


Product datasheet

Yeast Mitochondria Isolation Kit ab178779

[4 References](#) [1 Image](#)

Overview

Product name	Yeast Mitochondria Isolation Kit
Sample type	Suspension cells
Assay type	Direct
Species reactivity	Reacts with: <i>Saccharomyces cerevisiae</i> , <i>Pichia pastoris</i> Predicted to work with: <i>Schizosaccharomyces pombe</i> , <i>Candida albicans</i> 
Product overview	Isolated mitochondria are useful for studying mitochondrial respiration, assembly of respiratory complexes, apoptosis, mtDNA and mtRNA, and for protein profiling.

Abcam's Yeast Mitochondria Isolation Kit (ab178779) has been tested for *Pichia pastoris* and *Saccharomyces cerevisiae* and can be used to isolate mitochondria from other yeast.

Under fermentable media, the yield is ~ 150 - 200 µg of mitochondria, and under non-fermentable media (such as glycerol), # 200 - 250 µg of mitochondria from a culture growing at OD = 20.

ab178779 is only sufficient for 25 samples.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K259 Yeast Mitochondria Isolation Kit. K259-50 is the same size as the 50 test size of ab178779.

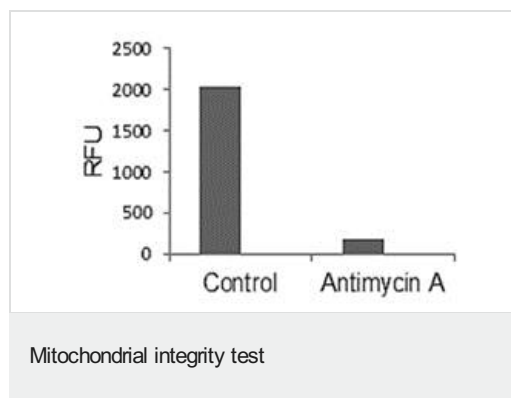
Properties

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

Components	50 tests
Buffer A	1 x 50ml
Buffer B	1 x 50ml
Dilution Buffer I	1 x 10ml
DTT IV	1 x 1ml

Components	50 tests
Lysis Enzyme Mix	1 x 200µl
Protease Inhibitor Cocktail I	1 vial
Yeast Homogenization Buffer	1 x 50ml

Images



Purified mitochondria were analyzed for intactness by using JC-1 dye, which tests the electrochemical proton gradient ($\Delta\Psi$) of the inner mitochondrial membrane. The intact purified mitochondria (control) show aggregation of JC-1 dye, whose signal can be measured at Ex/Em = 530/590 nm. Treatment with antimycin C (100 μ M), an inhibitor of the electron transport chain, dissipates the mitochondrial membrane potential resulting in reduced fluorescence signal.

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