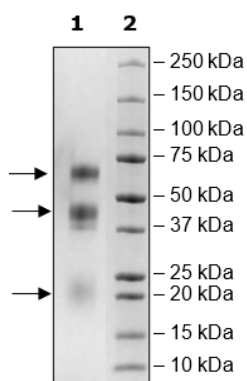


## Product Information

<b>Description:</b>	Recombinant human activated protein C (PROC), full length encompassing 1 – 461(end). This construct contains a C-terminal His-tag (10xHis). This recombinant protein was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	Activated Protein C (PROC) (1-461(end)-His)
<b>Concentration:</b>	0.52 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM Tris-HCl, pH 8.0, 138 mM NaCl, and 2.7 mM KCl
<b>MW:</b>	Reduced conditions: Heavy Chain: 60 kDa; Light Chain: 46 kDa; Propetide: 24 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_000312
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	Assay was done in 50 mM Tris-HCl, pH 8.8, 150 mM NaCl, 10 mM CaCl <sub>2</sub> 0.01% Brij35. Reactions were done in 50 µl duplicates with 100 µM Boc-VPR-AMC substrate (R&D ES011). PROC was titrated at various concentrations. Reaction was incubated for 10 minutes at room temperature. Fluorescence intensity was measured at Excitation 380 nm/ Emission 460 nm.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PROC Activity

