

## Product Information

<b>Description:</b>	KRAS (G12V), Isoform B, His-Tag (BPS Bioscience #100480), 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.
<b>Construct:</b>	KRAS (G12V) (His-2-185) – (GDP)
<b>Mutation:</b>	G12V
<b>Concentration:</b>	1.43 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	20 mM HEPES, pH 7.4, 150 mM NaCl, 1 mM DTT
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_004985
<b>MW:</b>	22 kDa
<b>Purity:</b>	≥90%

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

