

# Spike Trimer (S1+S2) (XBF, Omicron Variant), His-Tag (SARS-CoV-2) Recombinant

Catalog: 101702

Lot: 230215

## Product Information

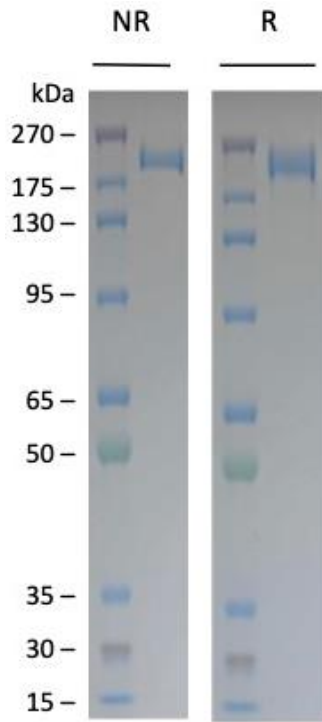
<b>Description:</b>	Recombinant SARS-CoV-2 Spike protein in its homotrimeric form, containing S1+S2 subunits. This protein corresponds to SARS-CoV-2 Omicron Variant XBF and contains the Omicron Spike mutations listed below. The construct also contains a C-terminal His-tag. The recombinant protein was affinity purified.
<b>Species:</b>	SARS-CoV-2
<b>Mutation:</b>	T19I, LPP24-26del, A27S, G142D, K147E, W152R, F157L, I210V, V213G, G257S, G339H, R346T, S371F, S374P, S375F, T376A, D405N, R408S, K417N, N440K, G446S, N460K, S477N, T478K, E484A, F486P, F490S, Q498R, N501Y, Y505H, D614G, H655Y, N679K, P681H, N764K, D796Y, Q954H, N969K
<b>Concentration:</b>	0.50 mg/ml
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	1X PBS, 0.09% NaN <sub>3</sub>
<b>MW:</b>	Apparent MW of >175 kDa in SDS-PAGE due to glycosylation
<b>Stability:</b>	At least 6 months at -80°C. This protein is highly sensitive to freeze/thaw cycles. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	The ELISA was performed by coating 100 µl of ACE2-Fc at 2 µg/ml in PBS at 4°C overnight. The plate was washed in PBS and blocked with 200 µl of 1% BSA in PBS for 1 hour at room temperature. Increasing amounts of Spike Trimer XBF were added to the plate (serially diluted in blocking buffer). The plate was washed in PBS before adding 100 µl of AffiniPure Peroxidase-conjugated Anti-His-Tag antibody (Jackson Immuno Research #300-035-240) diluted 1:5,000 in blocking buffer for 1 hour at room temperature. The plate was washed with PBS and 100 µl of MB/E Ultra-Sensitiv HRP Substrate (EMD Millipore #ES022-500ml) was added for 5 minutes at room temperature. The reaction was stopped with 100 µl of 0.6N H <sub>2</sub> SO <sub>4</sub> and read at 450 nm.
<b>Applications:</b>	Screen or titrate anti-spike antibodies or inhibitors; use as antigen control in assay design and optimization.

# Spike Trimer (S1+S2) (XBF, Omicron Variant), His-Tag (SARS-CoV-2) Recombinant

Catalog: 101702  
Lot: 230215

## Quality Control Data

### Coomassie Staining



### Binding Assay

