

Product datasheet

Mitochondria Isolation Kit For Tissue and Cultured Cells ab288084

[1 Image](#)

Overview

Product name	Mitochondria Isolation Kit For Tissue and Cultured Cells
Sample type	Adherent cells, Suspension cells, Tissue Homogenate
Product overview	<p>Mitochondria Isolation Kit For Tissue and Cultured Cells (ab288084), previously known as Mammalian Mitochondria Isolation Kit for Tissue and Cultured Cells (Cat# K288-50), offers two options for the simple and rapid isolation of intact mitochondria.</p> <p>The first option is utilizing a reagent-based method allowing parallel processing of multiple samples.</p> <p>The second option uses traditional dounce homogenization, which provides better mitochondrial yield.</p>

Properties

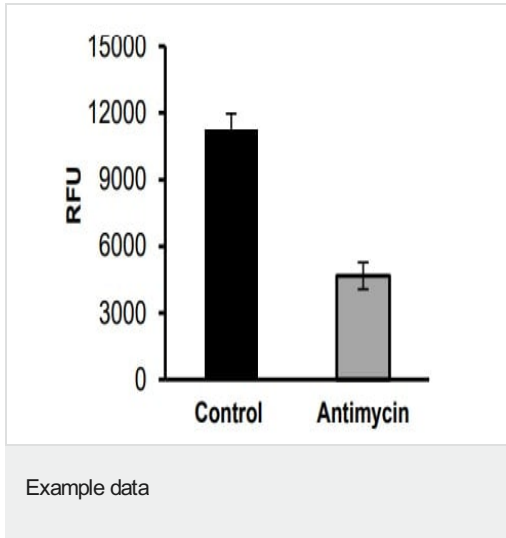
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	50 tests
Mitochondria Isolation Buffer	NM	2 x 100ml
Reagent A	Blue cap	1 x 1ml
Storage Buffer	NM	1 x 25ml

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

Images



Mitochondria Integrity Test:

Mitochondria (100 μg) were purified from the mouse liver. Purified mitochondria were analyzed for intactness by using JC-1 dye, which tests the electrochemical proton gradient ($\Delta\Psi$) of the inner mitochondrial membrane. Purified mitochondria was thawed on ice and incubated with JC-1 dye (1 $\mu\text{g/ml}$) for 15 minutes at 37°C. The intact purified mitochondria show aggregation of JC-1 dye whose signal can be measured at Ex/Em = 530/590nm. Treatment with Antimycin A (100 μM) 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