


Product datasheet

Yeast Mitochondria Isolation Kit ab178779

1 References 1 Image

Overview

Product name	Yeast Mitochondria Isolation Kit
Sample type	Suspension cells
Assay type	Direct
Species reactivity	<p>Reacts with: <i>Saccharomyces cerevisiae</i>, <i>Pichia pastoris</i></p> <p>Predicted to work with: <i>Schizosaccharomyces pombe</i>, <i>Candida albicans</i> </p>
Product overview	<p>Isolated mitochondria are useful for studying mitochondrial respiration, assembly of respiratory complexes, apoptosis, mtDNA and mtRNA, and for protein profiling.</p> <p>Abcam's Yeast Mitochondria Isolation Kit (ab178779) has been tested for <i>Pichia pastoris</i> and <i>Saccharomyces cerevisiae</i> and can be used to isolate mitochondria from other yeast.</p> <p>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.</p> <p>ab178779 is only sufficient for 25 samples.</p>

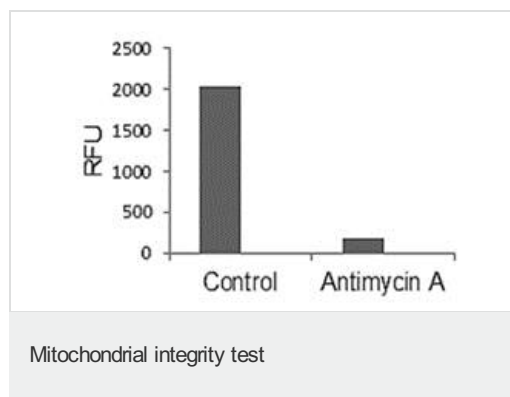
Properties

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

Components	Identifier	50 tests
Buffer A	Blue	1 x 50ml
Buffer B		1 x 50ml
DTT		1 x 1ml
Homogenization Buffer	Amber	1 x 50ml
Lysis Enzyme Mix	Orange	1 x 200µl
Protease Inhibitor Cocktail	Red	1 vial

Components	Identifier	50 tests
Storage Buffer		1 x 10ml

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