

BRD4 Bromodomain 2

Catalog No 15-0013
Lot No 22041005-01
Pack Size 100 µg

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

Product Description:

Recombinant human BRD4 Bromodomain 2 (HUNK1, accession O60885, amino acids 341-470) expressed in *E. coli* containing an N-terminal GST tag. BRD4 is a member of the BET sub-family of bromodomain-containing proteins, which also includes BRD2, BRD3 and BRDT. BET proteins have two bromodomains and an ET domain C-terminal to the second bromodomain. BRD4 serves as a cofactor for Myc and is crucial to Myc's role in acute myeloid leukemia. Inhibition of BRD4 binding suppresses Myc activity and leads to differentiation of AML cells and elimination of AML stem cells.

Formulation:

Recombinant GST-tagged protein at 1 mg/mL in 50 mM HEPES pH 7.3, 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 BRD4 Bromodomain 2 is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



EpiCypher®

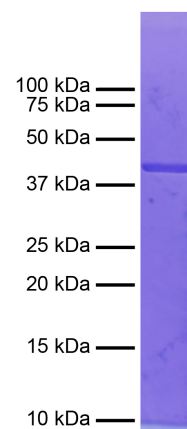


Figure 1: Protein gel data: BRD4 Bromodomain 2 (1 µg) was resolved via PAGE gel and stained with Coomassie blue to demonstrate the size and purity of the protein. The migration and molecular weight of the protein standards are indicated.

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