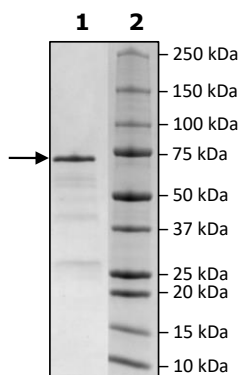


Product Information

Construct:	BRD3 (BD1+BD2) (GST-29-417)
Concentration:	0.52 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_007371
MW:	70 kDa
Purity:	77%
Assay Conditions:	Assay was done according to BRD3 (BD1+BD2) TR-FRET Assay Kit (BPS Bioscience #32615) with various amounts of BRD3 (BD1+BD2), GST-tag.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



BRD3 (BD1+BD2) Activity

