

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 1S (BPS Bioscience #78221): Thaw Medium 1 (BPS Bioscience # 60187) plus 200 µg/ml of Hygromycin B (Thermo Fisher, #10687010) and 5 µg/ml Blastidicin (Invivogen, #ant-bl-1) to ensure recombinant expression.<ul style="list-style-type: none">○ Thaw Medium 1 (BPS Bioscience #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Thermo Fisher, Cat. #26140079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 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.
Optimized for Use With:	CRISPRa (SAM) HeLa Cell Line (BPS Bioscience, #78193) CRISPRa (SAM) HepG2 Cell Line (BPS Bioscience, #78194)