

## Recombinant Histone H2BFWT

**Catalog No:** 31578, 31978

**Expressed In:** *E. coli*

**Quantity:** 100, 1000 µg

**Concentration:** 0.5 µg/µl

**Source:** Human

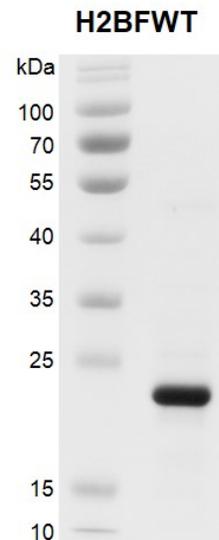
**Buffer Contents:** Recombinant Histone H2BFWT is supplied in 50 mM Tris pH 8.0, 150 mM NaCl, 5% glycerol. Please refer to product insert upon arrival for lot-specific concentration.

**Background:** H2BFWT (H2B histone family member W, testis specific) has been found in sperm nuclei and appears to be associated with the telomeric chromatin. In contrast to conventional H2B, H2BFWT is unable to recruit chromosome condensation factors and to participate in the assembly of mitotic chromosomes.

**Protein Details:** Recombinant Histone H2BFWT (accession number: NP\_001002916.3) was expressed in *E. coli* cells as full length protein without any tag. The molecular weight of H2BFWT is about 19.5 kDa. The recombinant histone is >92% pure by SDS-PAGE.

**Application Notes:** Recombinant histone H2BFWT 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 H2BFWT gel.** Histone H2BFWT was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.