

## Product Information

---

<b>Description:</b>	This liquid Cell Thawing Medium is optimized for thawing and plating select BPS Bioscience recombinant cell lines.
<b>Formulation:</b>	<ul style="list-style-type: none"><li>• <b>Thaw Medium 15 (BPS Bioscience #78412):</b> IMDM medium supplemented with 20% FBS, 1% Penicillin/Streptomycin.</li></ul>
<b>Quality:</b>	Sterile filtered (0.2 $\mu\text{m}$ ) and Mycoplasma-free. All components are low endotoxin. FBS is ISIA-certified and sourced from the United States of America.
<b>Storage:</b>	Upon receipt, the medium should be stored in the dark at 4°C.
<b>Stability:</b>	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.
<b>Culture Protocol:</b>	For detailed thawing and cell culture protocol, see cell line data sheet.