

PRODUCT INFORMATION



Citrullinated Histone H2B (human, recombinant)

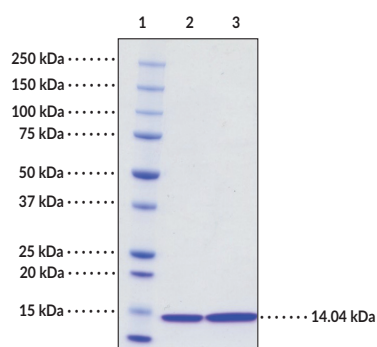
Item No. 30133

Overview and Properties

Synonyms: H2BC1, HIST1H2BA, hTSH2B, TH2B, TSH2B.1
Source: Recombinant human histone H2B type 1-A expressed in *E. coli*, citrullinated with human recombinant PAD4
Amino Acids: 2-127 (full length)
Uniprot No.: Q96A08
Molecular Weight: 14.04 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: Water
Protein Concentration: *batch specific* mg/ml

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

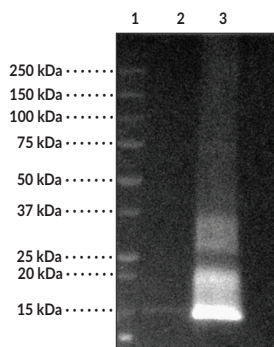
Images



Lane 1: MW Markers
Lane 2: Citrullinated Histone H2B (2 µg)
Lane 3: Citrullinated Histone H2B (4 µg)

SDS-PAGE analysis of Citrullinated Histone H2B.

Representative gel image shown; actual purity may vary between each batch.



Lane 1: MW Markers
Lane 2: Histone H2B (2 µg)
Lane 3: Citrullinated Histone H2B (4 µg)

Analysis of Histone H2B Citrullination. Histone H2B and citrullinated histone H2B were reacted with Cayman's Citrulline-specific Probe-biotin (Item No. 17450) and detected using Streptavidin:HRP (Item No. 16747).

Representative gel image shown; actual purity may vary between each batch.



Identification of modified sites in Citrullinated Histone H2B (Item No. 30133). Citrullinated Histone H2B was detected by LC-MS/MS and analyzed using Mascot and Scaffold PTM software. Deiminated arginines are indicated in teal.

Citrullination sites shown are representative of typical results. Batch-to-batch variations may occur.

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, 10/22/2020

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

Citrullinated Histone H2B Type 1-A is a citrullinated form of human histone H2B type 1-A (hTSH2B), a testis- and sperm-specific variant of the histone H2B protein that is encoded by *H2BC1*, previously known as *HIST1H2BA*.^{1,2} hTSH2B shares 85% identity with human somatic H2B.1, with the majority of amino acid differences between hTSH2B and somatic H2B occurring at the N-terminal end of the protein.¹ *H2BC1* is expressed in testis, and protein levels of hTSH2B are elevated in early round spermatids, decreasing over the course of spermiogenesis, with hTSH2B present in approximately 20% of mature spermatozoa.^{1,3} hTSH2B is enriched at promoter regions of genes important for sperm biology, capacitation, and fertilization and localizes to the basal nuclear area in mature spermatozoa.^{1,4} When incubated in *Xenopus* egg extracts, hTSH2B-positive human sperm cells decondense more rapidly and have increased long nuclear axis lengths compared with hTSH2B-negative sperm cells.⁵ Mice lacking the genes encoding the murine homolog of hTSH2B (TH2B), as well as the histone variant TH2A (*Th2a*^{-/-}*Th2b*^{-/-} mice), are sterile and have spermatogenesis defects.⁶ This product contains purified histone H2B type 1-A (human, recombinant) (Item No. 11081) that has been modified with PAD4 enzyme, which is subsequently depleted by affinity chromatography. Cayman's Citrullinated Histone H2B Type 1-A (human, recombinant) can be used for Western blot and ELISA applications.

References

1. Zalensky, A.O., Siino, J.S., Gineitis, A.A., *et al.* Human testis/sperm-specific histone H2B (hTSH2B). Molecular cloning and characterization. *J. Biol. Chem.* **277**(45), 43474-43480 (2002).
2. Marzluff, W.F., Gongidi, P., Woods, K.R., *et al.* The human and mouse replication-dependent histone genes. *Genomics* **80**(5), 487-498 (2002).
3. van Rooijen, H.J., Ooms, M.P., Spaargaren, M.C., *et al.* Immunoexpression of testis-specific histone 2B in human spermatozoa and testis tissue. *Hum. Reprod.* **13**(6), 1559-1566 (1998).
4. Hammoud, S.S., Nix, D.A., Zhang, H., *et al.* Distinctive chromatin in human sperm packages genes for embryo development. *Nature* **460**(7254), 473-478 (2009).
5. Singleton, S., Mudrak, O., Morshedi, M., *et al.* Characterisation of a human sperm cell subpopulation marked by the presence of the TSH2B histone. *Reprod. Fertil. Dev.* **19**(2), 392-397 (2007).
6. Shinagawa, T., Huynh, L.M., Takagi, T., *et al.* Disruption of *Th2a* and *Th2b* genes causes defects in spermatogenesis. *Development* **142**(7), 1287-1292 (2015).