

Data Sheet

Cas12, GST-tag (Lachnospiraceae)

Lachnospiraceae, Recombinant,
N-terminal GST-tag

Catalog #: 100740

Lot #: 200428

Conc: 0.10 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 1 mM DTT 10% glycerol.

Description: Lachnospiraceae Cas12, also known as LpCpf1, Genbank Accession No. ATB19154, a.a 1-1227, with an N-terminal GST-tag, expressed in an *E. coli* expression system. MW=150 kDa.

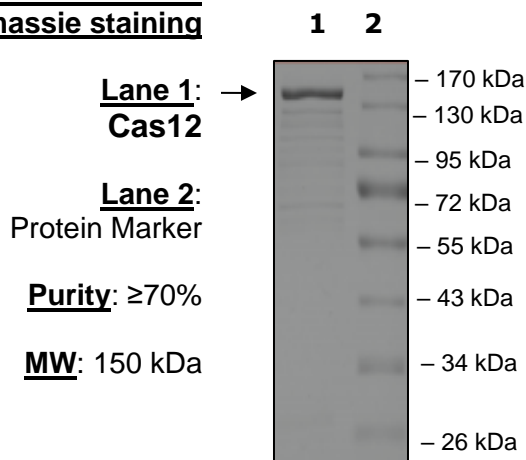
Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Zhou, P., *et al.*, *Nature*. 2020; **579**:270-289.
2. Stefano, S., *et al.*, *Biophysical Journal*. 2020; **118 (3)**: 223A-224A.

Quality Assurance

**10-20% SDS-PAGE
Coomassie staining**



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