EPIXTRACT® Total Histone Extraction Kit

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

EPIXTRACT[®] 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 10⁵cells or 1mg tissue

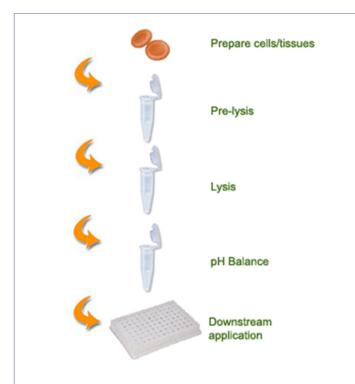


Fig 1. Schematic procedure for suing the $\mathsf{EPIXTRACT}^{\mathsf{®}}$ Total Histone Extraction Kit.

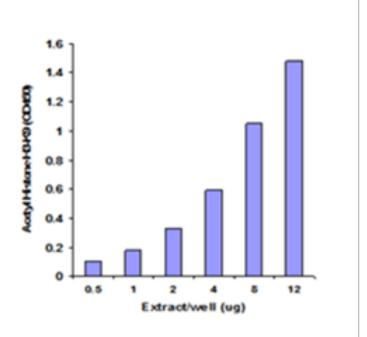


Figure 2. Histone extracts were prepared from MCF-7 cells using the EPIXTRACT[®] Total Histone Extraction Kit and acetyl histone H3-K9 was quantified.

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

Last modified: May 29, 2024



eu@enzolifesciences.com