

BRD7 (His)

CATALOG NO.: RD-11-471

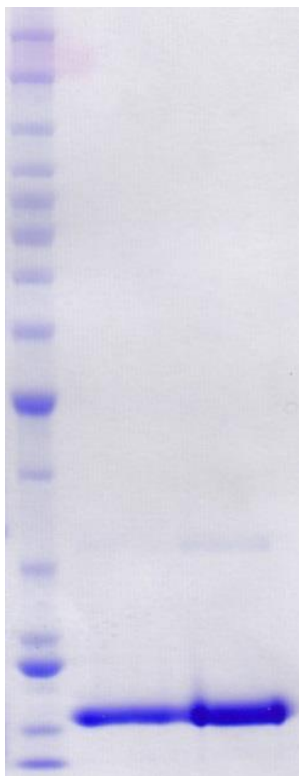
LOT NO.:

**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*.

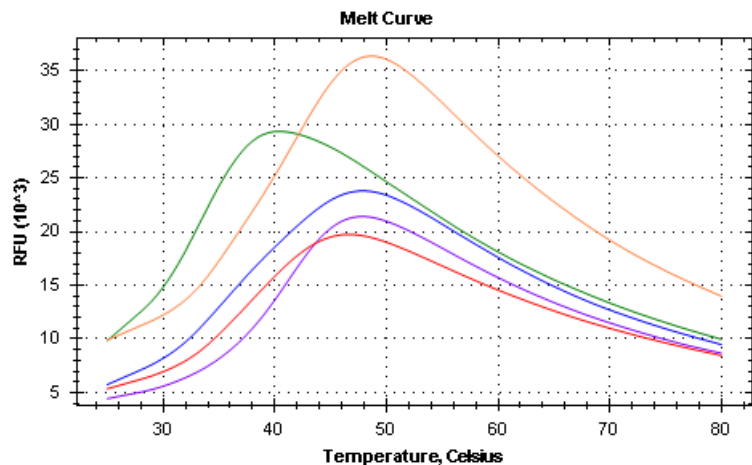
**PURITY:** >95% by SDS-PAGE

**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 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 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 $\mu\text{M}$  Bromosporine (blue), 25 $\mu\text{M}$  BI-9564 (red), 25 $\mu\text{M}$  NI-57 (orange) and 1% N-methyl-2-pyrrolidinone (purple) stabilizes the protein folding and shifts the T<sub>m</sub> (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.