

Data Sheet

Spike Trimer (S1+S2), His-tag, Eu-labeled (SARS-CoV-2)

SARS-CoV-2, Recombinant, C-terminal
 His-tag, Eu-labeled
Catalog #: 100894
Lot #: 200915 **Conc:** 1.35 mg/ml

Formulated in: 40 mM Tris pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole and 20% glycerol.

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

References:

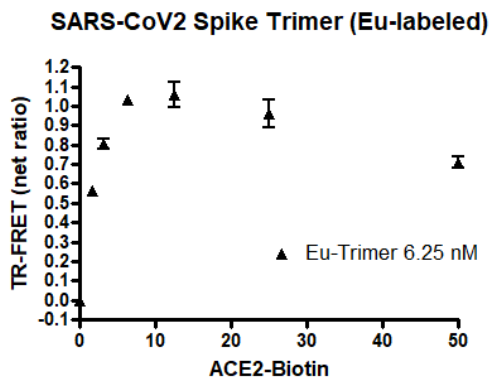
1. Klumpp-Thomas, C., *et al.*, *MedRxiv*. 2020 Jul; preprint doi: [10.1101/2020.07.08.20147371](https://doi.org/10.1101/2020.07.08.20147371)
2. Yurkovetskiy, L., *et al.*, *bioRxiv*. 2020 Jul; preprint doi: [10.1101/2020.07.04.187757](https://doi.org/10.1101/2020.07.04.187757)

Description: Severe acute respiratory Coronavirus Spike trimer (S1+S2), Genbank Accession No. MN908947, a.a. 16-1213, with a with C-terminal His-tag, Eu-labeled, expressed in a HEK293 expression system. MW=139 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Assay Conditions: SARS-CoV2-Trimer Eu-labeled protein (6.25 nM) was incubated with ACE2, His-Avi-Tag, Biotin-labeled HiP (BPS Bioscience, #100665) and the Dye-labeled acceptor in ACE2-Spike TR-FRET buffer for 1 hour at room temperature. TR-FRET signals were measured and the ratio was calculated as [Ex340nm/Em665nm] / [Ex340nm/Em620nm].

Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

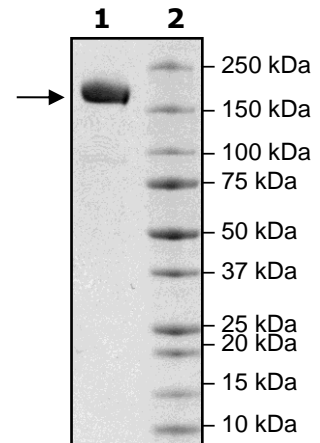


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3µg Spike Trimer
Lane 2:
Protein Marker

MW: 139 kDa +
glycans

Purity: ≥90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com