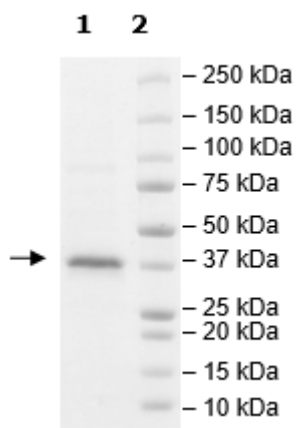


Product Information

Construct:	IRAK1 (His-186-530)
Concentration:	0.39 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 300 mM imidazole, 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_001569
MW:	40 kDa
Purity:	≥90%
Specific Activity:	≥40 pmol/min/μg
Assay Conditions:	Assay was done in Kinase buffer containing 1 mM DTT using MBP substrate (0.2 mg/ml) and 20 μM 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



IRAK1 Activity

