

Product Information

Description:	MCL-1 Peptide Ligand is useful as a substrate for screening services and optimized for use with the following TR-FRET assay kits: <ul style="list-style-type: none">• MCL-1 TR-FRET Assay kit (BPS Bioscience #79506)• MCL-1 TR-FRET Assay kit (Mouse) (BPS Bioscience #79928)
Format:	Lyophilized powder
MW:	3162 Da
Purity:	>99%
Sequence:	Biotinylated peptide
Storage/Stability:	-20°C or -80°C when diluted for 6 months. Avoid repeated freeze/thaw cycles.
Solubility:	Aqueous buffers