

Product datasheet

Lyophilized MCF7 Exosome Standards ab239691

★★★★★ [2 Abreviews](#) [1 Image](#)

Overview

Product name	Lyophilized MCF7 Exosome Standards
Species reactivity	Reacts with: Human
Product overview	Lyophilized MCF7 Exosome Standards (ab239691) are lyophilized exosomes (1×10^{12}) derived from human cancer cell line. MCF7 is a human breast cancer cell line.

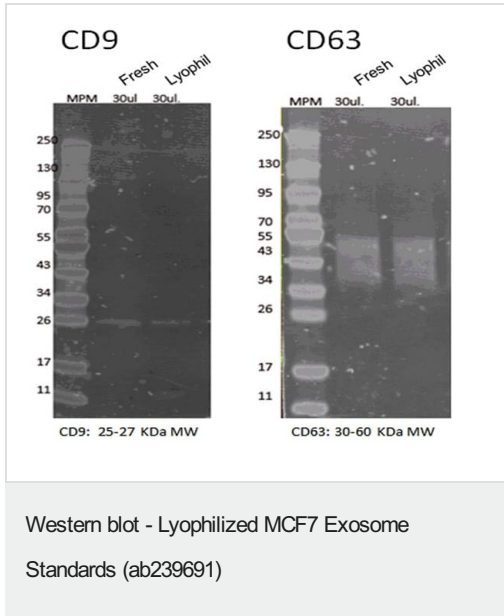
Properties

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

Components	100 µg
Lyophilized MCF7 Exosome Standards	1 x 100µg

Relevance Exosomes are best defined as extracellular vesicles that are released from cells upon fusion of an intermediate endocytic compartment, the multivesicular body (MVB), with the plasma membrane. This liberates intraluminal vesicles (ILVs) into the extracellular milieu and the vesicles thereby released are what we know as exosomes. They contain abundant proteins, DNA, mRNA, and non-coding RNAs. They are now considered as an additional mechanism for intercellular communication, allowing cells to exchange proteins, lipids and genetic material. Increasing studies have shown that exosomes play an important role in tumour initiation, growth, progression, metastasis, drug resistance and immune escape.

Images



Fresh and lyophilized Exosome batches were analyzed and compared by WB in native conditions for exosomal markers, by anti-CD9 (Clone VJ1/20) and anti-CD63 (Clon TEA3/18) antibodies at a 1:1000 dilution (0.1 mg/ml).

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