

HP1 β -[CHR] (His)

{CBX1-[CHR] (His)}

CATALOG NO.: RD-11-203

LOT NO.:

DESCRIPTION: Human recombinant HP1 β -chromodomain (residues 19-79; Genbank Accession # NM_006807; MW = 9.98 kDa) expressed with an N-terminal His-tag in *E. coli*.

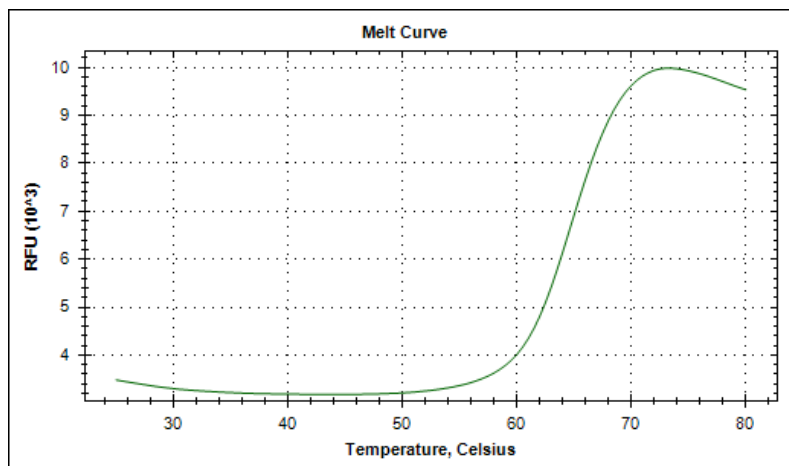
PURITY: >85% by SDS-PAGE

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 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 enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC HP1 β -[CHR] (His): MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Differential Scanning Fluorimetry of RBC HP1 β -[CHR] (His). Thermal denaturation of HP1 β -[CHR] (His) is detected (CFX384TM Touch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). HP1 β [CHR] (His) displays a T_m of 64.5°C.

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