



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Cas12, His-tag (Lachnospiraceae)

Lachnospiraceae, Recombinant,
C-terminal His-tag

Catalog #: 100741

Lot #: 200428

Conc: 0.10 mg/ml

Description: Lachnospiraceae Cas12, also known as LpCpf1, Genbank Accession No. ATB19154, a.a 1-1227, with a C-terminal His-tag, expressed in an *E. coli* expression system. MW=130 kDa.

Formulated in: 50 mM sodium phosphate, pH 7.5, 300 mM NaCl, 150 mM imidazole, 1 mM DTT 10% glycerol.

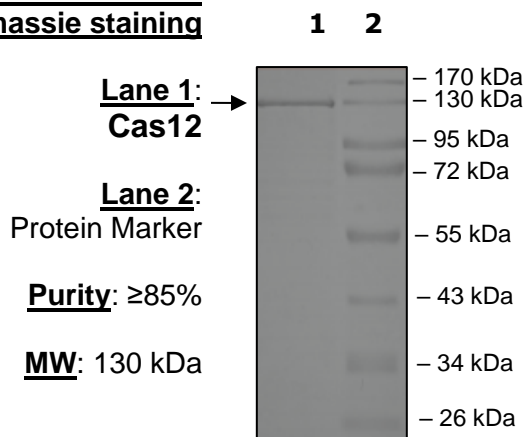
Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

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