

BRD4 Bromodomain 2, His-Tagged

Catalog No. 15-0047
Lot No. 15022001
Pack Size 100 µg

Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 16 kDa **Epitope Tag** His

Product Description:

Recombinant human BRD4 Bromodomain 2, His-Tagged (HUNK1, accession number O60885, amino acids 341 to 470) containing an N-terminal His-tag, expressed in *E. coli*. 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 6xHis-tagged protein at 1 mg/ml in 20 mM Tris-HCl pH 8.0, 200 mM NaCl, 1 mM DTT, 20% glycerol.

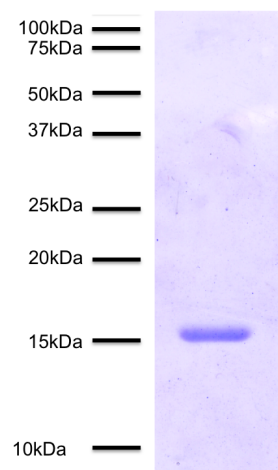
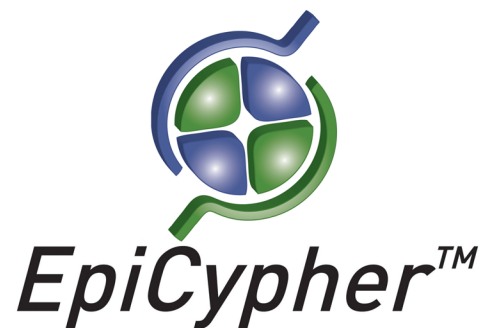
Storage and Stability:

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

Application Notes:

Recombinant BRD4 Bromodomain 2, His-Tagged is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: BRD4 Bromodomain 2, His-Tagged (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.