

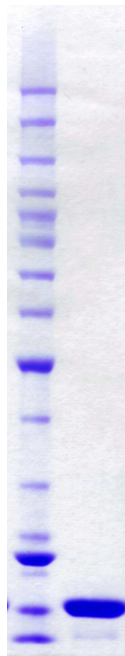
CECR2 (His) (Cat eye syndrome critical region protein 2; KIAA1740)

**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 formation<sup>1</sup>. 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)<sup>2</sup>. A crystal structure for the CECR2 bromodomain has been determined<sup>2</sup>.

**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 OD<sub>280</sub>.

**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  $\mu\text{l}$ ) 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  $\mu\text{g}$  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.