

Product datasheet

Plasma Membrane Isolation Kit ab284937

[2 Images](#)

Overview

Product name	Plasma Membrane Isolation Kit
Sample type	Cell culture media, Tissue Culture Media
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mammals
Product overview	<p>This Plasma Membrane Isolation Kit (ab284937) (previously known as OrgFrontier Plasma Membrane Isolation Kit K414) provides a proprietary set of reagents and buffers designed to enable optimized extraction and isolation of intact plasma membrane from mammalian tissues and cells.</p>

The protocol requires an ultracentrifuge to be performed and can be completed in a few hours.

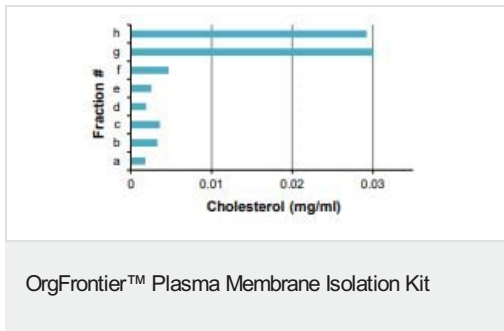
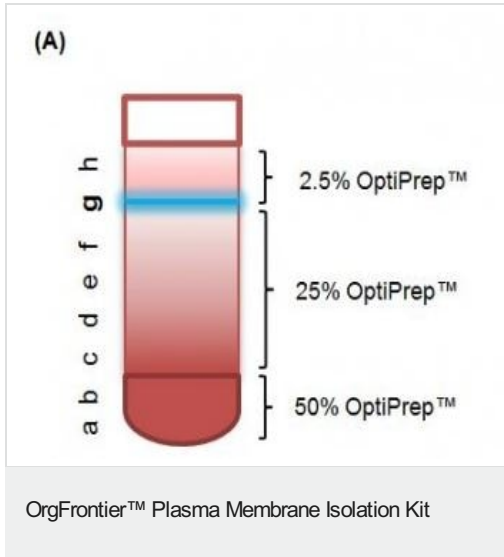
Notes This product is manufactured by BioVision, an Abcam company and was previously called K414 OrgFrontier™ Plasma Membrane Isolation Kit. K414-10 is the same size as the 10 test size of ab284937.

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	10 tests
Density Gradient Media II	1 x 85ml
Gradient Dilution Buffer II	1 x 100ml
Homogenization Buffer	1 x 50ml
Protease Inhibitor Cocktail II	1 x 1ml

Images



Cholesterol content by fraction, assayed using Cholesterol/Cholesteryl Ester Assay Kit - Quantitation (**ab65359**). Cholesterol is highly enriched in the plasma membrane (approximately 22% of the total lipid content of the plasma membrane), but is found in much lower concentration (2%-6% of total lipids) in other organelle membranes.

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