

## Product Information

<b>Description:</b>	This liquid Maintenance Medium is a complete medium optimized to maintain BPS Bioscience's induced pluripotent stem cells (iPSC)-derived cardiomyocytes.
<b>Formulation:</b>	<ul style="list-style-type: none"><li>• <b>Maintenance Medium 17 (BPS Bioscience #78509):</b> RPMI 1640 medium supplemented with 5% Cardiomyocyte Support Supplement and 1% Penicillin/Streptomycin.</li></ul>
<b>Quality:</b>	Sterile filtered (0.2 µm) and Mycoplasma-free. All components are low endotoxin.
<b>Storage:</b>	Upon receipt, the medium should be stored in the dark at 4°C.
<b>Stability:</b>	Equilibrate the medium at room temperature before use. Do <b>not</b> pre-warm at 37°C. The medium may be shipped with dry ice and arrive frozen; this does not affect performance. When stored in the dark at 4°C, the product is stable for <b>one month</b> after first thaw. Avoid repeated freeze/thaw cycles.
<b>Culture Protocol:</b>	For detailed thawing and cell maintenance protocols, see cell data sheet.