

PRODUCT INFORMATION



Histone H4 (human recombinant)

Item No. 10264

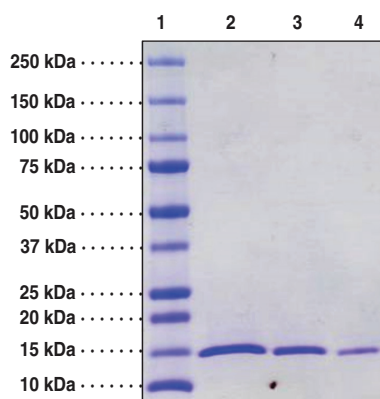
Overview and Properties

Synonyms: H4, Histone H4
Source: Recombinant protein, expressed in *E. coli*
Amino Acids: 1-103 (full-length)
Uniprot No.: P62805

Batch specific information can be found on the Certificate of Analysis or by contacting Technical Support

Molecular Weight: 11.5 kDa
Storage: -80°C (as supplied)
Stability: As supplied, 6 months from the QC date provided on the Certificate of Analysis, when stored properly
Purity: ≥95% (estimated by SDS-PAGE)
Supplied in: A lyophilized powder; recommended buffer for reconstitution is 50 mM sodium phosphate, pH 7.2, containing 100 mM sodium chloride and 20% glycerol.

Image(s)



Lane 1: MW Markers
Lane 2: Purified H4 (2 µg)
Lane 3: Purified H4 (5 µg)
Lane 4: Purified H4 (10 µg)

Representative gel image shown; actual purity may vary between each batch but protein will be ≥95% pure.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/02/2016

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

A nucleosome is the basic repeating unit of chromatin in which 146 base pairs of DNA wrap twice around a histone octamer consisting of two copies of each of the core histones, H2A, H2B, H3, and H4.¹ The combination of two H2A/H2B dimers and one H3/H4 tetramer create the nucleosome core.² Histone H4 undergoes many modifications which include acetylation, methylation, and phosphorylation that are important for regulation of gene transcription.¹

References

1. Bhaumik, S.R., Smith, E., and Shilatifard, A. Covalent modifications of histones during development and disease pathogenesis. *Nat. Struct. Mol. Biol.* **14(11)**, 1008-1016 (2007).
2. Tanaka, Y., Tawaramoto-Sasanuma, M., Kawaguchi, S., *et al.* Expression and purification of recombinant human histones. *Methods* **33**, 3-11 (2004).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM