



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

PLPro Substrate

Catalog #: 79997

Description: A peptide-AMC substrate for isopeptidase T based on the C-terminal residues of ubiquitin.

Sequence: Z-RLRGG-AMC

Molecular Weight: ~840 Da

Purity: ≥95%

Applications: Useful as a substrate in the Papain-like Protease (SARS-CoV-2) Assay Kit: Protease Activity (BPS Bioscience #79995), as well for screening services.

Storage: -20°C

Optimized for Use With:

Product	Catalog #	Size
Papain-like Protease (SARS-CoV-2) Assay Kit: Protease Activity	79995	96 reactions

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