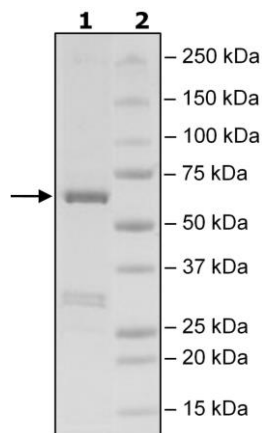


Product Information

Construct:	BRAF (V600E) (GST-His-417-766(end))
Mutation:	V600E
Concentration:	0.17 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
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_004333
MW:	69 kDa
Purity:	79%
Specific Activity:	166 pmol/min/g
Assay Conditions:	Assay was done in Kinase buffer containing 1 mM DTT, 0.1 mg/ml Raf substrate (Mek1 BPS Bioscience #40075 Lot 190508-3), and 10 μ M ATP. Reaction was done at 30°C for 45 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



BRAF Activity

