

BAZ1A (His)

CATALOG NO.: RD-11-225

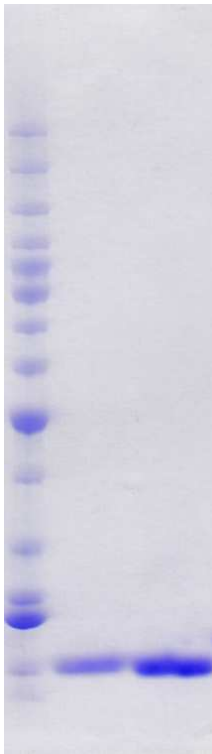
LOT NO.:

DESCRIPTION: Human recombinant BAZ1A bromodomain (residues 1423-1544; Genbank Accession # NM_013448; MW = 16.8 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

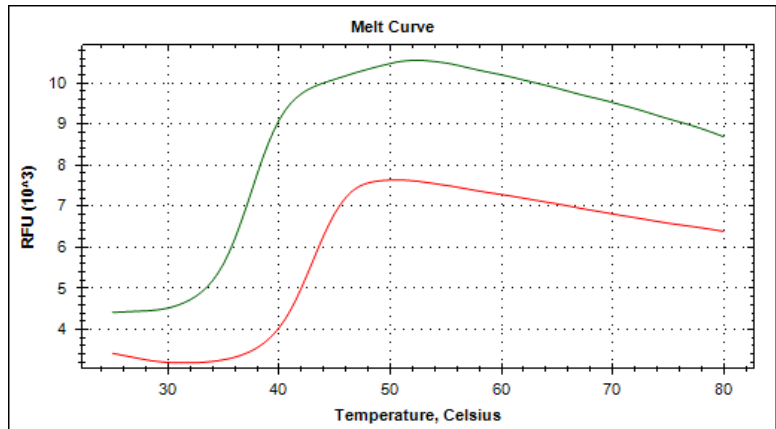
PURITY: >95% by SDS-PAGE

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 2 μg and 4 μg of RBC BAZ1A (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 BAZ1A (His).

Thermal denaturation of BAZ1A (His) is detected (CFX384TM Touch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). Addition of 25 μM GSK2801 (red) stabilizes the protein folding and shifts the T_m (inflection point) from 37.5°C to 43°C.

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