

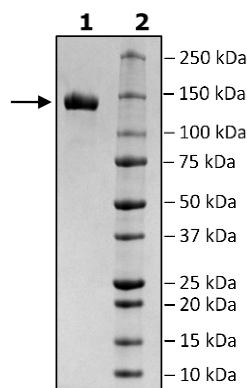
ACE2, Fc Fusion (IgG1), Avi-Tag

Product Information

Construct:	ACE2 (18-740-Fc(IgG1)-Avi)
Concentration:	1.76 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
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_021804
MW:	112 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Assay was done according to the ACE2 Inhibitor Screening Assay Kit (BPS Bioscience #79923) with various amounts of ACE2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



ACE2 Activity

