

ACF1 / BAZ1A Bromodomain

Catalog No. 15-0001
Lot No. 13182001
Pack Size 100 µg

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

Product Description:

Recombinant human ACF1 Bromodomain (BAZ1A, WCRF180, accession Q9NRL2, amino acids 1412 to 1551) containing an N-terminal GST tag, expressed in *E. coli*.

ACF1 is a component of the ACF complex, an ATP-dependent chromatin remodeling complex, that regulates spacing of nucleosomes. It is also part of the CHRAC complex which facilitates nucleosome sliding by the ACF complex and enhances ACF-mediated chromatin assembly. ACF1 may be involved in nuclear receptor-mediated repression of transcription through its interaction with NCOR1.

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.

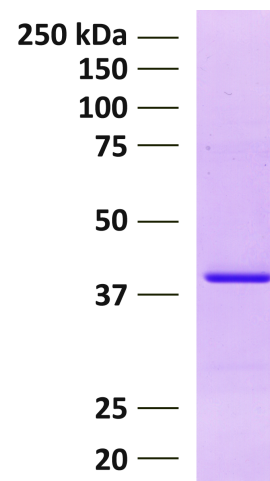
Storage and Stability:

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

Application Notes:

Recombinant ACF1 / BAZ1A Bromodomain is useful for protein binding and screening experiments examining acetylated protein substrates.

References:



Protein Gel Data: ACF1 / BAZ1A 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.