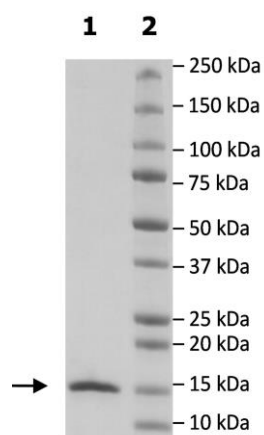


Product Information

Description:	Human Bromodomain adjacent to zinc finger domain, 2B or BAZ2B, GenBank Accession No. NM_013450, a.a. 2054 – 2168 (end) corresponding to single bromodomain with N-terminal His-tag, MW = 14.3 kDa, expressed in an <i>E. coli</i> expression system.
Species	Human
Construct:	BAZ2B (His-2054-2168(end))
Concentration:	1.34 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	14 kDa
Genbank Accession:	NM_013450
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was done according to BAZ2B Inhibitor Screening Assay Kit (BPS Bioscience #32600) with BAZ2B titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



BAZ2B Activity

