

EPIXTRACT® Nuclear Protein Isolation Kit

Rapid, reliable method to isolate nuclear proteins in mammalian cells and tissues.

EPIXTRACT® Nuclear Protein Isolation Kit II provides a simple and reliable method to isolate nuclear proteins. Extracts can be used for a variety of downstream applications including western blotting, protein-DNA binding assays, nuclear enzyme assays and others requiring nuclear proteins.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-45016-0100	100Reactions
----------------	--------------

Manuals, SDS & CofA

[View Online »](#)

- Rapid - extracts in 1 hour
- Simple - optimized protocol with easy steps
- Reliable - reproducible results
- High Yield - under optimal conditions, total yield up to 100µg

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For the extraction of nuclear proteins from mammalian cells or tissues.

Assay Time <1 hour

Contents Pre-extraction Buffer, Extraction Buffer, Protease Inhibitor Cocktail, DTT



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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