

Data Sheet

NSP7, His-tag (SARS-CoV-2)

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

Catalog #: 100829

Lot #: 200717-2

Conc.: 2.53 mg/ml

Formulated in: 20 mM HEPES pH 7.4, 150 mM NaCl, 4 mM MgCl₂, 250 mM imidazole, 10% glycerol.

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

References:

1. Kirchoderfer, R. N., *et al.*, *Nature Comm.* **10(2342)**: 1-9, 2019.
2. Zhai, Y. *et al.*, *Nature Struct. & Molec. Bio.* **12**: 980-986, 2005.

Description: SARS-CoV-2 nonstructural protein 7 (NSP7), Genbank Accession No.: YP_009725303.1, a.a. 1-84 (full length), with C-terminal His-tag, expressed in an *E. coli* expression system. MW= 10 kDa.

Quality Assurance

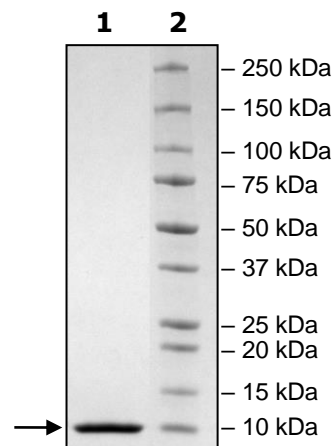
4-20% SDS-PAGE **Coomassie staining**

Lane 1:
2µg SARS-CoV-2
NSP7

Lane 2:
Protein Marker

MW: 10 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

Please visit our website at: www.bpsbioscience.com