

Mitochondrial DNA Isolation Kit

To isolate intact mitochondrial DNA from a variety of cells and tissue types without contamination from genomic DNA.

Mitochondria are membrane-bound organelles found in all eukaryotic cells. They are often referred to as the powerhouses of the cell. Their primary function is to produce ATP necessary for various cellular processes, including metabolism, growth, and movement. Dysfunction in mitochondria is associated with mitochondrial disorders including muscle weakness, fatigue, metabolic strokes, seizures, cardiomyopathy, arrhythmias, developmental or cognitive disabilities, diabetes mellitus etc. The Mitochondrial DNA Extraction Kit provides a simple, fast and reliable tool for isolating intact mitochondrial DNA from a variety of cells and tissue types without contamination from genomic DNA. The recovered mitochondrial DNA is compatible with various downstream applications including PCR, cloning etc.

Ordering Information

[Order Online »](#)

AKR-PI-0105	50preparations
-------------	----------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Species Reactivity Species independent

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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