## **EPIXTRACT**<sup>®</sup> Total Histone Extraction Kit (High Throughput)

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

FOIXTPACT® That district in the district in the coughput 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



- 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 3x10<sup>4</sup>cells

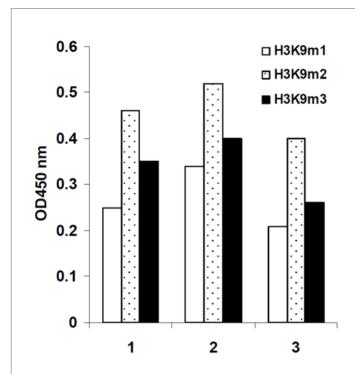


Figure 2. Histone extracts were prepared from MCF-7 cells (30,000 cells/well) using the EPIXTRACT<sup>®</sup> Total Histone Extraction Kit (High Throughput) and acetyl histone H3 was quantified.

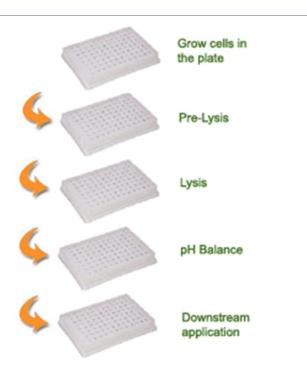


Figure 1. Schematic procedure for using the EPIXTRACT® Total Histone Extraction Kit (High Throughput)

## **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



info-

eu@enzolifesciences.com