

Data Sheet

2019-nCoV Spike RBD, Fc-tag

SARS-CoV-2, Recombinant, C-terminal Fc-tag
Catalog #: 100699
Lot #: 200331 **Conc.:** 0.10 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 90 mM Glycine, 01 mM PMSF

Stability: At least 6 months at -80°C. *Avoid freeze/thaw cycles.*

References:

1. Zhou, P. *et al.*, *Nature*. **272**: 270-289, 2020.
2. Xiao, X. *et al.*, *Cell Mol Life Sci*. **61(19-20)**: 2428-2430, 2004.

Description: SARS-CoV-2 2019-nCoV Spike protein S1 subunit, receptor binding domain (RBD), also known as SARS-CoV-2 spike RBD, novel coronavirus spike RBD and nCoV spike RBD, GenBank Accession No. QHD43416.1, a.a. 319-541, with C-terminal Fc-tag, expressed in a CHO cell expression system. MW=50 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2019-nCoV Spike
 RBD

Lane 2:
 Protein Marker

Purity: ≥90%

MW: 50 kDa+ glycans

