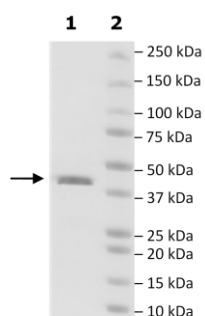


Product Information

Construct:	IDO2 (His-15-420(end))
Concentration:	3.68 mg/ml
Species:	Human
Formulated In:	20 mM Tris-HCl pH 7.5, 250 mM NaCl, 20% glycerol
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_194294
MW:	46 kDa
Purity:	≥90%
Assay Conditions:	IDO2 activity was confirmed using the IDO2 Inhibitor Screening Assay Kit (BPS Bioscience, #72022). Reaction mixture was incubated for 2 hours at 30°C, and absorbance was measured at 320 nm.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



IDO2 Activity

