

Recombinant Histone TH2B

Catalog No: 31577, 31977

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

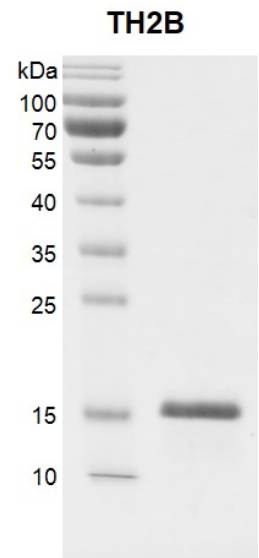
Buffer Contents: Recombinant Histone TH2B is supplied in 50 mM Tris pH 8.0, 150 mM NaCl, 5% glycerol. Please refer to product insert upon arrival for lot-specific concentration.

Background: TH2B (testis-specific H2B) is originally identified as a testis-specific variant of histone H2B. TH2A is a testis-specific variant of histone H2A. They are abundant in the testis, oocytes and fertilized eggs. TH2A and TH2B are controlled by a shared promoter that is located between them. Disruption of Th2a and Th2b genes causes defects in spermatogenesis. TH2A/TH2B may enhance reprogramming by introducing processes that normally operate in zygotes and during SCNT. TH2A and TH2B induce nucleosome instability. TH2B controls the chromatin-to-nucleoprotamine transition.

Protein Details: Recombinant Histone TH2B (accession number: NP_733759.1) was expressed in *E. coli* cells as full length protein without any tag. The molecular weight of TH2B is about 14.2 kDa. The recombinant histone is >95% pure by SDS-PAGE.

Application Notes: Recombinant histone TH2B is suitable for use as substrate for histone modification enzymes, or to generate chromatin *in vitro*.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant Histone TH2B gel. Histone TH2B was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.