

Mitochondria/Cytosol fractionation kit

The core kit contains unique formulations of buffers and reagents that allow isolation of mitochondria and cytosol in only two easy steps. No ultracentrifugations are required. No toxic chemicals are involved.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-2760	1Kit
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Protect from light. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

For the effective isolation of a highly enriched mitochondrial fraction from the cytosolic fraction of various mammalian cells including both apoptotic and non-apoptotic cells. The enriched mitochondrial and cytosolic fractions can be used to study apoptosis and signal transduction pathways by detecting translocations of factors of interest between the two fractions by ELISA, Western blot or other assays.

Contents

10ml Mitochondria Extraction Buffer, 20ml 5X Cytosol Extraction Buffer, 100µl DTT (1M), 1 vial Protease Inhibitor Cocktail.

Quantity

~100 tests.

Species Reactivity

Mammalian

Last modified: February 17, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com