

# Certificate of Analysis

## Histone H3

**Expression System:** (Recombinant protein expressed in E. coli cells)

**Catalog Numbers:** 12-800, 12-800-K, 12-800M

**Parent Lot Number:** D22AP010N

## Product Description

N-terminal 6His-tagged, recombinant, human Histone H3 amino acids 2-end expressed in E. coli cells. Purified using immobilized metal affinity chromatography. Molecular Weight = 17.5 kDa.

## Purity

95% by SDS-PAGE and Coomassie blue staining.

## Formulation

2.63 mg/ml of protein in 10 mM Sodium phosphate buffer, pH 7.4, 130 mM NaCl, 10% glycerol, 0.2mM PMSF, 0.1% 2-mercaptoethanol. Frozen solution.

## Storage and Stability

On receipt of material store at -70°C. Unopened reagent is stable for a minimum of 1 year from date of shipment when stored at recommended storage temperature. Avoid repeat freeze/thaw cycles. For maximum recovery of product, centrifuge original vial prior to removing the cap.

## Handling Recommendations

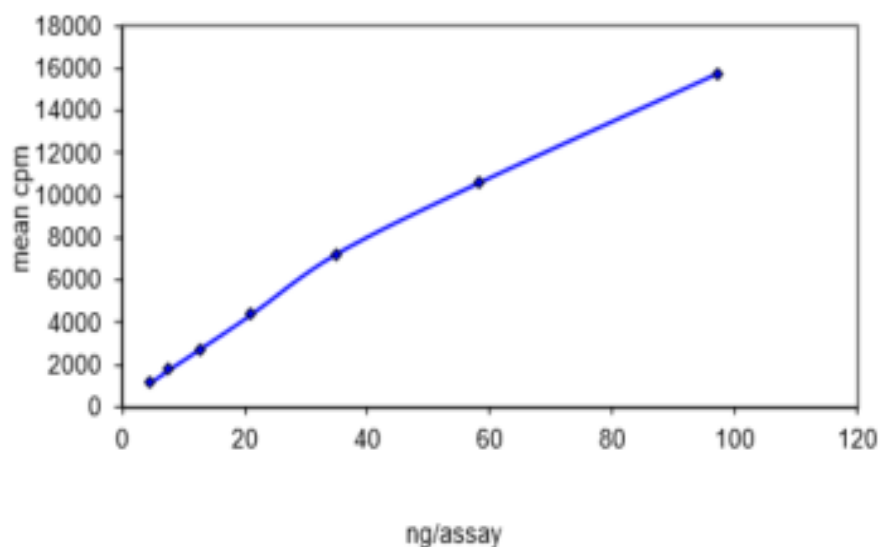
Rapidly thaw the vial under cold water and immediately place on ice. Aliquot unused material into pre-chilled micro-centrifuge tubes and immediately snap-freeze the vials in liquid nitrogen prior to re-storage at -70°C.

## Quality Control Testing

### Kinase Assay:

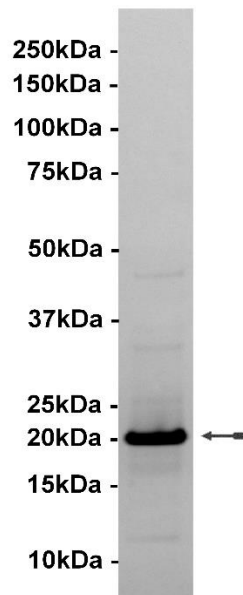
At 230  $\mu$ M, this lot was successfully phosphorylated by GAK (Catalogue # 16-055).

**Histone H3 data**



MS Tryptic Fingerprint: Confirmed identity as Histone H3 with the translated sequence.

SDS-PAGE and Coomassie Stain: Purity was assessed by SDS-PAGE and Coomassie blue staining using 3 µg of Histone H3.



The data presented in this document apply to the parent lot shown above and to all pack sizes derived from subsequent vialling runs of this parent lot. An alphabetical suffix after the parent lot number is used to denote each vialling run.

## Sequence Information

<b>Protein</b>	Human Histone H3
<b>Tags</b>	N-terminal 6His
<b>Native sequence</b>	A21 of the recombinant protein is equivalent to A2 of human Histone H3
<b>Accession number</b>	GenBank BC031333.1

### ***For Research Use Only***

Unless otherwise stated in our catalogue or other company documentation accompanying the product(s), our products are intended for research use only and are not to be used for any other purpose, which includes but is not limited to, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses or any type of consumption or application to humans or animals.

**Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)**  
**Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)**  
**General product information: [www.discoverx.com](http://www.discoverx.com)**