

HP1 α -[CHR] (GST) {CBX5-[CHR] (GST)}

CATALOG NO.: RD-11-420

LOT NO.:

DESCRIPTION: Human recombinant HP1 α -[CHR] (residues 18-78; Genbank Accession # NM_001127322.1; MW = 34.5 kDa) expressed as an N-terminal GST-fusion 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 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 α -[CHR] (GST). 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.