

EPIXTRACT[®] Total Histone Extraction Kit (High Throughput)

**Rapid, reliable, high throughput method to
isolate histone proteins in mammalian cells**

EPIXTRACT[®] Total Histone Extraction Kit (High Throughput) provides a simple and reliable method to isolate histone proteins in 1 hour. Extracts can be used for a variety of downstream applications such as post-translational modification assays.

Ordering Information

[Order Online »](#)

ENZ-45013-0002	2x96Reactions
----------------	---------------

Manuals, SDS & CofA

[View Online »](#)

- Rapid - extracts in 1 hour
- Simple - optimized protocol with easy steps
- Reliable - reproducible results that keep post-translational modifications intact
- High Efficiency - extract histones from as low as 3×10^4 cells

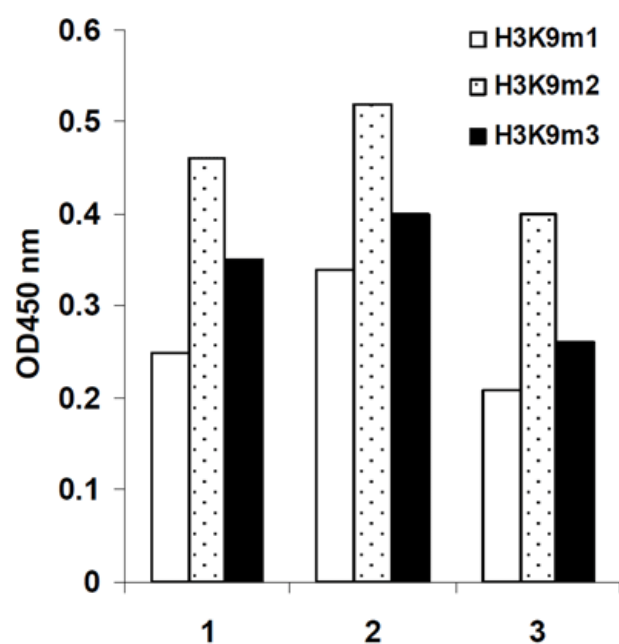


Figure 2. Histone extracts were prepared from MCF-7 cells (30,000 cells/well) using the EPIXTRACT® Total Histone Extraction Kit (High Throughput) and acetyl histone H3 was quantified.

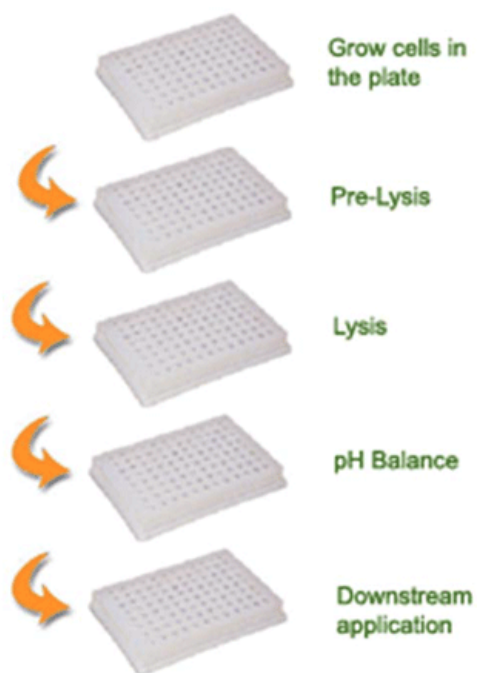


Figure 1. Schematic procedure for using the EPIXTRACT® Total Histone Extraction Kit (High Throughput)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

For the extraction of total histone proteins from mammalian cells in a 96 well plate format.

Assay Time

1 hour

Contents

Pre-Lysis Buffer, Lysis Buffer, Balance Buffer, DTT

Last modified: May 29, 2024



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)