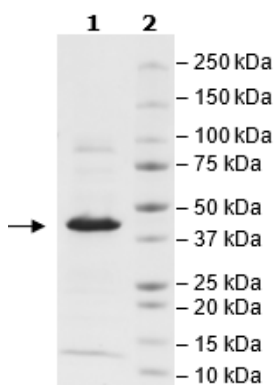


Product Information

Construct:	TDO (His-2-406(end))
Concentration:	0.35 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, 250 mM imidazole
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_005651
MW:	49 kDa
Purity:	85%
Assay Conditions:	Assay was performed according to the TDO Inhibitor Screening Assay Kit (BPS Bioscience #72023) with TDO enzyme titrated from 2,000 ng/rxn – 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



TDO Activity

