

## Recombinant BRDT (257-382), GST-tag

---

**Catalog No:** 31483, 31883

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

**Quantity:** 100, 1000 µg

**Concentration:** 1 µg/µl

**Source:** Human

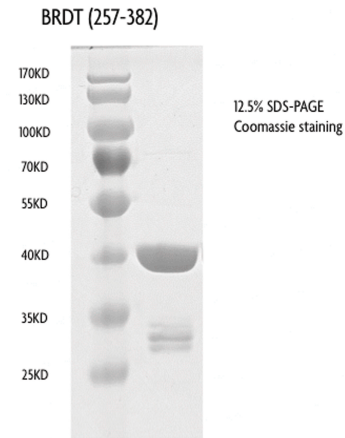
**Buffer Contents:** Recombinant BRDT (257-382), GST-tag protein expressed in *E. coli* at a concentration of 1 µg/µl in 25 mM Tris pH 7.4, 150 mM NaCl and 5% glycerol.

**Background:** BRDT is similar to the RING3 protein family. It possesses two bromodomain motifs and a PEST sequence (a cluster of proline, glutamic acid, serine, and threonine residues), characteristic of proteins that undergo rapid intracellular degradation. The bromodomain functions as a 'reader' of epigenetic histone marks and regulates chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets.

**Protein Details:** The peptide corresponding to amino acids 257-382 that contains the second bromodomain sequences of BRDT (accession number NP\_997072.1) was expressed in *E. coli* and contains an N-terminal GST tag with an observed molecular weight of 42.8 kDa. The recombinant protein is >90% pure by SDS-PAGE.

**Application Notes:** Recombinant BRDT (257-382), GST-tag is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

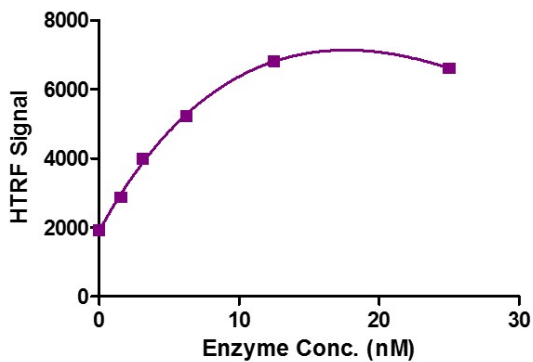
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product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



**Recombinant BRDT (257-382), GST-tag protein gel.**

BRDT (257-382), GST-tag protein was run on an SDS-PAGE gel and stained with Coomassie blue.

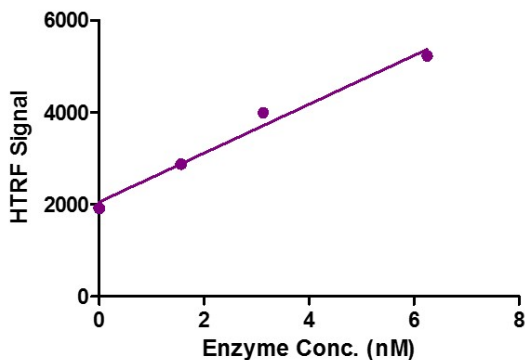
**GST-BRDT (257-382) Titration**



**Recombinant BRDT (257-382) HTRF activity assay.**

3.3  $\mu$ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRDT (257-382) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-GST antibody was used to detect reaction products.

**GST-BRDT (257-382) Titration**



**Recombinant BRDT (257-382) HTRF activity assay.**

3.3  $\mu$ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRDT (257-382) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-GST antibody was used to detect reaction products.