

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 1H (BPS Bioscience #79546): Thaw Medium 1 (BPS Bioscience, #60187) plus 5 µg/ml of Blasticidin (ThermoFisher, #R21001).<ul style="list-style-type: none">○ Thaw Medium 1 (BPS Bioscience #60187): 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).
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.
Optimized for Use With:	T-Rex HEK293 Recombinant Cell Line (BPS Bioscience, #71227)