

Product Information

Description:	Recombinant human KRAS (GTPase), isoform a, encompassing amino acids 2-186 (end). In this construct the N-terminal end of the protein contains a 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). The protein also contains mutation of interest G12D. The protein was affinity purified.
Species	Human
Construct:	KRAS (G12D) (His-TEV-2-186(end))
Mutation:	G12D
Concentration:	1.39 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	20 mM HEPES, pH 7.4, 150 mM NaCl, 10% glycerol, and 1 mM DTT
MW:	23 kDa
Genbank Accession:	NM_033360
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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

