

Product datasheet

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 | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 10% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| Overall | | | | < 13% |

Sample type Cell culture supernatant

Assay type Sandwich

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

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)

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.

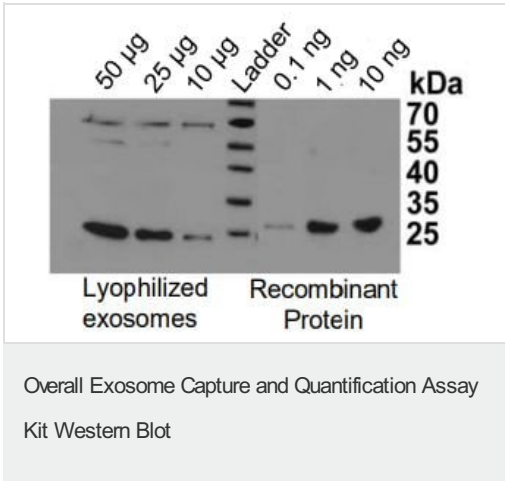
Post-translational modifications

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

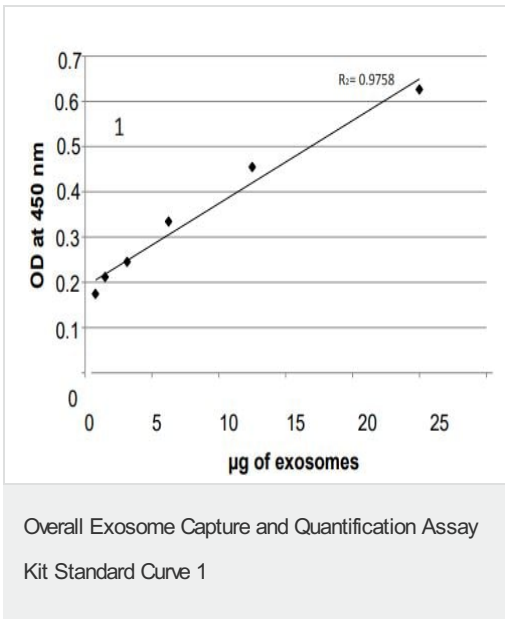
Cellular localization

Membrane.

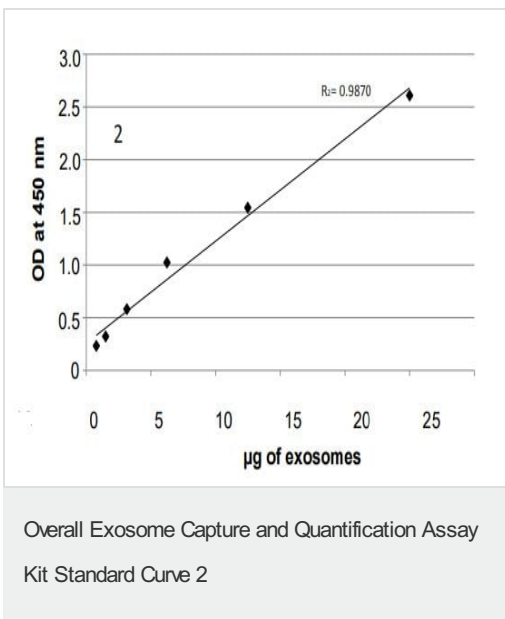
Images



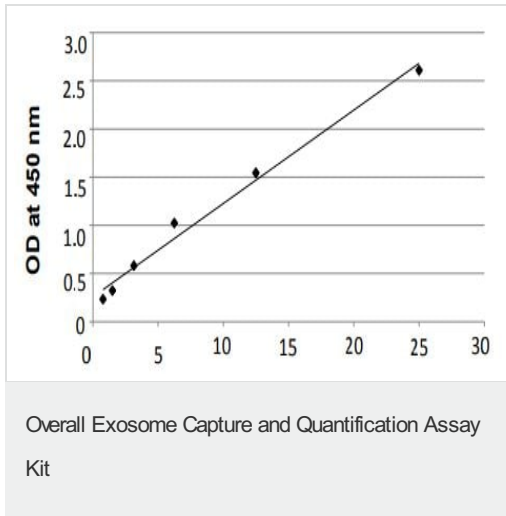
CD9 marker detection by Western Blotting on lyophilized exosomes from human plasma and the recombinant CD9 protein.



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