

Product Information

| | |
|---------------------------|--|
| Construct: | Spike Trimer (B.1.351 Variant) (16-1213-His) (SARS-CoV-2) |
| Mutation: | L18F, D80A, D215G, R246I, K417N, E484K, N501Y, D614G, A701V |
| Concentration: | 0.51 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: | 136 kDa |
| Purity: | ≥90% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

