

Product Data Sheet

Research use only

Covid19 Spike Peptide library - 96well format - 159 peptides - 1 library

Catalog # : EG-COV19-LIB1

Spike Protein 96 well Overlapping Peptide Library

Coronavirus Spike glycoproteins mediate the virus attachment to host cell surface receptors and facilitate the virus entry by assisting fusion between viral and host cell membranes. They are composed of two subunits, S1 and S2. The Spike protein of SARS-CoV-2 harbors furin cleavage sites, one being located at the boundary between S1 and S2.

Spike glycoproteins are extensively studied as targets for antibodies, vaccine candidates or protease inhibitors, as examples.

We developed a peptide library spanning the entire Spike glycoprotein from SARS-CoV-2. The library is provided in a convenient 96-well plate format adapted for screening purposes.

Spike glycoprotein from SARS-CoV-2: NCBI reference sequence YP_009724390 (1273 aa)

Number of peptides: 159

Peptide length: 15 aa, 9aa for the C-terminal peptide

Peptide sequences: provided on CTDS

Offset number: 7 aa

Counter Ion: TFA

Synthesis scale: 5 µmol

Purity: crude

Gross peptide weight: not quantified

Molecular Mass/ Weight: 1100 to 2072 Da

QC validation: MALDI-TOF QC on 10 % of peptides

Format: lyophilized in 96well plate (P-DW-20-C Axygen®) – 2 plates

Storage Conditions: - 20 °C

For further information please contact our Customer Help Desk

EU : info@eurogentec.com +32 4 372 76 65 – **US :** service@anaspec.com +1 510 791 9560