

Product Information

Description:	Forster resonance energy transfer substrate peptide for 3CL Protease (MERS-CoV). Optimal excitation wavelength 485 nm, emission wavelength 528 nm.
Supplied as:	Powder
Size:	78021-2: 1 mg; 78021-3: 10 mg
Solubility	DMSO
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles. Protect from light.
Storage:	-20°C
MW:	2291.6
Purity:	≥95% by HPLC
Applications:	Suitable for studying enzyme kinetics and screen small molecule inhibitors of 3CL Protease (MERS-CoV) for drug discovery and HTS applications.