

Recombinant BRDT (21-137) protein

Catalog No: 31450, 31850

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant BRDT (21-137) protein is supplied at a concentration of 1 µg/µl in 25 mM Tris pH 8.0, 500 mM NaCl, 20% 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.

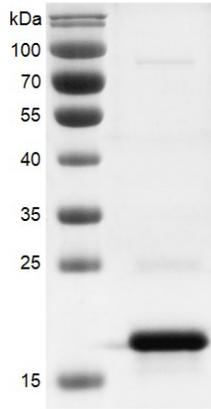
Protein Details: Recombinant BRDT (21-137) protein that includes amino acids 21-137 of human BRDT protein (accession number NP_001717.2) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with a molecular weight of 20 kDa. The purity of the protein is >85% by SDS-PAGE. 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.

Application Notes: Recombinant BRDT (21-137) 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 guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

BRDT (21-137)



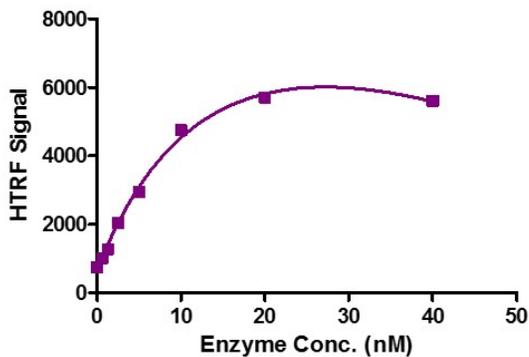
Recombinant BRDT (21-137) protein gel.

BRDT (21-137) protein was run on a 12% SDS-PAGE gel and stained with Coomassie Blue.

MW: 20 kDa

Purity: > 85%

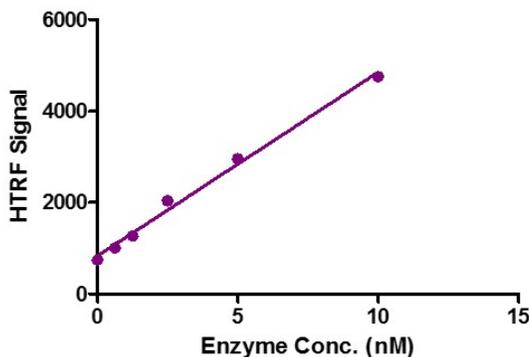
BRDT (21-137) Titration



HTRF Assay for Recombinant BRDT (21-137) protein activity.

3 μ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRDT (21-137) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA overnight at room temperature. Anti-Flag antibody was used to detect reaction products.

BRDT (21-137) Titration



HTRF Assay for Recombinant BRDT (21-137) protein activity.

3 μ M histone peptide H4K5/8/12/16(4Ac) was incubated with BRDT (21-137) protein in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA overnight at room temperature. Anti-Flag antibody was used to detect reaction products.