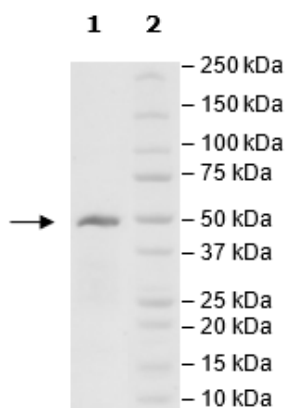


Product Information

Construct:	BRD4 (BD1+BD2) (His-49-460)
Concentration:	1.49 mg/ml
Species:	Human
Formulated In:	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, and 10% 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:	47 kDa
Purity:	85%
Assay Conditions:	BRD4 (BD1+BD2), His-tag (BPS Bioscience #31045) protein is incubated with biotinylated acetyl-histone peptide substrate at 50 nM (BPS Bioscience #33000) in 3x BRD Homogeneous Assay Buffer 1 (BPS Bioscience #33001) in a 10 µl reaction for 1 hour at RT. 20 µl of 1:1 mixture AlphaLISA® Nickle Chelate acceptor beads (PerkinElmer AL108C) diluted 1:250 fold and Alphascreen® Streptavidin -conjugated donor beads (PerkinElmer 6760002S) diluted 1:125 fold was added and incubated for 1 hour at room temperature. After 1 hour incubation at RT, Alphacounts are measured.
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

