

BRD7 (GST)

CATALOG NO.: RD-11-254

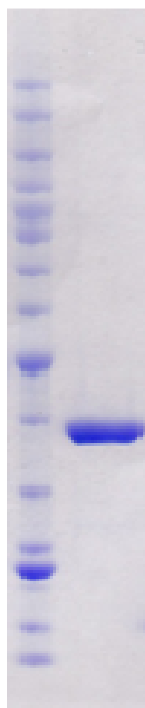
LOT NO.:

DESCRIPTION: Human recombinant BRD7 bromodomain (residues 129-254; Genbank Accession # NM_013263; MW = 41.83 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

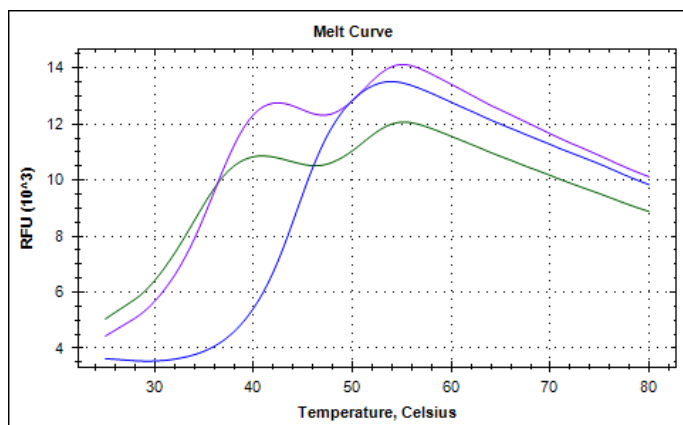
PURITY: >95% by SDS-PAGE

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 BRD7 (GST). 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 (GST) in presence or absence of common bromodomain ligands.

Thermal denaturation of BRD7 (GST) 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) and RVX-208 (purple) stabilizes the protein folding and shifts the T_m (inflection point) from 34°C to 44.5°C and 36°C, respectively.

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