

Product Information

Description:	This liquid assay medium is optimized for use in cellular assays in select BPS Bioscience cell lines.
Formulation:	<ul style="list-style-type: none">• Assay Medium 6A (BPS Bioscience #82211): Phenol red-free DMEM, 10% Charcoal Stripped FBS, and 1% Penicillin/Streptomycin.
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.