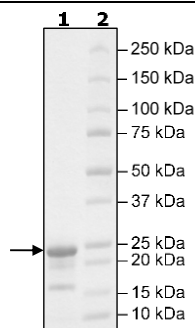


Product Information

Construct:	KRAS (His-2-185)-(BODIPY-GDP)
Concentration:	0.55 mg/ml
Species:	Human
Formulated In:	20 mM HEPES, pH 7.4, 150 mM NaCl and 1 mM DTT
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_033360
MW:	23 kDa
Purity:	72%
Assay Conditions:	Assay was performed by following the KRAS Nucleotide Exchange Assay Kit (BPS Bioscience, #79859) protocols with BODIPY-GDP-loaded KRAS(WT) concentrated at 0.12 mg/ml and GTP titrated 3-fold from 100 μ M – 0 μ M.
Applications:	Useful for KRAS(WT) GDP/GTP nucleotide exchange reaction.
Notes:	KRAS (WT) His-Tag (BPS Bioscience, #11308), is loaded with BODIPY-GDP. Unbound BODIPY-GDP was removed by spin column and it is ready for the GDP/GTP exchange assay.

Quality Control Data

4-20% SDS-Page Coomassie Staining



KRAS BODIPY-GDP Activity

