

BRD2 Bromodomain 1

Catalog No. 15-0009
Lot No. 13197001
Pack Size 100 µg

Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 42 kDa **Epitope Tag** GST

Product Description:

Recombinant human BRD2 Bromodomain 1 (RING3, accession number P25440, amino acids 57 to 200) containing an N-terminal GST tag, expressed in *E. coli*. BRD2 is a member of the BET sub-family of bromodomain-containing proteins, which also includes BRD3, BRD4 and BRDT. BET proteins have two bromodomains and an ET domain C-terminal to the second bromodomain. BRD2 has been reported to bind multiple sites of acetylation on histone H4. BRD2 is essential for mammalian development, and has been linked to Juvenile Myoclonic Epilepsy.

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.

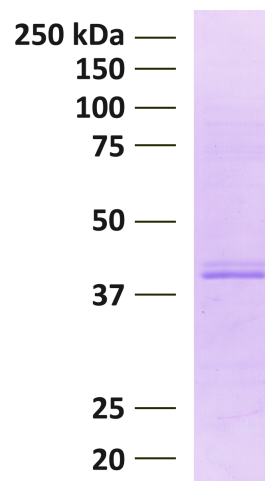
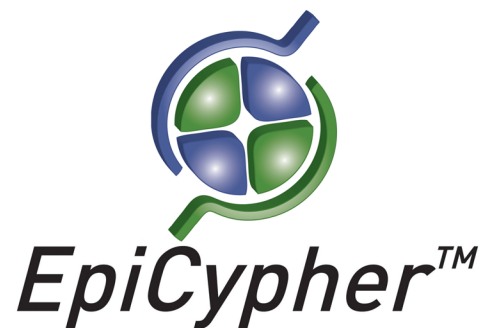
Storage and Stability:

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

Application Notes:

Recombinant BRD2 Bromodomain 1 is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: BRD2 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.