

## Product Information

---

<b>Applications:</b>	Study enzyme kinetics, and screen small molecular inhibitors of 3CL MERS-CoV protease for HTS applications in drug discovery.
<b>Supplied As:</b>	Liquid
<b>Format</b>	Aqueous buffer solution
<b>Stability:</b>	Store at -20°C. Avoid freeze/thaw cycles. Stable at least 6 months from date of receipt, when stored as directed.
<b>Storage:</b>	-20°C