

# PRODUCT INFORMATION



## Histone H2B Type 1-A (human, recombinant)

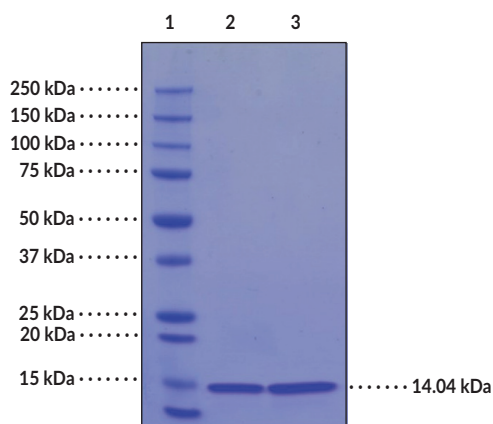
Item No. 11081

### Overview and Properties

**Synonyms:** H2BC1, HIST1H2BA, hTSH2B, TH2B, TSH2B.1  
**Source:** Recombinant human histone H2B type 1-A expressed in *E. coli*  
**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.

### Image



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

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

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, 01/22/2020

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

---

Human histone H2B type 1-A (hTSH2B) is a testis- and sperm-specific variant of the histone H2B protein that is encoded by *H2BC1*, previously known as *HIST1H2BA*.<sup>1,2</sup> 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.<sup>1</sup> *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.<sup>1,3</sup> 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.<sup>1,4</sup> 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.<sup>5</sup> Mice lacking the genes encoding the murine homolog of hTSH2B (TH2B), as well as the histone variant TH2A (*Th2a*<sup>-/-</sup>*Th2b*<sup>-/-</sup> mice), are sterile and have spermatogenesis defects.<sup>6</sup> Cayman's 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).

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