

BRD3 Bromodomain 1

Catalog No. 15-0011
Lot No. 13197001
Pack Size 100 µg

Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 41 kDa **Epitope Tag** GST

Product Description:

Recombinant human BRD3 Bromodomain 1 (RING3L, accession number Q15059, amino acids 25 to 160) containing an N-terminal GST tag, expressed in *E. coli*. BRD3 is a member of the BET sub-family of bromodomain-containing proteins, which also includes BRD2, BRD4 and BRDT. BET proteins have two bromodomains and an ET domain C-terminal to the second bromodomain. BRD3 interacts with acetylated histone and non-histone proteins, including acetylated GATA1. BRD3 is found fused to the NUT protein in a subset of carcinomas, contributing to the oncogenic phenotype by suppressing

Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

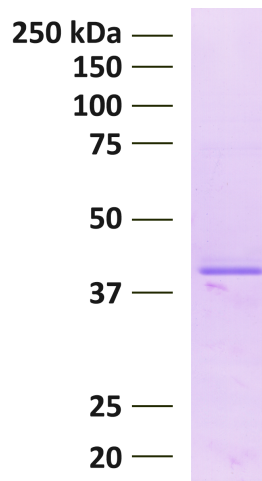
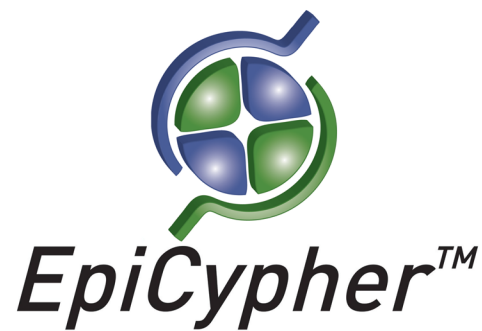
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Recombinant BRD3 Bromodomain 1 is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: BRD3 Bromodomain 1 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.

This product is for *in vitro* research use only and is not intended for use in humans or animals.