

Product Information

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:	$C_{21}H_{30}NaO_8S$
Amount:	50 μ g
Molecular Mass:	507.5 Da
Format:	Powder
Solubility:	H ₂ O. To obtain a 500 μ M solution, resuspend in 200 μ l 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.

