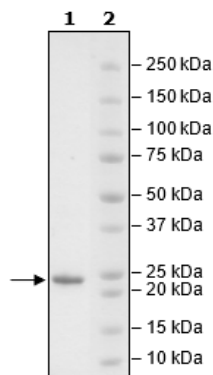


Product Information

Description:	KRAS, His-Tag (BPS Bioscience #11308), is loaded with GDP. Unbound GDP was removed by spin column. Ready for use in the GDP/GTP exchange assay or as a negative control for KRAS-RAF interaction
Construct:	KRAS (His-2-185) – (GDP)
Concentration:	1.35 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_004985
MW:	22 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed according to KRAS (G12C) Coupled Nucleotide Exchange Assay Kit (BPS Bioscience #78004) with KRAS, Isoform B, His-Tag, GDP-loaded Recombinant (BPS Bioscience #101522) titrated from at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



KRAS Activity

