

Product Information

Construct:	KRAS (G12R) (His-2-185)
Mutation:	G12R
Concentration:	3.47 mg/ml
Species:	Human
Formulated In:	20 mM HEPES pH 7.4, 150 mM NaCl, 1 mM DTT, and 10% glycerol.
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_004985
MW:	22 kDa
Purity:	>90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

