

Mitochondria Isolation Kit for Cultured Cells ab110170

★★★★★ [2 Abreviews](#) [67 References](#) [1 图像](#)

概述

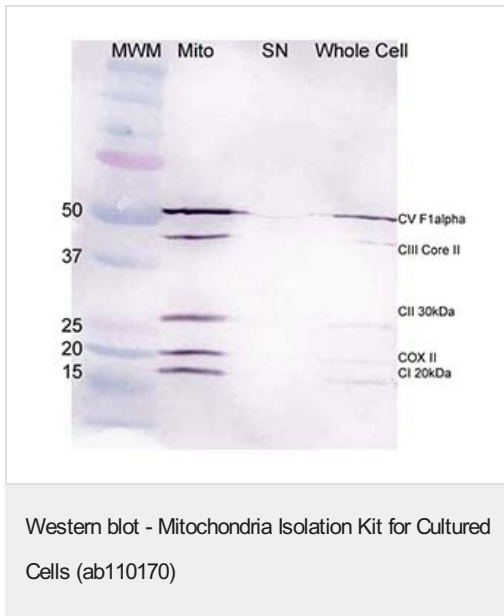
产品名称	Mitochondria分离试剂盒for Cultured Cells
检测时间	1h 00m
种属反应性	与反应: Mammals
产品概述	<p>ab110170 contains sufficient reagents for up to 20 benchtop mitochondria isolations from mammalian cultured cells. Isolated mitochondria can be analyzed by Western Blotting, Immunocapture, or Activity assay.</p> <p>Available with a dounce homogenizer as kit ab110171.</p> <p>Mitochondrial isolation protocol summary:</p> <ul style="list-style-type: none">- suspend cells in reagent A and incubate for 10 min- homogenize using dounce homogenizer- spin at 1,000 g for 10 min and retain supernatant, repeat- spin at 12,000 g for 15 min and retain pellet- resuspend pellet in reagent C and analyze or freeze

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 kit
Reagent A	1 x 50ml
Reagent B	1 x 50ml
Reagent C	1 x 10ml

图片



Isolated mitochondria show enriched signal when compared to the whole cell extract. In lane 1, MRC5 mitochondria isolated with Abcam's ab110170 (MS852) Mitochondria Isolation Kit for Cultured Cells were loaded at 20 µg. In lanes 2 and 3, post-spin supernatant (SN) and whole cell extract were loaded at 20 µg.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors