

Data Sheet

GC376

Catalog #: 78013

Description: GC376 a 3CL protease inhibitor is optimized for use with the 3CL Protease (SARS-CoV-2) Assay Kit (BPS Bioscience #79955) and 3CL Protease (SARS-CoV) Assay Kit (BPS Bioscience #78015).

Molecular Formula: C₂₁H₃₀NaO₈S

Amount: 50 µg

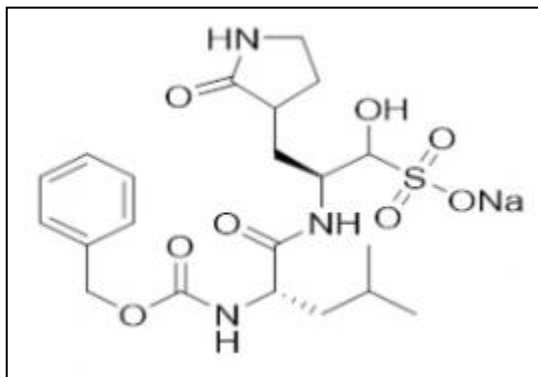
Molecular Mass: 507.5 Da

Supplied as: powder

Solubility: H₂O. To obtain a 500 µM solution, resuspend in 200 µl water.

Applications: Study enzyme kinetics and HTS applications in drug discovery.

Storage: Store at -20°C. Avoid freeze/thaw cycles. Stable at least 12 months from date of receipt, when stored as directed.



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