

BAZ2B Bromodomain

Catalog No. 15-0042
Lot No. 13259001
Pack Size 100 µg

Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 39 kDa **Epitope Tag** GST

Product Description:

Recombinant human BAZ2B Bromodomain (WALp4, accession Q9UIF8, amino acids 2041 to 2168) containing an N-terminal GST tag, expressed in *E. coli*. BAZ2B is an bromodomain protein that has been shown to interact with acetylated lysines on histone H3. BAZ2B also contains a PHD finger and an MBD (methylated DNA binding) domain.

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.

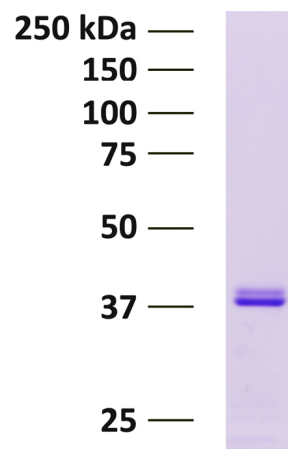
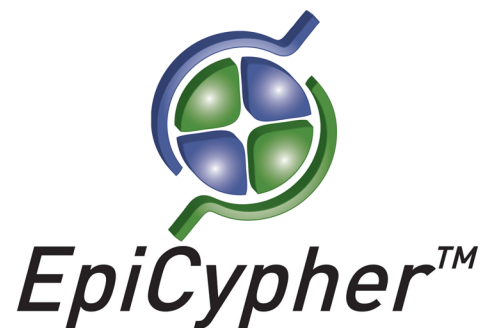
Storage and Stability:

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

Application Notes:

Recombinant BAZ2B Bromodomain is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: BAZ2B Bromodomain (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.