

# EPIXTRACT<sup>®</sup> Total Histone Extraction Kit

**Rapid, reliable method to isolate histone proteins in mammalian cells and tissues for histone modification applications.**

EPIXTRACT<sup>®</sup> Total Histone Extraction Kit 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-45014-0100	100Reactions
----------------	--------------

## 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 10<sup>5</sup> cells or 1mg tissue



Fig 1. Schematic procedure for using the EPIXTRACT® Total Histone Extraction Kit.

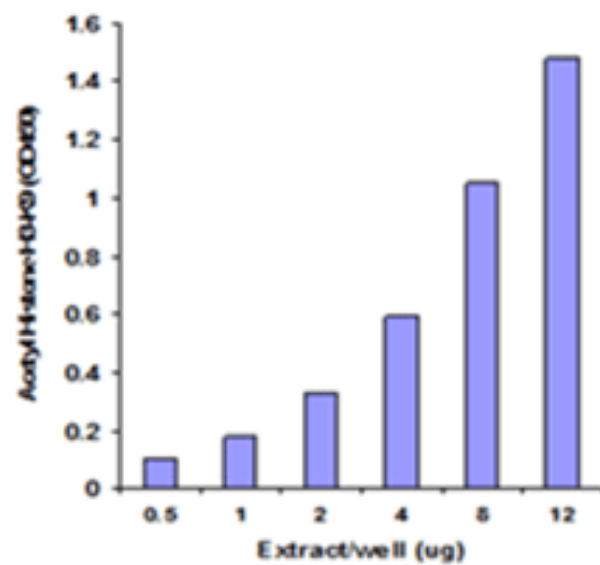


Figure 2. Histone extracts were prepared from MCF-7 cells using the EPIXTRACT® Total Histone Extraction Kit and acetyl histone H3-K9 was quantified.

## 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 or tissues.

Assay Time                    1 hour

Contents                        10x Pre-Lysis Buffer, Lysis Buffer, Balance Buffer, 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)