

Product Information

Description:	This liquid Medium is a complete medium optimized to culture select BPS Bioscience cell lines.
Formulation:	BPS Growth Media contain antibiotics for maintenance of selective markers in stable cell lines. <ul style="list-style-type: none">• Growth Medium 4C (BPS #79631): DMEM/F12 (1:1) supplemented with 10% FBS, 1% Penicillin/Streptomycin. Plus 1 µg/ml Doxycycline and 3 mM Na-butyrate.
Quality:	Sterile filtered (0.2 µm) and Mycoplasma-free. 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.