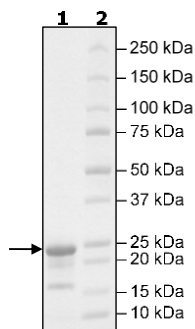


## Product Information

<b>Description:</b>	KRAS, 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.
<b>Label:</b>	BODIPY™ FL GDP ((ThermoFisher Scientific, Guanosine 5'-Diphosphate, BODIPY™ FL 2'-(or-3')-O-(N-(2-Aminoethyl) Urethane), Bis (Triethylammonium) Salt), is a mixed isomer in which the BODIPY® FL fluorophore has been attached to the 2' or 3' position of the ribose ring via a linker. Invitrogen BODIPY® FL dye is a green-fluorescent dye with similar excitation and emission to fluorescein (FITC) or Alexa Fluor 488. It is characterized by a high extinction coefficient and high quantum yield and is relatively insensitive to pH changes. Its hydrophobic properties make it useful with lipophilic compounds. The dye has an excited-state lifetime of 5 nanoseconds or longer, making it useful for fluorescence polarization-based assays.
<b>Construct:</b>	KRAS (His-2-185)-(BODIPY-GDP)
<b>Concentration:</b>	0.55 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	20 mM HEPES, pH 7.4, 150 mM NaCl and 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>	72%
<b>Assay Conditions:</b>	Assay was performed by following the KRAS Nucleotide Exchange Assay Kit (BPS Bioscience #79859) protocols with BODIPY-GDP-loaded KRAS concentrated at 0.12 mg/ml and GTP titrated 3-fold from 100 μM – 0 μM.
<b>Applications:</b>	Useful for KRAS(WT) GDP/GTP nucleotide exchange reaction.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



KRAS BODIPY-GDP Activity

