

Data Sheet

NSP10/NSP16 Complex (SARS-CoV-2)

SARS-CoV-2, Recombinant

Catalog #: 100747

Lot #: 200428-1 **Conc.:** 0.12 mg/ml

Formulated in: 20 mM HEPES pH 7.4, 150 mM NaCl, 4 mM MgCl₂, 100 µg/mL FLAG, 10% glycerol.

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

References:

1. Rosas-Lemus, M., *et. al.*, *Europe PMC*. DOI: 10.1101/2020.04.17.047498, 2020.
2. Krafcikova, P. *et. al*, *BioRxiv*. DOI: 10.1101/2020.05.15.097980, 2020.

Description: Complex of SARS-CoV-2 nonstructural protein 10 (**NSP10**), GenBank Accession No. YP_009742617/YP_009725306, a.a. 1-139(full length), with N-terminal FLAG-tag, MW=16 kDa and SARS-CoV-2 nonstructural protein 16 (**NSP16**), Genbank Accession No. YP_009725311, a.a. 1-298(full length), with N-terminal His-tag, MW=34 kDa, co-expressed in a HEK293 cell expression system.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:

4µg SARS-CoV-2

NSP10/NSP16

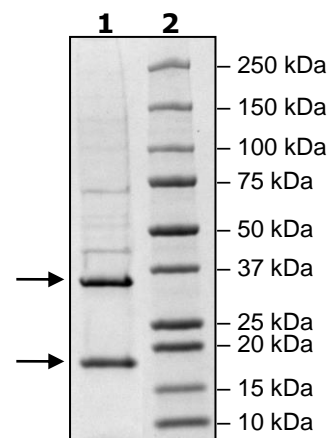
Lane 2:

Protein Marker

MW: NSP16 34 kDa

NSP10 16 kDa

Purity: ≥85%



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