

Data Sheet

3CL Protease (SARS-CoV-1 / SARS-CoV-2) Substrate

Catalog #: 79952

Description: Sensitive internally quenched fluorogenic (FRET) substrate for SARS main protease with a K_m value of 17 μM and a k_{cat} value of 1.9 s^{-1} .

Synonyms: DABCYL-Lys-HCoV-SARS Replicase Polyprotein 1ab (3235-3246)-Glu-EDANS trifluoroacetate salt and DABCYL-KTSAVLQSGFRKME-EDANS

Molecular Formula: $\text{C}_{95}\text{H}_{141}\text{N}_{25}\text{O}_{24}\text{S}_2$

Molecular Mass: 2081.45

Purity: >95% by HPLC

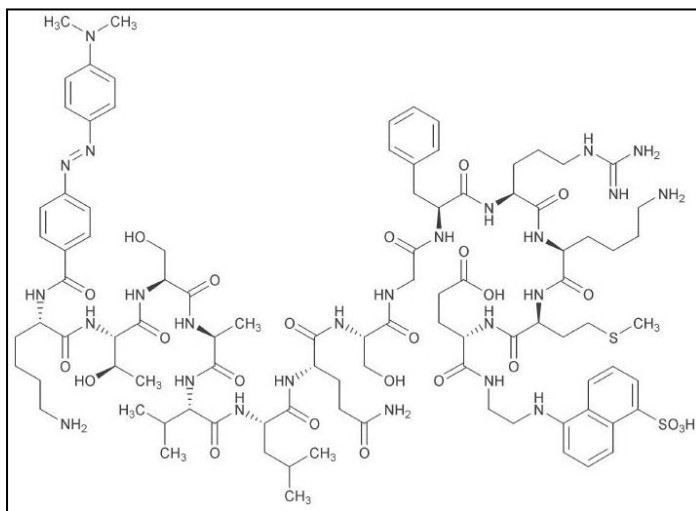
Concentration: 10 mM

CAS Registry No.: 730985-86-1

Supplied as: DMSO, dark red solution

Formulation: DMSO

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



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