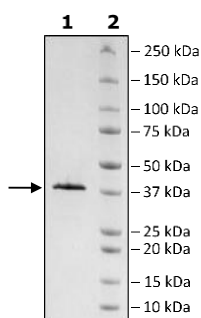


## Product Information

<b>Construct:</b>	BRD2 (GST-65-187)
<b>Concentration:</b>	0.78 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_005104
<b>MW:</b>	41 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was done according to the BRD2 (BD2) Inhibitor Screening Assay Kit (BPS Bioscience #32522) with BRD2, GST-Tag titrated from 1000 ng/rxn – 15.6 ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## BRD2 Activity

