

Product Information

Description:

The PLPro (Papain-Like Protease) substrate is a small 5-mer fluorogenic peptide, Z-RLRGG-AMC, used to measure the activity of SARS family PLPro proteases. Release of the C-terminal AMC dye (7-amido-4-methylcoumarin) by proteolysis generates fluorescence, which can be monitored at an excitation wavelength of 345 nm and an emission wavelength of 445 nm.

The peptide corresponds to a short sequence of ubiquitin and may also be used as substrate of USP5 (isopeptidase T) and ubiquitin C-terminal hydrolases (UCH family, for example UCHL1).

Sequence:

Z-Arg-Leu-Arg-Gly-Gly-AMC

Molecular Formula:

$C_{40}H_{56}N_{12}O_9$

Supplied As:

DMSO, solution

Formulation:

DMSO

Stability:

At least 6 months at -80°C . Avoid freeze/thaw cycles. Protect from light.

Storage:

-80°C

MW:

~840 Da

Cas No.

167698-69-3

Purity:

$\geq 95\%$

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.

Optimized for Use With:

Papain-like Protease (SARS-CoV-2) Assay Kit: Protease Activity (BPS Bioscience #79995)

Chemical Structure

