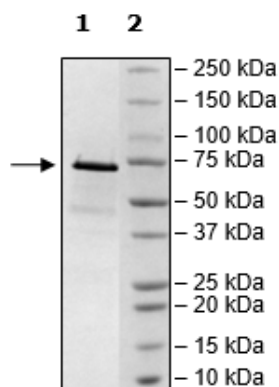


Product Information

| | |
|---------------------------|---|
| Construct: | BRD4 (BD1+BD2) (GST-49-460) |
| Concentration: | 0.32 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, and 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_058243 |
| MW: | 73 kDa |
| Purity: | ≥90% |
| Assay Conditions: | Assay was performed according to BRD4 (BD1+BD2) TR-FRET Assay Kit (BPS Bioscience #32612) with BRD4 (BD1+BD2) (BPS Bioscience #31044) titrated from 72 ng/rxn – 1.125 ng/rxn. |
| 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

