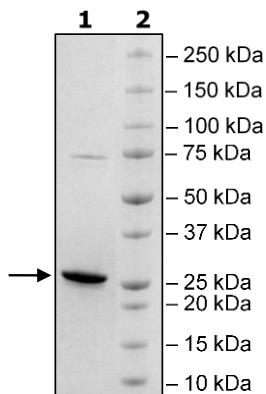


Product Information

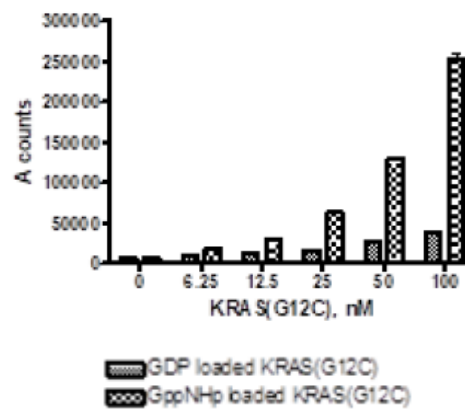
Description:	KRAS (G12C) His-Tag (BPS Bioscience #100413), is loaded with GDP. Unbound GDP was removed by spin column. Ready for use in the GDP/GTP exchange assay or as a negative control for KRAS-RAF interaction.
Construct:	KRAS (G12C) (His-TEV-2-186(end)) - (GDP)
Mutation:	G12C
Concentration:	0.38 mg/ml
Species:	Human
Formulated In:	20 mM HEPES, pH 7.4, 150 mM NaCl, and 1 mM DTT
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_033360
MW:	23 kDa + GDP
Purity:	≥90%
Assay Conditions:	KRAS (G12C) GDP-loaded was tested for binding to the RBD of RAF1. KRAS (G12C) GDP-loaded was compared with KRAS (G12C) GppNHp-loaded (BPS Bioscience #100641) for RBD-RAF1 (BPS Bioscience #100519) binding by Bio-layer Interferometry (GATOR, Probe Life) and AlphaScreening (PerkinElmer).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



RBD-RAF1 (2.5 nM)/KRAS (G12C)



Bio-layer Interferometry (GATOR)

