

Recombinant CECR2 (425-538) protein

Catalog No: 31374, 31774

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1.5 µg/µl

Source: Human

Buffer Contents: Recombinant CECR2 (425-538) protein was expressed in *E. coli* cells at a concentration of 1.5 µg/µl in 25 mM Tris-HCl pH 8.0, 500 mM NaCl, 0.04% Triton X-100, 20% glycerol, 1mM TCEP.

Background: CECR2 (Cat Eye Syndrome Chromosome Region, Candidate 2) protein is the bromodomain-containing subunit of the CERF (CECR2-containing-remodeling factor) complex. Bromodomains recognize acetylated histone lysine residues and function as 'readers' of these epigenetic histone marks to regulate chromatin structure and gene expression by linking associated proteins to the recognized acetylated nucleosomal targets. The CERF complex functions to facilitate the perturbation of chromatin structure in an ATP-dependent manner. In addition to its chromatin remodeling function, CECR2 interaction with LRPPRC may be involved in the integration of the cytoskeletal network with processes such as vesicular trafficking, nucleocytoplasmic shuttling, transcription, chromosome remodeling and cytokinesis.

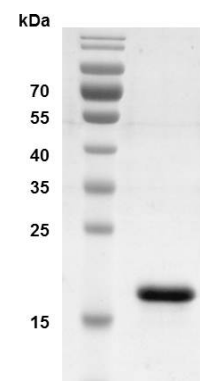
Protein Details: The peptide corresponding to amino acids 425 - 538 that contains the bromodomain sequences of CECR2 (accession number NM_031413.3) was expressed in *E. coli* and contains an N-terminal His tag and C-terminal FLAG tag with an observed molecular weight of 19.7 kDa. It shows binding specificity for acetylated H3K9, H3K14, H3K9/K14 and H3K9/S10/K14. The recombinant protein is >95% pure by SDS-PAGE.

Application Notes: Recombinant CECR2 (425-538) is suitable for use in binding assays, inhibitor screening, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

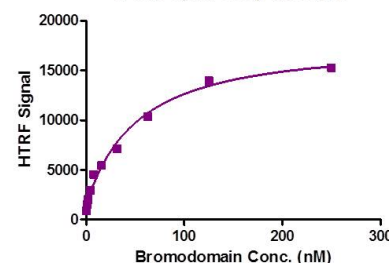
CECR2 (425-538)



Recombinant CECR2 (425-538) protein gel.

CECR2 (425-538) protein was run on a 12.5% SDS-PAGE gel and stained with Coomassie Blue.

CECR2 (425-538) Titration



Recombinant CECR2 (425-538) activity assay.

3 µM histone peptide H4K5/8/12/16 (ac4) was incubated with CECR2 (425-538) in reaction buffer including 50 mM HEPES-NaOH pH 7.0, 0.1% BSA for 1 hour at room temperature. Anti-FLAG antibody was used to detect reaction products.