

## Product Information

|                           |   |
|---------------------------|---|
| <b>Description:</b>       | KRAS (G12V), Isoform B, His-Tag Recombinant (BPS Bioscience, #100480), 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 (G12V) (His-2-185)-(BODIPY-GDP)  |
| <b>Mutation:</b>          | G12V  |
| <b>Concentration:</b>     | 0.42 mg/ml  |
| <b>Species:</b>           | Human   |
| <b>Formulated In:</b>     | 20 mM HEPES, pH 7.4, 150 mM NaCl, and 1.0 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

