

CytoFrac Cell Fractionation Kit

To provide reproducible separation of four distinct subcellular fractions: **Cytoplasmic, Nuclear, Subcellular Membrane/Particulate and Cytoskeletal fractions from a single**

mammalian sample. The CytoFrac Cell Fractionation system is designed to provide reproducible separation of four distinct subcellular fractions: Cytoplasmic, Nuclear, Subcellular Membrane/Particulate and Cytoskeletal fractions from a single mammalian sample. The procedure is fast and easy, requiring only 2 hours with no ultracentrifugation required. All four fractions obtained are suitable for diverse downstream applications such as 1 or 2D gel electrophoresis, enzyme activity assays, gel shift assay, and/or Western blotting.

Ordering Information

[Order Online »](#)

AKR-PI-0106	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