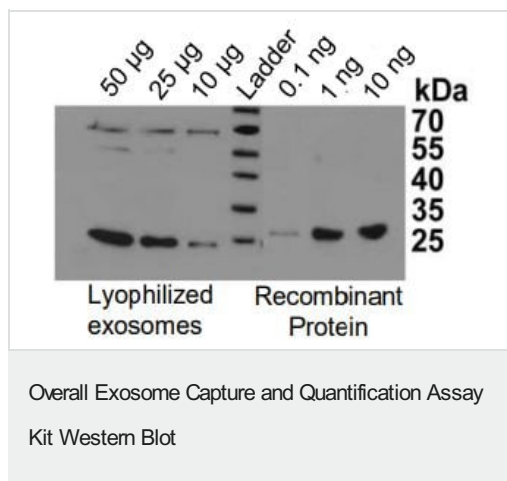
**Post-translational modifications**

Protein exists in three forms with molecular masses between 22 and 27 kDa, and is known to carry covalently linked fatty acids.

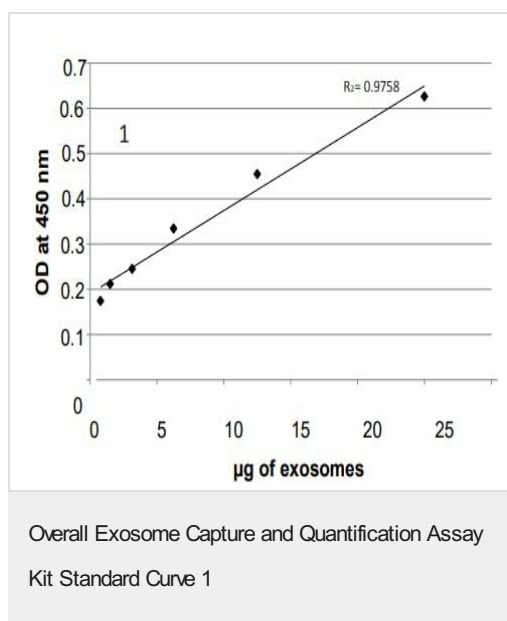
**Cellular localization**

Membrane.

**Images**

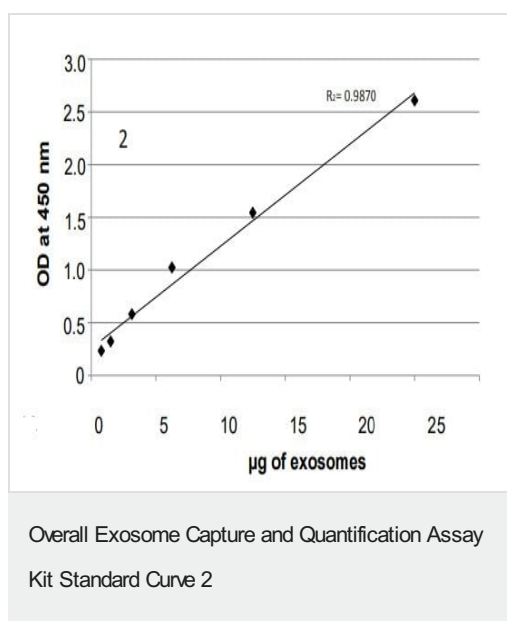


CD9 marker detection by Western Blotting on lyophilized exosomes from human plasma and the recombinant CD9 protein.



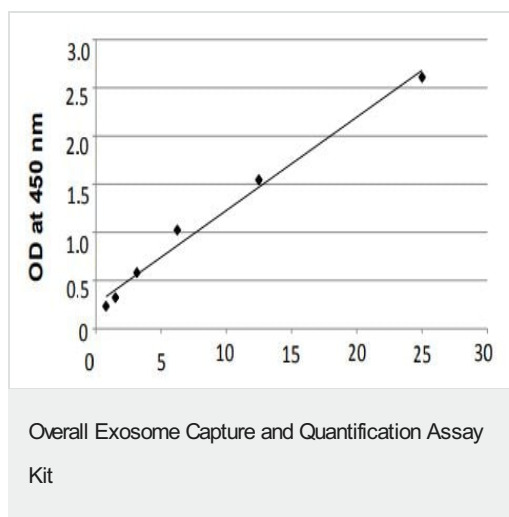
Example of standard curves obtained. Exosome standard preparation derived from COLO1 Cell Culture Supernatant

This standard curve is for demonstration only. For quantification purposes, a standard curve must be obtained with every assay.



Example of standard curves obtained. Exosome standard preparation derived from Plasma of healthy donors.

This standard curve is for demonstration only. For quantification purposes, a standard curve must be obtained with every assay.



CD9 titration of plasma healthy donor exosome standards.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

#### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors