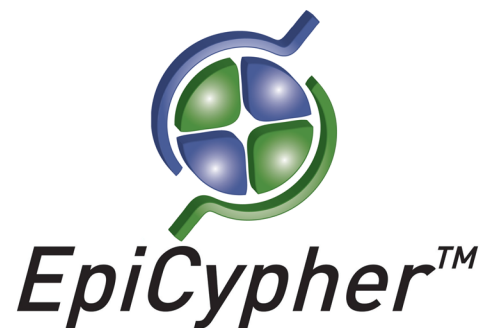


BATZFD Bromodomain

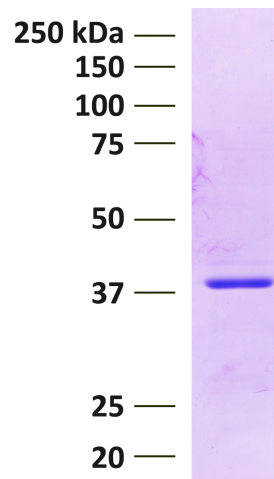
Catalog No. 15-0004
Lot No. 13182001
Pack Size 100 µg



Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 38 kDa **Epitope Tag** GST

Product Description:

Recombinant human BATZFD Bromodomain (Bromodomain adjacent to zinc finger domain, accession number Q59G54, amino acids 1021 to 1150) containing an N-terminal GST tag, expressed in *E. coli*.



Protein Gel Data: BATZFD 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.

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

References:

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