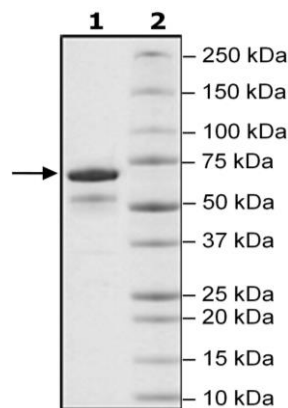


Product Information

Description:	Human Peptidyl-prolyl cis-trans isomerase D (PPID), also known as peptidyl-prolyl cis-trans isomerase Cyclophilin-40 (CYP-40, CyP-D), corresponding to amino acids 2-370(end). The construct contains an N-terminal GST-tag. The recombinant protein was affinity purified.
Species	Human
Construct:	PPID (GST -2-370(end))
Concentration:	1.09 mg/ml
Expression System:	<i>E. coli</i>
Purity:	77%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 16 mM glutathione
MW:	68 kDa
Genbank Accession:	NM_005038
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was performed according to HSP90α C-Terminal Domain TR-FRET Assay Kit (BPS Bioscience #50289) with HSP90α titrated from 48 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



HSP90α: PPID Activity

