

BAZ1B (GST) (Bromodomain adjacent to zinc finger domain protein 1B; WSTF)

CATALOG NO.: RD-11-252

LOT NO.:

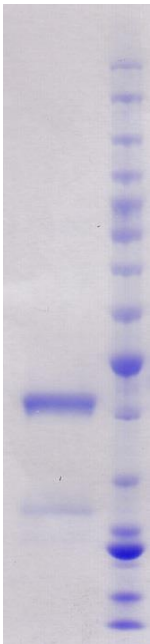
**DESCRIPTION:** Human recombinant BAZ1B bromodomain (residues 1340-1457; Genbank Accession # NM\_032408.2; MW = 44.1 kDa) as an N-terminal GST-fusion protein with C-terminal His- and Strep-tags expressed in *E. coli*. BAZ1B (WSTF) is a key subunit of several ATP-dependent chromatin remodeling complexes with activities in nucleosome and chromatin assembly, in transcriptional regulation (at vitamin D-receptor (VDR) target genes) and in DNA replication and repair (see review<sup>1</sup>). 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<sup>2,3</sup>. The BAZ1B bromodomain displays affinity for several acetylated histone tail peptides, particularly ones incorporating H3K14(Ac)<sup>4</sup>. Interaction between BAZ1B and the gene promoter for the vitamin D synthetic enzyme 25(OH)D<sub>3</sub> 1 $\alpha$ -hydroxylase, most likely mediated by bromodomain binding to acetylated histones, is necessary for ligand-dependent transcriptional repression by VDR<sup>4</sup>. By virtue of its regulatory role in aromatase gene expression, BAZ1B may hold potential as a target for therapy in estrogen-dependent breast cancer<sup>5</sup>.

**PURITY:** >80% by SDS-PAGE

**SUPPLIED AS:**  $\mu$ g/ $\mu$ L in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v)

**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  $\mu$ 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 (12% acrylamide) of 2 $\mu$ g of RBC BAZ1B (GST)**

MW markers (right) 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.