

Product Information

| | |
|--------------------------|---|
| Description: | This liquid Medium is a complete medium optimized to culture select BPS Bioscience cell lines. |
| Formulation: | BPS Growth contain antibiotics for maintenance of selective markers in stable cell lines. <ul style="list-style-type: none">• Growth Medium 4A (BPS #79535): DMEM/F12 (1:1) supplemented with 10% FBS, 1% Penicillin/Streptomycin. Plus 10 µg/ml Blasticidin, and 200 µg /ml Zeocin. |
| 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. |