

## BAZ2A (GST)

**CATALOG NO.:** RD-11-272

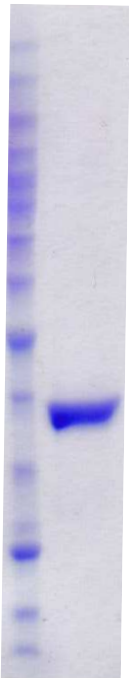
**LOT NO.:**

**DESCRIPTION:** Human recombinant BAZ2A bromodomain (residues 1787-1901; Genbank Accession # NM\_013449; MW = 40.7 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

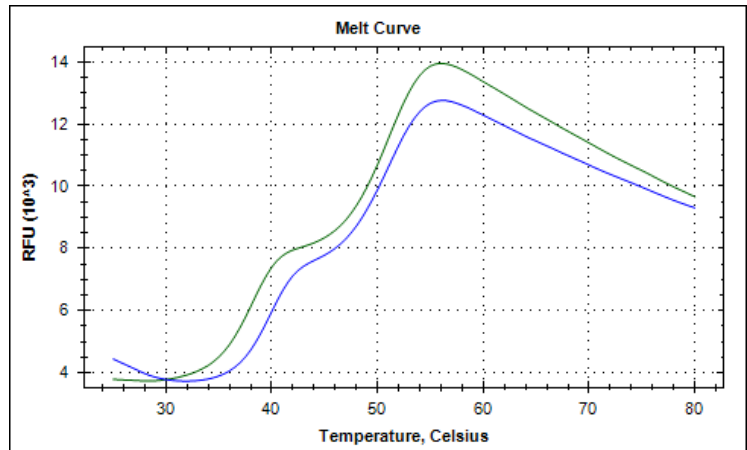
**PURITY:** >95% by SDS-PAGE

**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 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  $\mu\text{g}$  of RBC BAZ2A (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 BAZ2A (GST) in presence or absence of common bromodomain ligands.**

Thermal denaturation of BAZ2A (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  $\mu\text{M}$  Bromosporine (blue) stabilizes the protein folding and shifts the T<sub>m</sub> (inflection point) from 38°C to 40°C.

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