

## Product Information

<b>Description:</b>	This liquid Growth Medium is a complete medium optimized to culture select BPS Bioscience cell pools.
<b>Formulation:</b>	<ul style="list-style-type: none"><li>• <b>Growth Medium 1R (BPS Bioscience #78180):</b> Thaw Medium 1 (BPS Bioscience, #60187) plus 0.25 µg/mL Puromycin (Invivogen, #ant-pr-1) to ensure recombinant expression.<ul style="list-style-type: none"><li>○ <b>Thaw Medium 1 (BPS Bioscience #60187):</b> MEM medium (Gibco, #11095-080) supplemented with 10% FBS (Gibco, #26140-079), 1X MEM Non-essential Amino Acids (Corning, #25-025-Cl), 1% Sodium Pyruvate (Corning, #25-000-Cl), and 1% Penicillin/Streptomycin (Gibco, #15140-122).</li></ul></li></ul>
<b>Quality:</b>	Sterile filtered (0.2 µm) and Mycoplasma-free. Each lot is tested for growth of recombinant cells. 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.
<b>Culture Protocol:</b>	For detailed thawing and cell culture protocol, see cell pool data sheet.
<b>Optimized for Use With:</b>	Cas9-Expressing MCF7 Cell Pool (BPS Bioscience, #78179)