

BRWD3-2 (GST)

CATALOG NO.: RD-11-301

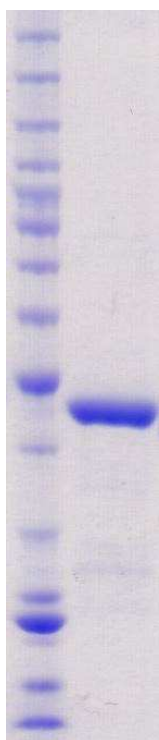
LOT NO.:

DESCRIPTION: Human recombinant BRWD3-2 bromodomain (residues 1292-1444; Genbank Accession # NM_153252; MW = 44.8 kDa) expressed in *E. coli* as an N-terminal GST fusion protein.

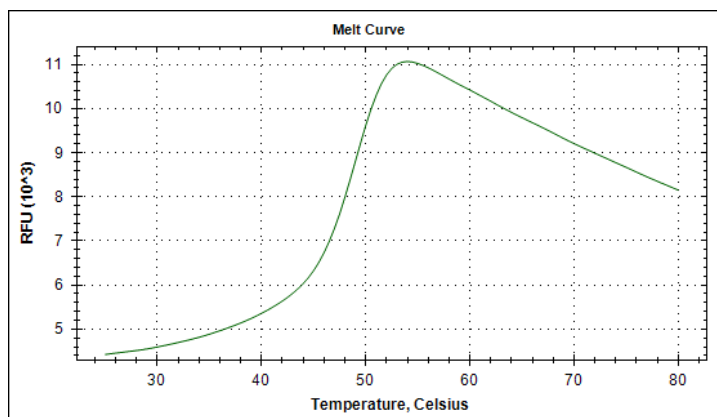
PURITY: >95% by SDS-PAGE

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 BRWD3-2 (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 BRWD3-2 (GST)

Thermal denaturation of BRWD3-2 (GST) is detected (CFX384TM Touch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies).

Apo form of BRWD3-2 (GST) displays a T_m of 49°C and is not stabilized in the presence of various known bromodomain ligands (JQ1, PF11, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208, GSK2801 and PF13; all tested at 25 μM).

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