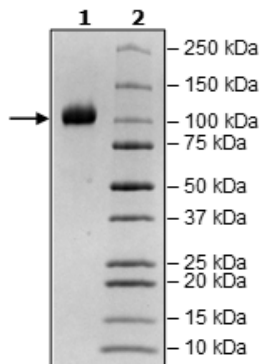


Product Information

Construct:	ACE2 (18-740-His)-(trFluor-Eu)
Label:	trFluor™ Eu (Ex/Em = 346nm / 617 nm) is a fluorophore with large Stokes shifts and an extremely long emission half-life. It is ideal for applications that require high sensitivity and for time-resolved fluorometry because a delay between excitation and the detection of emission minimizes interference. EU-labeled probes are characterized by a relatively high stability and are often insensitive to quenching. trFluor™ Eu is covalently and randomly conjugated to the target protein.
Concentration:	0.55 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:	85 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Assay was performed according to the Spike S1-Biotin (SARS-CoV-2): ACE2 TR-FRET Assay Kit (BPS Bioscience #79949-1) with Spike S1-Biotin (BPS Bioscience #100679) enzyme titrated from 200 ng/rxn – 16 ng/rxn and ACE-Eu at 12.5 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



ACE2-Eu Activity

