

BAZ1B (His)

(Bromodomain adjacent to zinc finger domain protein 1B; WSTF)

CATALOG NO.: RD-11-208

LOT NO.:

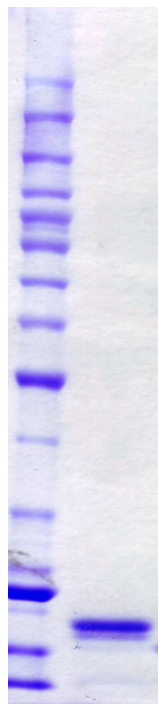
DESCRIPTION: Human recombinant BAZ1B bromodomain (residues 1340-1457; Genbank Accession # NM_032408; MW = 17.1 kDa) expressed in *E. coli* with C-terminal His- and Strep-tags. BAZ1B (WSTF) is a key subunit of several ATP-dependent chromatin remodeling complexes with activities in nucleosome and chromatin assembly, in transcriptional regulation (at vitamin D-receptor (VDR) target genes) and in DNA replication and repair (see review¹). The full-length protein comprises multiple domains, including the N-terminal domain which confers the atypical tyrosine kinase activity responsible for H2A.X Tyr142 phosphorylation, a key regulatory mark in the DNA damage response^{2,3}. The BAZ1B bromodomain displays affinity for several acetylated histone tail peptides, particularly ones incorporating H3K14(Ac)⁴. Interaction between BAZ1B and the gene promoter for the vitamin D synthetic enzyme 25(OH)D₃ 1 α -hydroxylase, most likely mediated by bromodomain binding to acetylated histones, is necessary for ligand-dependent transcriptional repression by VDR⁴. By virtue of its regulatory role in aromatase gene expression, BAZ1B may hold potential as a target for therapy in estrogen-dependent breast cancer⁵.

PURITY: >85% by SDS-PAGE

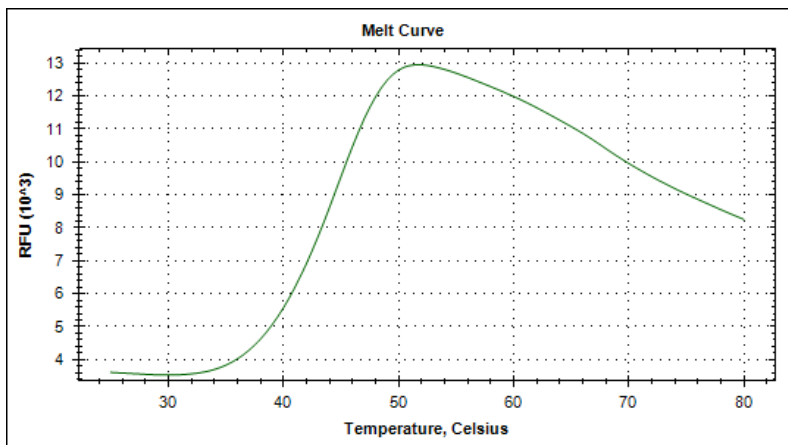
SUPPLIED AS: μ g/ μ L in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v) as determined by OD₂₈₀

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 μ l) or storage of diluted protein is not recommended.

REFERENCES: 1) C. Barnett & J.E. Krebs *Biochem. Cell Biol.* 2011 **89** 12; 2) A. Xiao *et al. Nature* 2009 **457** 57; 3) N. Singh *et al. Proc. Natl. Acad. Sci. USA* 2012 **109** 14381; 4) R. Fujiki *et al. EMBO J.* 2005 **24** 3881; 5) J. Lundqvist *et al. Biochim. Biophys. Acta* 2013 **1833** 40



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μ g of RBC BAZ1B (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 BAZ1B (His): Thermal denaturation of BAZ1B (His) is detected (CFX384TM Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). The apo form of BAZ1B (His) displays a T_m of 44.5°C and is not stabilized in the presence of various known bromodomain ligands (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151 and RVX-208; all tested at 25 μ M; not shown).

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