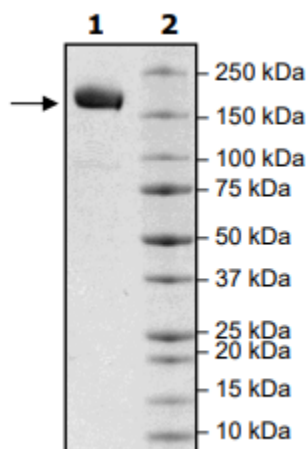


## Product Information

<b>Construct:</b>	Spike Trimer (S1+S2)(SARS-CoV-2) (16-1213-His)-( trFluor-Eu)
<b>Label:</b>	trFluor™ Eu (Ex/Em = 346nm / 617 nm) is a fluorophore with large Stokes shifts and an extremely long emission half-live. 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.
<b>Concentration:</b>	1.35 mg/ml
<b>Species:</b>	SARS-CoV-2
<b>Formulated In:</b>	40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole and 20% glycerol.
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	MN90894
<b>MW:</b>	139 kDa + glycans
<b>Glycosylation:</b>	This protein runs at higher M.W. by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	SARS-CoV2-Trimer Eu-labeled protein (6.25 nM) was incubated with ACE2, His-Avi-Tag, Biotin-labeled HiP (BPS Bioscience, #100665) and the Dye labeled acceptor in ACE2-Spike TR-FRET buffer for 1 hour at room temperature. TRFRET signals were measured and the ratio was calculated as $[\text{Ex}340\text{nm}/\text{Em}665\text{nm}] / [\text{Ex}340\text{nm}/\text{Em}620\text{nm}]$ .
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



SARS-CoV-2 Spike Trimer (Eu-labeled)

