

## Product Information

---

<b>Description:</b>	Buffer optimized for use in BPS Bioscience METTL Assay Kits.
<b>Applications:</b>	Study enzyme kinetics, screen drugs in high throughput applications.
<b>Supplied As:</b>	Liquid
<b>Format</b>	Tris-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
<b>Optimized for Use With:</b>	METTL3/METTL14 Complex Chemiluminescent Assay Kit (BPS Bioscience #79614)