

## CHD4-[CHR] (His)

**CATALOG NO.:** RD-11-440

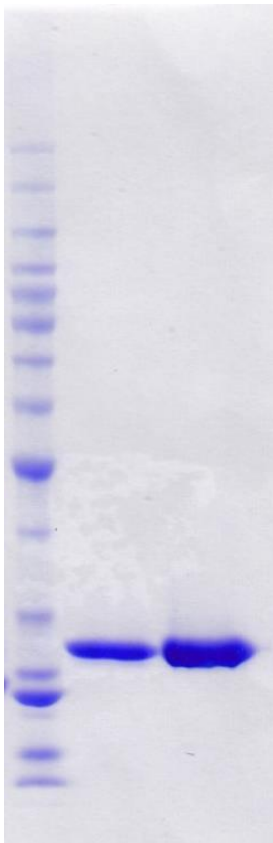
**LOT NO.:**

**DESCRIPTION:** Human recombinant CHD4-[CHR] (residues 497-680; Genbank Accession # NM\_001273.3; MW = 23.7 kDa) expressed as a C-terminal His-tag protein in *E. coli*.

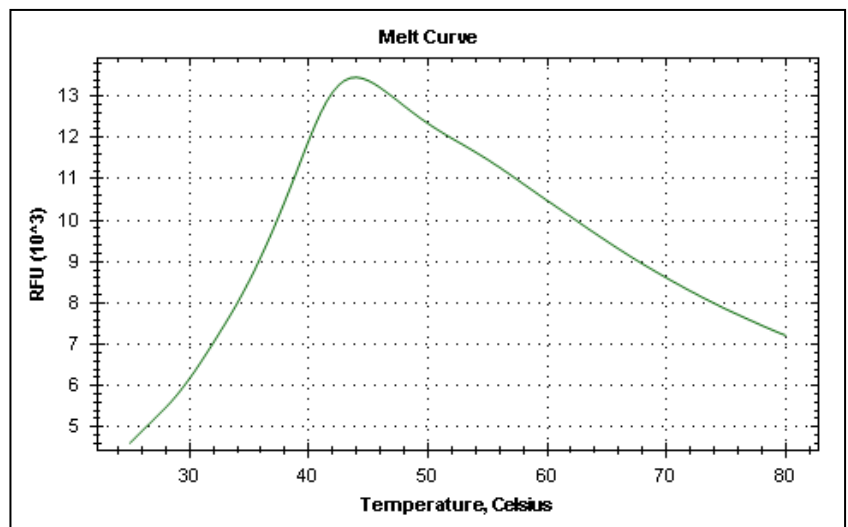
**PURITY:** >95% by SDS-PAGE

**SUPPLIED AS:**  $\_ \mu\text{g}/\mu\text{L}$  in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol 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 enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 and 10  $\mu\text{g}$  of RBC CHD4-[CHR] (His).** MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Differential Scanning Fluorimetry of RBC CHD4-[CHR] (His).** Thermal denaturation of CHD4-[CHR] (His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). The apo form of CHD4-[CHR] (His) displays a T<sub>m</sub> of 39.0°C.

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