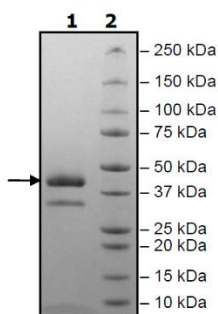


Product Information

Construct:	Renin (67-406-His)
Concentration:	2.00 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 250 mM imidazole
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_000537
MW:	39 kDa
Purity:	80%
Specific Activity:	255 pmol/min/μg
Assay Conditions:	Assay was done according to the Renin Assay Kit (BPS Bioscience #80211) with Renin titrated from 8 ng/rxn -0.125 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

