

## Product Information

<b>Description:</b>	Recombinant human KRAS (GTPase), isoform a, encompassing amino acids 2-186(end). This construct contains an N-terminal His-tag (6xHis) followed by an optimal TEV protease target sequence that allows cleavage of the His-tag from the protein (TEV: tobacco etch virus cysteine protease BPS Bioscience #50308) and mutation of interest (G12C). The recombinant protein was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	KRAS (G12C) (His-TEV-2-186)
<b>Mutation:</b>	G12C
<b>Concentration:</b>	1.27 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	20 mM HEPES, pH 7.4, 150 mM NaCl, 1 mM DTT, and 10% Glycerol
<b>MW:</b>	23 kDa
<b>Genbank Accession:</b>	NM_033360
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

## Quality Control Data

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

