

## Data Sheet

### **Spike Trimer (S1+S2) Mutant, His-tag (SARS-CoV-2)**

SARS-CoV-2, Recombinant, C-terminal  
His-tag

**Catalog #:** 100728

**Lot #:** 200429

**Conc:** 0.17 mg/ml

**Formulated in:** 8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

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

**Description:** Severe acute respiratory Coronavirus Spike trimer (S1+S2), with 682RRAR685>A, K986P, and V987P mutations, Genbank Accession No. MN908947, a.a. 1-1213, with a C-terminal His-tag, expressed in a HEK293 expression system. MW=139 kDa.

#### **References:**

1. Yan, R. *et al.*, *Science*. 2020 Mar; **367(6485)**:1444-1448.
2. Amanant, F., *et al.*, *MedRxiv*. 2020 Apr; preprint doi: [10.1101/2020.03.17.20037713](https://doi.org/10.1101/2020.03.17.20037713)

---

### **Quality Assurance**

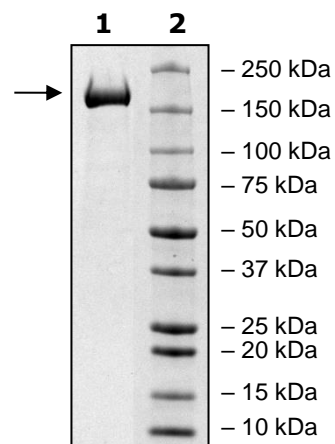
#### **4-20% SDS-PAGE Coomassie staining**

**Lane 1:**  
2µg SARS-CoV-2  
Trimer

**Lane 2:**  
Protein Marker

**MW:** 139 kDa

**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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)