



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Assay Medium 3A

Catalog #: 79816

Description

This liquid assay medium is optimized for use in cellular assays in select BPS Bioscience cell lines.

Formulation

F-12K Medium (Kaighn's Modification of Ham's F-12 Medium) supplemented with 1% FBS (Thermo Fisher, #26140079), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

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.

Protocol

For detailed cellular assay protocol, see cell line data sheet.

Optimized for Use With

- CD155 / TCR Activator - CHO Recombinant Cell Line #60548
- TIGIT / NFAT Reporter - Jurkat Cell Line #60538

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com