

## Data Sheet

### RNA Polymerase, FLAG-tag (SARS-CoV-2)

SARS-CoV-2, Recombinant, N-terminal  
FLAG-tag

**Catalog #:** 100729

**Lot #:** 200428-1

**Conc:** 0.23 mg/ml

**Formulated in:** 20 mM Tris-HCl pH 8.0, 150 mM NaCl, 4 mM MgCl<sub>2</sub>, 100 µg/ml FLAG, 10% glycerol.

**Description:** Severe acute respiratory Coronavirus 2 RNA dependent RNA polymerase, GenBank Accession No. YP\_00972530, a.a. 1-932(full length), with an N-terminal FLAG-tag expressed in a HEK293 expression system. MW=108 kDa.

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

#### References:

1. Cheng, A. *et al.*, *Virology*. 2005 May; **335(2)**:165-176.
2. Wanchao, Y., *et al.*, *Science*. 2020 May; DOI: 10.1126/science.abc1560

### Quality Assurance

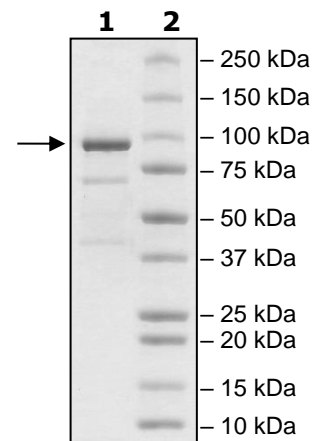
#### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2µg SARS-CoV-2  
RNA polymerase

**Lane 2:**  
Protein Marker

**MW:** 108 kDa

**Purity:** ≥82%



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)