

CECR2 (His)

(Cat eye syndrome critical region protein 2; KIAA1740)

CATALOG NO.: RD-11-210

LOT NO.:

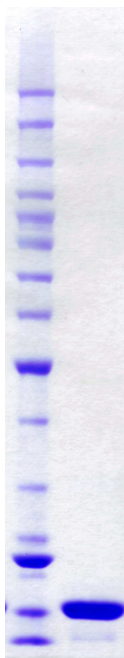
DESCRIPTION: Human recombinant CECR2 bromodomain (residues 425-538; Genbank Accession # NM_031413 ; MW = 16.4 kDa) expressed in *E. coli* with an N-terminal His-tag. CECR2, along with SMARCA1 (SNF2L) forms the ATP-dependent chromatin remodeling complex CERF and plays an essential role in neural tube formation¹. The CECR2 bromodomain displays binding affinity for multiple histone H2A and H3 Lys(Ac) residues in singly acetylated histone peptide microarrays (H2A: K15Ac, K36Ac, K75Ac; H3: K9Ac, K14Ac, K18Ac & others), with the H3K9Ac and H3K14Ac interactions confirmed by isothermal titration calorimetry (ITC)². A crystal structure for the CECR2 bromodomain has been determined².

PURITY: >95% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris-HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v) as determined by OD₂₈₀.

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 μl) or storage of diluted protein is not recommended.

REFERENCES: 1) G.S. Banting *et al. Hum. Mol. Genet.* 2005 **14** 513; 2) P. Filippakopoulos *et al. Cell* 2012 **149** 214



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 5 μg CECR2 (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355
 requests@reactionbiology.com www.reactionbiology.com