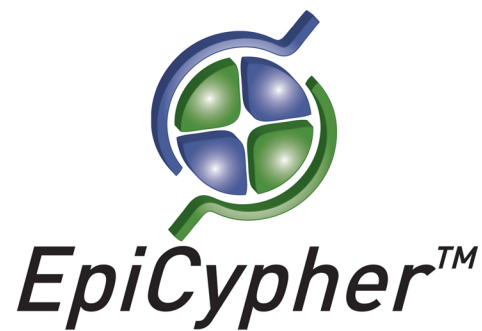


CECR2 Bromodomain

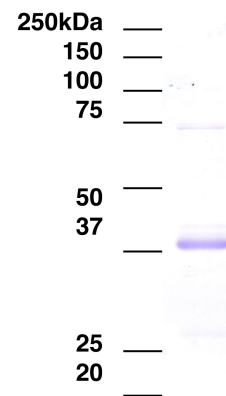
Catalog No. 15-0024
Lot No. 13228001
Pack Size 100 µg



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

Product Description:

Recombinant human CECR2 Bromodomain (Cat eye syndrome critical region protein 2, accession number Q9BXF3, amino acids 421 to 550) containing an N-terminal GST tag, expressed in *E. coli*. CECR2 is a member of the CERF chromatin remodeling complex, also containing SNF2L / SMARCA1. The bromodomain from CECR2 binds several sites of histone acetylation on H2A, H2B and H3. The CECR2 gene a candidate for Cat Eye Syndrome, a developmental defect resulting from a duplication of the region of chromosome 22 that contains CECR2.



Protein Gel Data: CECR2 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 CECR2 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.