

Product Information

| | |
|---------------------------|--|
| Construct: | Spike S1 RBD (K417T, E484K, N501Y) (SARS-CoV-2) (319-541-Avi-His) |
| Mutation: | K417T, E484K, N501Y |
| Concentration: | 2.66 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: | 28 kDa +glycans |
| Glycosylation: | This protein runs at higher M.W. by SDS-PAGE due to glycosylation. |
| Purity: | >90% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

