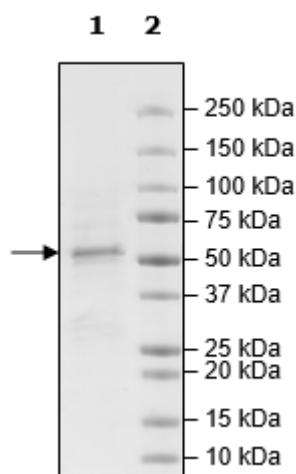


Product Information

Construct:	aRAF (GST-282-606(end)-His)
Mutations:	YY301, 302DD
Concentration:	0.08 mg/ml
Species:	Human
Formulated In:	40mM Tris, pH8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween 20, 3mM DTT, and 20% 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_001654
MW:	63 kDa
Purity:	61%
Specific Activity:	6 pmol/min/μg
Assay Conditions:	Assay was done according to the BRAF (V600E) Kinase Assay Kit (BPS Bioscience #48688) with aRAF titrated from 800 ng/rxn – 12.5 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

