

Product Information

- Description:** This liquid Cell Thawing Medium is optimized for use with tetracycline-inducible systems.
- Formulation:**
- **Thaw Medium 16 (BPS Bioscience #78647):** MEM medium supplemented with 10% TET-free FBS, 1% non-essential amino acids, 1 mM Na pyruvate, 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.