

Product Information

Description:	This liquid Medium is a complete medium optimized to culture select BPS Bioscience cell lines.
Formulation:	<ul style="list-style-type: none">• Assay Medium 7B (BPS #79718): Opti-MEM I + 0.5% FBS + 1% Non-essential amino acids + 1 mM sodium pyruvate + 1% penicillin/streptomycin.
Quality:	Sterile filtered (0.2 μ m) 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.