

## Data Sheet

### **Spike Trimer (S1+S2), His-tag (HCoV-NL63)**

HCoV-NL63, Recombinant, C-terminal  
His-tag

**Catalog #:** 100788

**Lot #:** 200610

**Conc:** 1.71 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:** Human Coronavirus NL63 Spike trimer (S1+S2) (HCoV-NL63 S protein), Genbank Accession No. APF29071, a.a. 1-1296(full length), with a C-terminal His-tag, expressed in a HEK293 expression system. MW=146 kDa. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

#### **References:**

1. Arden, K.E., *et al.*, *Med. Virology*. 2005 Mar; **75(3)**: 455-462
2. Vabret, A., *et al.*, *Emerg. Infect. Dis.* 2005 Aug; **11(8)**: 1225-1229.

---

### **Quality Assurance**

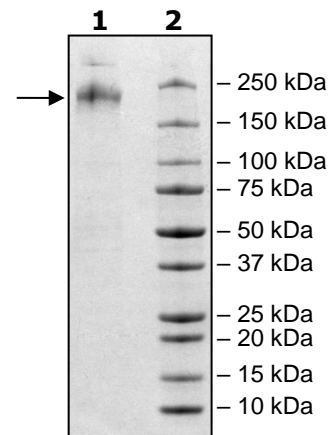
#### **4-20% SDS-PAGE Coomassie staining**

**Lane 1:**  
4µg HCoV NL63 S  
Trimer

**Lane 2:**  
Protein Marker

**MW:** 146 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](mailto:info@bpsbioscience.com)

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