

BRD7 (His)

CATALOG NO.: RD-11-471

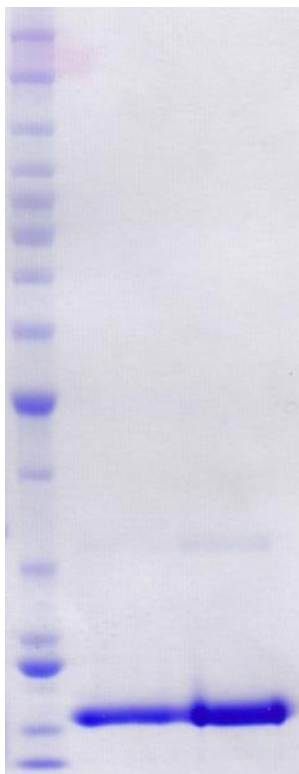
LOT NO.:

DESCRIPTION: Human recombinant BRD7 bromodomain (residues 129-254; Genbank Accession # NM_013263; MW = 17.5 kDa) expressed as an N-terminal 6xHis-fusion protein in *E. coli*.

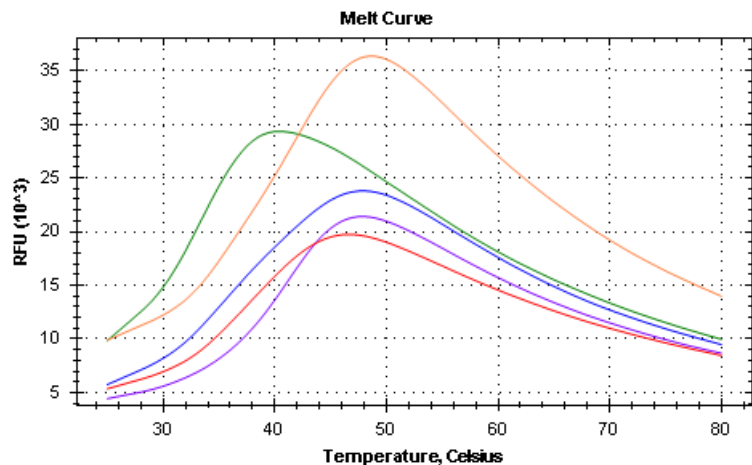
PURITY: >95% by SDS-PAGE

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 and 10 μg of RBC BRD7 (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 BRD7 (His) in presence or absence of common bromodomain ligands.

Thermal denaturation of BRD7 (His) is detected (CFX384 TMTouch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). Addition of 25 μM Bromosporine (blue), 25 μM BI-9564 (red), 25 μM NI-57 (orange) and 1% N-methyl-2-pyrrolidinone and (purple) stabilizes the protein folding and shifts the T_m (inflection point) from 33.5°C to 35.5°C, 38°C, 42°C and 41°C, respectively.

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