

PB1-6 (His) (Polybromo-1, PBRM1 bromodomain-6)

CATALOG NO.: RD-11-233

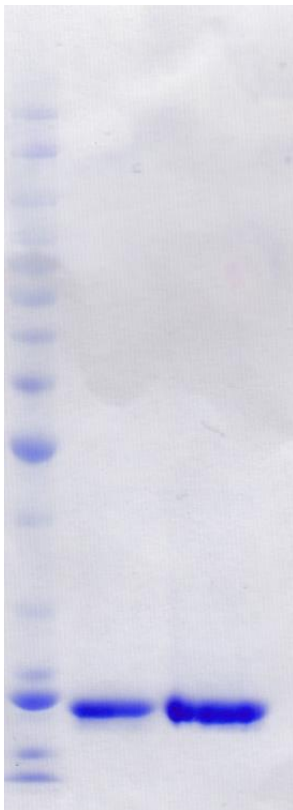
LOT NO.:

DESCRIPTION: Human recombinant PB1-6 bromodomain protein (residues 773-917; Genbank Accession # NM_018313; MW = 20.0 kDa) expressed with an N-terminal 6xHis-tag in *E. coli*.

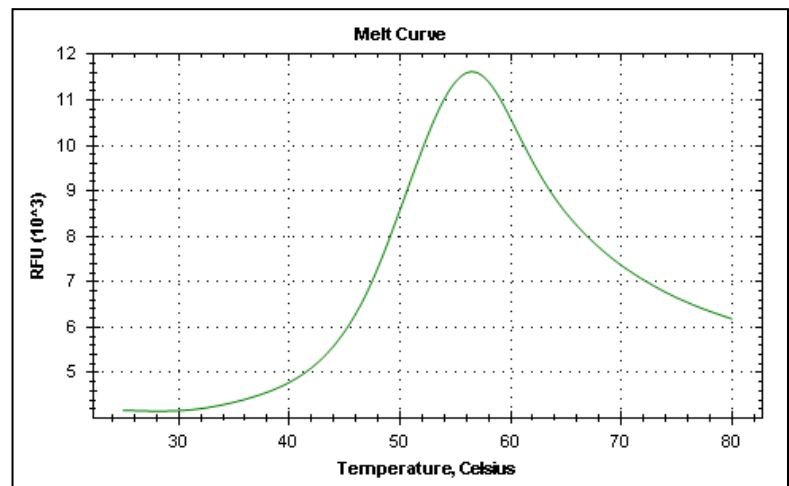
PURITY: >95% by SDS-PAGE

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 and 10 μg of RBC PB1-6 (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 PB1-6 (His). Thermal denaturation of PB1-6 (His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). Apo form of PB1-6 (His) displays a T_m of 50.5°C.

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