

Data Sheet

NSP8, His-tag (SARS-CoV-2)

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

Catalog #: 100830

Lot #: 200717-1

Conc.: 0.54 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 8 (NSP8), Genbank Accession No.: YP_009725304.1, a.a. 1-198(full length), with C-terminal His-tag, expressed in an *E. coli* expression system. MW= 23 kDa.

Quality Assurance

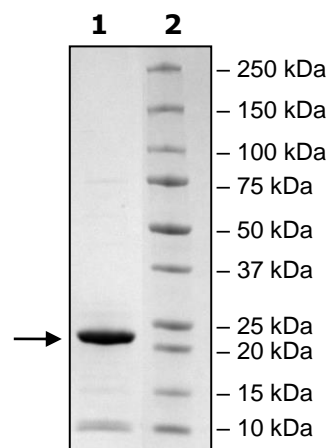
4-20% SDS-PAGE **Coomassie staining**

Lane 1:
2µg SARS-CoV-2
NSP8

Lane 2:
Protein Marker

MW: 23 kDa

Purity: 84%



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