

## Product Information

---

<b>Description:</b>	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"><li>• MCL-1 TR-FRET Assay kit (BPS Bioscience #79506)</li><li>• MCL-1 TR-FRET Assay kit (Mouse) (BPS Bioscience #79928)</li></ul>
<b>Format:</b>	Lyophilized powder
<b>MW:</b>	3162 Da
<b>Purity:</b>	>99%
<b>Sequence:</b>	Biotinylated peptide
<b>Storage/Stability:</b>	Store at -20°C. Aqueous solutions may be stored in aliquots at -80°C for up to 6 months. Avoid freeze/thaw cycles of solutions.
<b>Solubility:</b>	Aqueous buffers