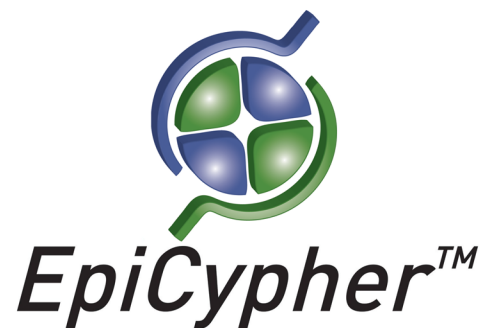


BRG1 / SMCA4 Bromodomain

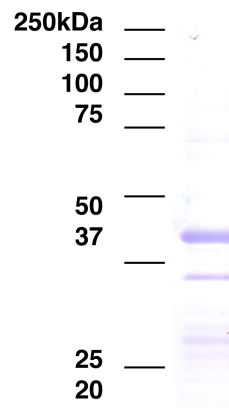
Catalog No. 15-0018
Lot No. 13228001
Pack Size 100 µg



Type Bromodom **Expressed In** *E. coli*
Mol. Wgt. 41 kDa **Epitope Tag** GST

Product Description:

Recombinant human BRG1 Bromodomain (Brahma-related gene 1, SMC4A, SMARCA4, BAF190A, BRG1, SNF2B, SNF2L4, accession number P51532 amino acids 1441 to 1580) containing an N-terminal GST tag, expressed in *E. coli*. BRG1 / SMCA4 is an ATP-dependent chromatin remodeling protein and a component of several protein complexes involved in transcriptional activation, including CREST and BAF. BRG1 is a component of an embryonic stem cell-specific BAF complex, and BRG1 is essential for ESC self-renewal and proliferation. Mutations in the BRG1 gene have been described in a



Protein Gel Data: BRG1 / SMCA4 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 BRG1 / SMCA4 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.