

Mitochondrial Fractionation Kit

reproducible mitochondrial and cytosolic fractionation for apoptosis research

Active Motif's Mitochondrial Fractionation Kit isolates highly enriched mitochondrial and cytosolic fractions from mammalian cell lines, simplifying the study of protein translocation events that occur during apoptosis and in many signal transduction pathways. The procedure is simple to perform and does not require ultra-centrifugation or use of any toxic chemicals. Isolated fractions can be used in Western blots, ELISAs and in other assays.

MITOCHONDRIAL EXTRACT KIT ADVANTAGES

- Eliminates cross-contamination of fractions
- Quality-controlled reagents ensure reproducibility
- No need to optimize your own reagents and protocol

The Mitochondrial Fractionation Kit's high-quality reagents and optimized protocol eliminate cross-contamination and produce high yields of properly segregated mitochondrial and cytosolic fractions (Figure 1). This ensures that you get accurate, reproducible results.

CONTENTS

Cytosol and Mitochondria Buffers, DTT and Protease Inhibitor. Sufficient reagents are included to perform 100 fractionations from a 100 mm dish.

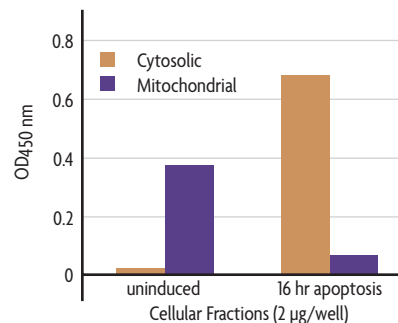


Figure 1: Location of cytochrome c in HeLa cells.

HeLa cells grown to 90-95% confluence were treated with 10 μ M Actinomycin D to induce apoptosis. Samples were harvested before as well as 16 hours after treatment. The Mitochondrial Fractionation Kit was used to isolate cytosolic and mitochondrial fractions. Two μ g of each fraction was assayed using FunctionELISA[™] Cytochrome c, which has since been discontinued, to monitor changes in mitochondrial and cytosolic cytochrome c levels.

Product	Format	Catalog No.
Mitochondrial Fractionation Kit	100 rxns	40015