

Mitochondria/Cytosol Fractionation Kit ab65320

96 References

Overview

Product name	Mitochondria/Cytosol Fractionation Kit
Sample type	Adherent cells, Suspension cells
Assay type	Direct
Assay time	4h 00m
Product overview	<p>Abcam's Mitochondria/Cytosol Fractionation Kit provides unique formulations of reagents for effective isolation of a highly enriched mitochondrial fraction from cytosolic fraction of mammalian cells including both apoptotic and nonapoptotic cells. The enriched mitochondrial and cytosolic fractions can be used for studying apoptotic and signal transduction pathways to detect translocation of factors interested between the two fractions by Western blotting, ELISA, or other assays. Procedures are simple and easy to perform, no ultracentrifugations and toxic chemicals are involved.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

Notes This product is manufactured by BioVision, an Abcam company and was previously called K256 Mitochondria/Cytosol Fractionation Kit. K256-100 is the same size as the 100 test size of ab65320.

Properties

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

Components	25 tests	100 tests	100 tests	25 tests
5X Cytosol Extraction Buffer I	1 x 5ml	1 x 20ml	1 x 20ml	1 x 5ml
DTT II	1 x 100µl	1 x 100µl	1 x 110µl	1 x 110µl
Mitochondria Extraction Buffer I	1 x 2.5ml	1 x 10ml	1 x 10ml	1 x 2.5ml
Protease Inhibitor Cocktail I	1 vial	1 vial	1 vial	1 vial

Relevance Mitochondria/Cytosol Fractionation is the process for effective isolation of a highly enriched mitochondrial fraction from the cytosolic fraction of mammalian cells.

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