

## Product Information

---

<b>Description:</b>	This liquid Assay Medium is optimized for use with BPS Bioscience cell-based assay.
<b>Formulation:</b>	<ul style="list-style-type: none"><li>• <b>Assay Medium 1C (BPS Bioscience #78674):</b> MEM medium supplemented with 0.5% charcoal stripped FBS, 1% non-essential amino acids, 1 mM Na pyruvate</li></ul>
<b>Quality:</b>	Sterile filtered (0.2 µm) and Mycoplasma-free. All components are low endotoxin. FBS is ISIA-certified and sourced from the United States of America.
<b>Storage:</b>	Upon receipt, the medium should be stored in the dark at 4°C.
<b>Stability:</b>	Do not exceed 37°C. When stored in the dark at 4°C, the product is stable for six months after first thaw. Avoid repeated freeze/thaw cycles.