

DESCRIPTION

<i>Description</i>	Recombinant Human Coronavirus SARS-CoV-2 Spike Protein S1 subunit, Receptor-Binding Domain (RBD), NCBI Accession Number: MN908947
<i>Sequence</i>	Arg319-Phe541
<i>Expression system</i>	CHO cells
<i>Tag</i>	HIS Tag C-Terminus
<i>Purification</i>	Affinity chromatography
<i>Extinction coefficient</i>	33850 M <sup>-1</sup> .cm <sup>-1</sup> Abs 0.1% (=1 g/l) 1.29 assuming all pairs of Cys residues form cystines
<i>Predicted Molecular Weight</i>	26.1 kDa

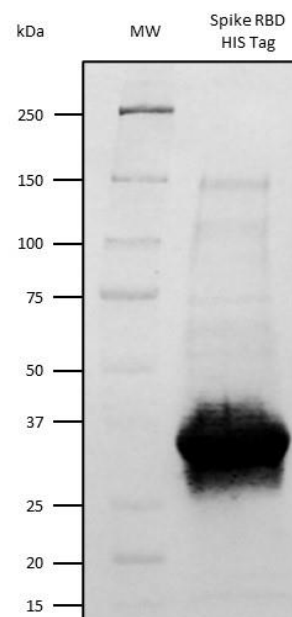
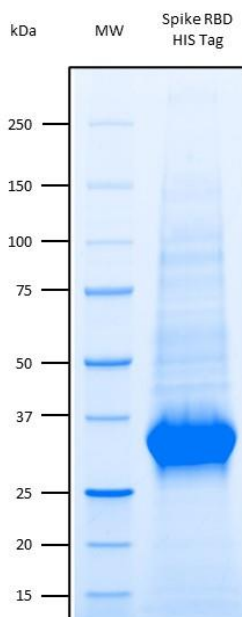
SPECIFICATIONS

<i>SDS Page</i>	Approx. 30 kDa
<i>Concentration</i>	1 mg/ml
<i>Size</i>	100 µg
<i>Purity</i>	>95% by SDS PAGE gel
<i>Formulation</i>	PBS
<i>Endotoxins</i>	Negative on cell-based assay
<i>Activity</i>	Active form: positive binding with recombinant ACE2 Human. Recognized by CR3022 anti RBD recombinant antibody and Covid-19 positive patients serum samples.
<i>Stability and Storage</i>	Store at -20°C. Avoid repeated freeze-thaw cycles Stable one week at 2 to 8°C after reconstitution under sterile conditions Stable 12 months -20°C to -80°C after reconstitution under sterile conditions

DATA

SDS-PAGE 4-15% under reducing conditions and visualized by Coomassie blue staining showing a band at approx. 30 kDa

Western Blot: Spike RBD HIS Tag detected with Anti-HIS-Tag antibody HRP conjugate showing a band at approx. 30 kDa



SARS-CoV-2 Spike Glycoprotein (S1) RBD

Spike protein (S protein) is one of four structural proteins of Coronavirus (SARS-Cov, SARS-Cov-2, MERS amongst other), S protein plays the most important role in viral attachment, fusion and entry, and it serves as a target for development of antibodies, entry inhibitors and vaccines.

In the S protein, the Receptor Binding Domain (RBD) mediates viral entry of SARS-Cov and SARS-Cov-2 C into host cells by its interaction with the membrane receptor ACE2 (angiotensin-converting enzyme 2).

RELATED PRODUCTS

Code	Products	Quantity
715-H16-0BU	Spike RBD His Tag produced in HEK293 cells	100 µg
705-H17-0BU	Nucleoprotein His Tag produced in CHO cells	100 µg
715-H17-0BU	Nucleoprotein His Tag produced in HEK293 cells	100 µg
725-H17-0BU	Nucleoprotein His Tag produced in <i>E. coli</i>	100 µg
857.140.000	Anti S1 Domain Spike protein monoclonal antibody (SARS-CoV-2) – Clone CR3022- Azide free	200 µg/200 µl
857.140.005	Anti S1 Domain Spike protein monoclonal antibody (SARS-CoV-2) – Clone CR3022 Azide free	500 µg/500 µl
857.150.000	Anti Nucleoprotein monoclonal antibody (SARS-CoV-2) – Clone CR3018- Azide free	200 µg/200 µl
857.150.005	Anti Nucleoprotein monoclonal antibody (SARS-CoV-2) – Clone CR3018- Azide free	500 µg/500 µl

Products Manufactured and Distributed by:

Diacclone SAS

6 Rue Dr Jean-François-Xavier Girod

BP 1985, 25020 Besançon Cedex

France

Tel +33 (0)3 81 41 38 38 Fax +33 (0)3 81 41 36 36

Email: [info@diacclone.com](mailto:info@diacclone.com) [www.diacclone.com](http://www.diacclone.com)



**EAGLE BIOSCIENCES**

EAGLEBIO.COM | INFO@EAGLEBIO.COM

P: 617-419-2019 | F: 617-419-1110