

Product datasheet

Mitochondrial Protein IP Kit ab239710

Overview

Product name Mitochondrial Protein IP Kit

Product overview Mitochondrial Protein IP Kit (ab239710) is optimized for immunoprecipitation (IP and co-IP) using mitochondria and mitochondrial extracts. The buffer is a gentle formulation, which maintains the stability of mitochondrial complexes. The Mitochondrial Protein IP kit is provided with different choices of detergents like n-Dodecyl-beta-D-maltoside, Triton X-100 and digitonin to achieve different stringency conditions for protein-protein interaction studies. Triton X-100 is the most commonly used detergent especially for membrane protein solubilization. However, in case of fragile complexes digitonin or n-Dodecyl-beta-D-maltoside is the choice of detergents.

Notes Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Tested applications **Suitable for:** IP

Properties

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

Components	50 tests
10% Digitonin	1 x 1ml
10% n-Dodecyl-beta-D-maltoside	1 x 1ml
10% Triton X-100	1 x 1ml
Mitochondria Protein IP Buffer	1 x 50ml
Protease Inhibitor Cocktail	1 vial
Wash Buffer (5X)	1 x 50ml

Relevance Mitochondria are the power house of the cell. They are distinct organelles with two membranes. Usually they are rod shaped, however they can be round. The outer membrane limits the organelle and the inner membrane is thrown into folds or shelves that project inward and are called "cristae"

mitochondriales".

Cellular localization

Mitochondrial

Applications

Our [Abpromise guarantee](#) covers the use of **ab239710** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
IP		Use at an assay dependent concentration.

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