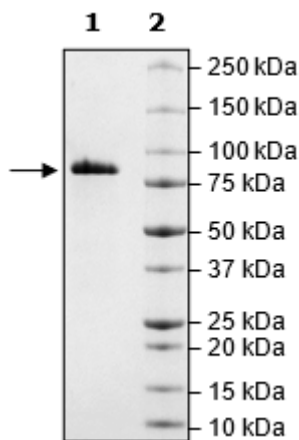


Product Information

Description:	Human Fibroblast Activation Protein also known as Integral Membrane Serine Protease) encompassing amino acids 29-760(end), This construct contains a C-terminal His-tag (6xHis). The recombinant protein was affinity purified.
Species:	Human
Construct:	FAP (29-760(end)-His)
Concentration:	0.89 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 mM Imidazole, 3 mM DTT, and 20% glycerol
MW:	86 kDa
Genbank Accession:	NM_004460
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.
Specific Activity:	≥83 pmol/min/μg
Assay Conditions:	Assay was done according to FAP Assay Kit (BPS Bioscience #80210) with various amounts of FAP.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



FAP Activity

