



## Certificate of Analysis

### Human Type I IFN Neutralizing Antibody Mixture

**Catalog No:** 39000-1

**Lot No:**

**Expiration:**

**Volume:** 0.2 ml

**Description:** Mixture of monoclonal and polyclonal antibodies directed against human Type I interferon receptor and several Type I interferons.

**Activity:** At a working dilution of 1:50, this mixture will neutralize 50% of the protective effect of the following interferon amounts in an antiviral assay:

<b>Hu-IFN Alpha, 2a</b>	ug/ml
<b>Hu-IFN Beta, 1a</b>	ug/ml
<b>Hu-IFN Omega</b>	ug/ml

**Buffer:** PBS

**Endotoxin:** < 1 EU/ g

**Purification Method:** A combination of Protein G affinity chromatography and other conventional chromatographical methods.

**Specificity:** Blocks biological activity of human Type I interferons (alpha, beta, omega, kappa, epsilon). At this working dilution, no cross reactivity is observed with human interferon gamma, human interferon lambda 1, human interferon lambda 2, human interferon lambda 3, mouse interferon alpha, mouse interferon beta or mouse interferon gamma.

**Assay Used to Measure Bioactivity:** Encephalomyocarditis virus (EMCV) was used to challenge A549 cells [Budd et al. (1985) Canc. Chem. Pharm. 12:39] which were incubated with a constant concentration of the antibody mixture and the titrated interferon.

**Tested Applications:** Neutralization

*Optimal dilutions should be determined by each laboratory for each application*

**Shipping Conditions:** Dry Ice

**Physical State of Product During Shipping:** Frozen

**Storage Conditions/Comments:** After receipt, the product should be kept at -70°C or below for retention of full activity. Thaw product vial by incubation in cold tap water until just thawed . the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA). For more information on protein handling, visit our Resource Library at [www.pbl assaysci.com](http://www.pbl assaysci.com).

### Authorization

Released by: \_\_\_\_\_

Date:

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.