

Product Information

Description:	This liquid Maintenance Medium is a complete medium optimized to maintain BPS Bioscience's induced pluripotent stem cells (iPSC)-derived cardiomyocytes.
Formulation:	<ul style="list-style-type: none">• Maintenance Medium C17 (BPS Bioscience #78509): RPMI 1640 medium supplemented with 5% Cardiomyocyte Support Supplement and 1% Penicillin/Streptomycin.
Quality:	Sterile filtered (0.2 µm) and Mycoplasma-free. All components are low endotoxin.
Storage:	Upon receipt, the medium should be stored in the dark at 4°C.
Stability:	Equilibrate the medium at room temperature before use. Do not 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 one month after first thaw. Avoid repeated freeze/thaw cycles.
Culture Protocol:	For detailed thawing and cell maintenance protocols, see cell data sheet.