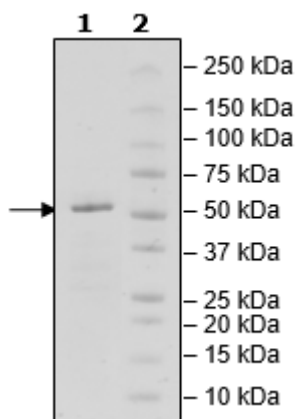


Product Information

Description:	Human Bromodomain containing 4 (also known as BRD4, HUNK1, or MCAP), encompassing amino acids 49-460 corresponding to bromodomains 1 and 2. This construct does not contain a tag.
Species	Human
Construct:	BRD4 (BD1+BD2) (49-460)
Concentration:	0.20 mg/ml
Expression System:	<i>E. coli</i>
Purity:	85%
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:	47 kDa
Genbank Accession:	NM_058243
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 performed according to BRD4 (BD1+BD2) TR-FRET Assay Kit (BPS Bioscience #32612) with BRD4 (BD1+BD2) Recombinant titrated at various concentrations as a competitive antagonist to BRD4 (BD1+BD2) GST-Tag Recombinant (BPS Bioscience #31044).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



BRD4 (BD1+BD2) Activity

