



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Thaw Medium 10 Catalog #: 79704

Description

This liquid medium is optimized to culture select BPS Bioscience primary cells and cell lines.

Formulation

RPMI 1640 medium (Thermo Fisher, #A1049101) supplemented with 10% FBS (Thermo Fisher, #26140079), 1mM Sodium pyruvate (Corning, #25-000-CL), 1% Non-essential amino acids (Corning, #25-025-CL), 1% Penicillin/streptomycin (Thermo Fisher, #15140122).

Quality

Sterile filtered (0.2 μ m) and Mycoplasma-free. Each lot is tested for growth of recombinant cells. All components are low endotoxin. FBS is ISIA-certified and sourced from the United States of America.

Storage and Stability

Upon receipt, the medium should be stored in the dark at 4°C. Avoid repeated freeze/thaw cycles. Do not exceed 37°C. When stored in the dark at 4°C, the product is stable for six months after first thaw.

Culture Protocol

For detailed thawing and cell culture protocol, see cell line data sheet.

Optimized for Use With

Product	Cat #	Size
LentiLight GFP Lentivirus	#79703	2 x 1 ml
CD4+ T cells, Negatively Selected (Human)	#79752	10 million cells
CD8+ T cells, Negatively Selected (Human)	#79753	10 million cells
Normal Human Peripheral Blood Mononuclear Cells	#79059	10 or 30 million cells
Firefly Luciferase–RPMI 8226 Recombinant Cell Line	#79834	2 vials
Anti-BCMA CAR Lentivirus	#79701	500 μ l x 2

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com