



Certificate of Analysis

iLite[™] Human Type I Interferon Assay Ready Cells

Catalog No: 51100-7

Lot No: 7498

Volume: 2.5 ml/vial

Expiration: March 25, 2023

Instructions on Thawing and Plating for Use in Human Interferon Alpha Assay:

1. Rapidly thaw the vial of cells using a 37°C water bath. Thawing should be accomplished within 3-4 minutes. Gently invert vial a minimum of ten times to ensure a uniform cell suspension. Transfer cells to a sterile reservoir. **Do not vortex cells.**
2. Immediately transfer 25µl of the cell suspension to each well of a sterile, 96-well microtiter plate using a multichannel micropipette with sterile tips. Cells must be used within 30 minutes of thawing.

Shipping Information

Shipping Conditions: Dry Ice

Physical State of Product During Shipping: Frozen

Storage Conditions/Comments: After receipt, store the product at -80°C only for retention of full activity.

For more information on protein handling, visit our Resource Library at www.pbl assaysci.com.

Authorization

Released by: _____

Date: March 2, 2022

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.