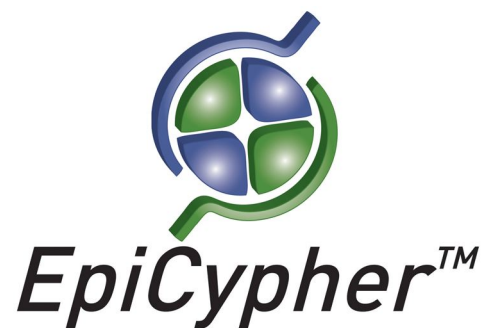


BRD9 Bromodomain, Recombinant Human

Catalog No. 15-0068
Lot No. 17041001
Pack Size 100 µg



Type Bromo **Expressed In** *E. coli*
Mol. Wgt. 37 kDa **Epitope Tag** GST

Product Description:

BRD9 Bromodomain, Recombinant Human (Bromodomain-containing protein 9, Rhabdomyosarcoma antigen MU-RMS-40.8, accession number Q9H8M2, amino acids 130 - 250) containing an N-terminal GST tag, expressed in *E. coli*. In addition to binding to acetylated histone proteins, the bromodomain (BD) of BRD9 has been reported to bind butyryllysine as well. BRD9 is reported to be part of SWI/SNF chromatin remodeling complexes in leukemic cells and inhibition of BRD9 BD binding has an anti-proliferative effect on these cells.

Formulation:

Recombinant GST-tagged protein at 1 mg/ml in 50 mM TRIS pH 8.4, 250 mM NaCl, 1 mM DTT, 20% glycerol.

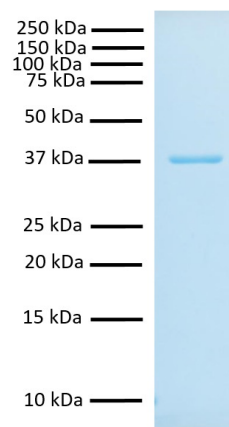
Storage and Stability:

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

Application Notes:

BRD9 Bromodomain, Recombinant Human is useful for protein binding and screening experiments examining acylated (acetyl and butyryl) protein and peptide substrates.

References:



Protein Gel Data: BRD9 Bromodomain, Recombinant Human (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.