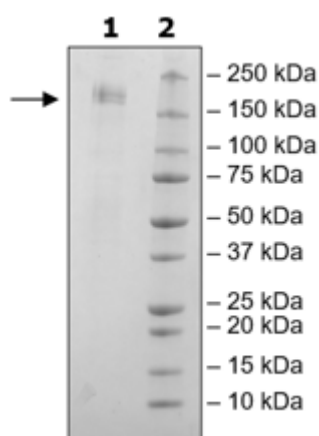


Product Information

Construct:	Spike Trimer (S1+S2) (B.1.1.529 BA.1 Variant) (16-1213-His) (SARS-CoV-2)
Mutations:	A67V, Δ69-70, T95I, G142D, Δ143-145, Δ211, L212I, Ins214EPE, G339D, S371L, S373P, S375F, K417N, N440K, G446S, S477N, T478K, K484A, Q493R, G496S, Q498R, N501Y, Y505H, T547K, D614G, H655Y, N679K, P681H, N764K, D796Y, N856K, Q954H, N969K, L981F
Concentration:	0.48 mg/ml
Species:	SARS-CoV-2
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:	MN908947
MW:	137 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	80%
Assay Conditions:	Assay was done according to the Spike Trimer (S1+S2) (B.1.1.529 BA.1, Omicron Variant) (SARS-CoV-2): ACE2 Inhibitor Screening Colorimetric Assay Kit (BPS Bioscience #78365) with serial dilutions of ACE2, His-Avi-Tag, Biotinlabeled HiP™ (BPS Bioscience #100665). The dilutions ranged from 0 to 1 μM in 3-fold dilutions.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Spike Trimer (S1+S2) (B.1.1.529 BA.1, Omicron Variant) Activity

