

BRD8-1 (His)

(Bromodomain containing protein 8, bromodomain 1)

CATALOG NO.: RD-11-302

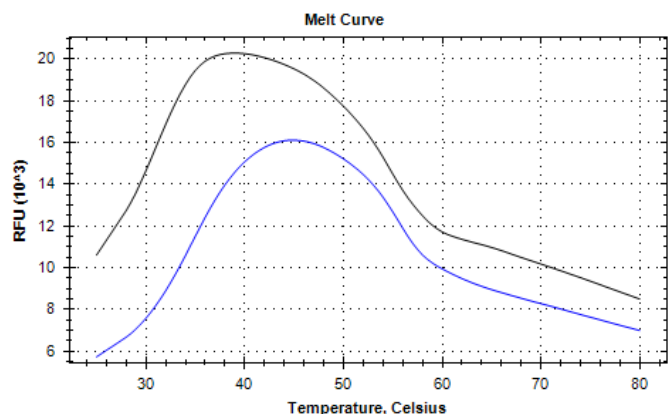
LOT NO.:

DESCRIPTION: Human recombinant BRD8, bromodomain 1 (residues 701-819; Genbank Accession # NM_139199.1; MW = 16.6 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

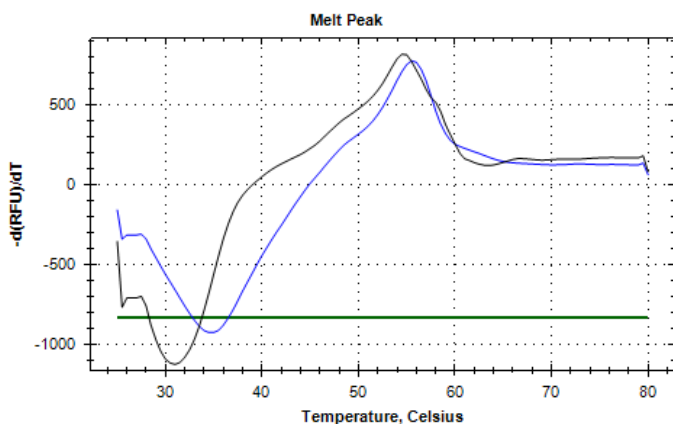
PURITY: >80% by SDS-PAGE

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.



Differential Scanning Fluorimetry of RBC BRD8-1 (His). Thermal denaturation of BRD8-1 (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 1% 1-methyl-2-pyrrolidone stabilizes the protein folding and shifts the T_m (inflection point) from 31°C to 34.5°C.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC BRD8-1 (His). 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.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com