

## Product Information

---

<b>Description:</b>	Forster resonance energy transfer substrate peptide for 3CL Protease (MERS-CoV). Optimal excitation wavelength 485 nm, emission wavelength 528 nm.
<b>Concentration:</b>	50 $\mu$ l at 1 mM
<b>Supplied as:</b>	DMSO solution
<b>Size:</b>	50 $\mu$ l
<b>Formulation:</b>	DMSO
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
<b>Storage:</b>	-80°C
<b>Purity:</b>	$\geq$ 95% by HPLC
<b>Applications:</b>	Suitable for studying enzyme kinetics and screen small molecule inhibitors of 3CL Protease (MERS-CoV) for drug discovery and HTS applications.