

Product Information

Description:	This liquid Growth Medium is a complete medium optimized to culture select BPS Bioscience cell lines.
Formulation:	<ul style="list-style-type: none">• Growth Medium 2M (BPS Bioscience #78181): Thaw Medium 2 (BPS Bioscience #60184) plus 1 µg/ml of Puromycin (InvivoGen, #ant-pr-1).<ul style="list-style-type: none">○ Thaw Medium 2 (BPS Bioscience #60184): RPMI 1640 medium (Thermo Fisher, Cat. #A1049101) supplemented with 10% FBS (Thermo Fisher, Cat. #26140079), 1% Penicillin/Streptomycin (Hyclone #SV30010.01).
Quality:	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.
Storage:	Upon receipt, the medium should be stored in the dark at 4°C.
Stability:	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.
Culture Protocol:	For detailed thawing and cell culture protocol, see cell line data sheet.