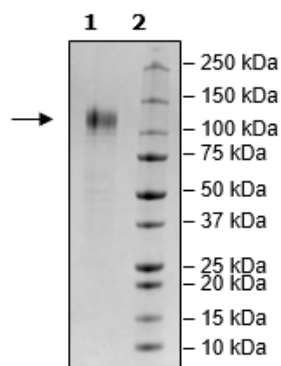


Product Information

Construct:	Spike S1 (B.1.1.529 BA.1 Variant) (16-685-Avi-His) - (Biotin) (SARS-CoV-2)
Mutations:	A67V, Δ 69-70, T95I, G142D, Δ 143-145, Δ 211, I212I, Ins214EPE, G339D, S371L, S373P, S375F, K417N, N440K, G446S, S477N, T478K, E484A, Q493R, G496S, Q498R, N501Y, Y505H, T547K, D614G, H655Y, N679K, P681H
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$.
Concentration:	1.13 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:	78 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	$\geq 90\%$
Assay Conditions:	Assay run according to the Spike S1 (B.1.1.529 BA.1, Omicron Variant) (SARS-CoV-2): ACE2 TR-FRET Assay Kit (BPS Bioscience #78358). 5 ng of ACE2-Eu, 400-fold diluted Dye-labeled acceptor, and various amounts of Spike S1[B] incubated for 1 hour at room temperature. TR-FRET measured at $\text{ex}340 \pm 20\text{nm}$, $\text{em}620 \pm 10\text{nm}$; and $\text{ex}340 \pm 20\text{nm}$, $\text{em}665 \pm 10\text{nm}$.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



ACE2-Eu: Spike S1 [Biotinylated] (Omicron Variant) Binding

