

Recombinant BRDT (21-137) protein, GST-Tag

Catalog No: 81156, 81856

Lot No: 13818001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.5 µg/µl

Source: Human

Buffer Contents: Recombinant BRDT (21-137) protein, GST-Tag is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol.

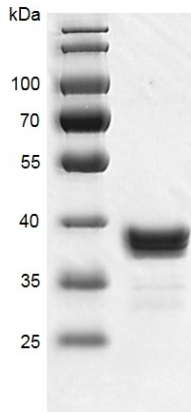
Background: BRDT is similar to the RING3 protein family. It possesses two bromodomain motifs and a PEST sequence (a cluster of proline, glutamic acid, serine, and threonine residues), characteristic of proteins that undergo rapid intracellular degradation. The bromodomain functions as a 'reader' of epigenetic histone marks and regulates chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. It shows binding specificity for acetylated H2BK12/15, H3K9/14/18/23, H4K12, H4K5/8, H4K8/12, H4K12/16, H4K5/12, H4K5/8/12/16 and H4K5/8/12/16.

Protein Details: Recombinant BRDT (21-137) protein, GST-Tag that includes amino acids 21-137 (Bromo domain) of human BRDT protein (accession number NP_001229735.2) was expressed in *E. coli* and contains an N-terminal GST Tag with a molecular weight of 40.2 kDa.

Application Notes: This protein 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.

GST-BRDT (21-137)



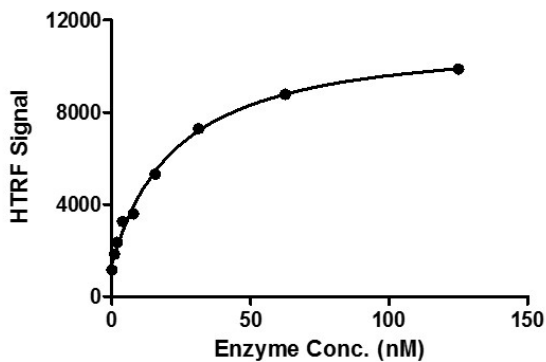
Recombinant BRDT (21-137) protein, GST-Tag protein gel

10% SDS-PAGE Coomassie staining

MW: 40.2 kDa

Purity: >80%

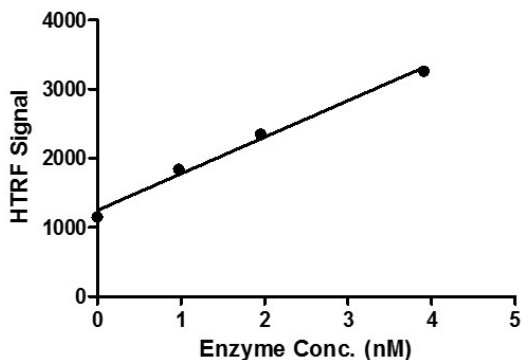
GST-BRDT (21-137)



HTRF assay for recombinant BRDT (21-137) protein, GST-Tag activity

1 μ M H4K5/8/12/16(ac4) peptide was incubated with different concentrations of BRDT (21-137) protein, GST-tag in a 10 μ l reaction system containing 50 mM HEPES-NaOH pH 7.4, 0.1% BSA for 1 hour, then 10 μ l GST antibody and SA-XL665 mixture (each 1:100 dilution in the reaction buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.

GST-BRDT (21-137)



HTRF assay for recombinant BRDT (21-137) protein, GST-Tag activity

1 μ M H4K5/8/12/16(ac4) peptide was incubated with different concentrations of BRDT (21-137) protein, GST-tag in a 10 μ l reaction system containing 50 mM HEPES-NaOH pH 7.4, 0.1% BSA for 1 hour, then 10 μ l GST antibody and SA-XL665 mixture (each 1:100 dilution in the reaction buffer) was added to each reaction system and incubated for 30 min. All the operations and reactions were performed at room temperature. HTRF assay was used for detection.