



## Certificate of Analysis

### Anti-Human Interferon Alpha, Rabbit Serum (PAb)

**Catalog No:** 31101-1

**Lot No:** 7357R

**Size:**  $\geq 1 \times 10^5$  Neutralizing Units/vial

**Description:** Rabbit Polyclonal Antibody against Human Interferon Alpha

**Volume:** 0.100 ml

**Activity:**  $4.64 \times 10^6$  NU/ml

**Concentration:** 0.67 mg/ml (estimate)

**Buffer:** Phosphate buffered saline (PBS)

**Antigen:** Recombinant human interferon alpha 2 (alpha 2b) (expressed in *E. coli*)

**Assay Used to Measure Bioactivity:** One neutralization unit is the amount of antiserum required to neutralize one unit of human interferon alpha (Hu-IFN- $\alpha$ ) to a 50% endpoint. Interferon was titrated with the use of the cytopathic effect inhibition assay [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) *J. Virol.* 37, 755-758] using A549 cells and EMCV was used as the challenge virus [Budd *et al.* (1985) *Canc. Chem. Pharm.* 12:39]. In this antiviral assay for interferon, about 1 unit/ml of interferon is the quantity necessary to produce a cytopathic effect of 50%. The units are determined by use of a laboratory standard calibrated with respect to the international reference standard for Hu-IFN- $\alpha$  provided by the National Institutes of Health [Gxa01-901-535]. This material is prepared specifically for effective neutralization of Hu-IFN- $\alpha$

**Tested Applications:** Neutralization

**Suggested Applications:** ELISA; Western blot; immunoprecipitation; immunohistochemistry

*Please note that these applications are presented for suggested use only and have not been fully evaluated by PBL.*

**Product Information:** The most common use of this product has been in a pool with Rabbit Anti-Human Interferon Beta (PBL # 31410) and Mouse Anti-IFNAR2 (PBL# 21385) to block the activity of all Type I IFNs. Krug *et al* [*Eur. J. Immunol.* (2001) 31:2154] used a pool of PBL #31101, #31410 and #21385 to demonstrate PBMC-derived Type I IFN is responsible for increased IFN-gamma secretion after stimulation with CpG ODN. Ito *et al.* [*J. Exp. Med.* (2002) 195:1507] used a similar pool to determine that TLR7 stimulated pDC but not mDC induce Th1 development through Type I IFN. Lande *et al.* [*J. Immunol.* (2003) 170:1174] used a pool of PBL#31410 and #31101 to show a blockade of autocrine Alpha/Beta action leads to decreased CXCL-10 secretion from *Mycobacterium*-infected dendritic cells. Jago *et al.* [*Immunity* (2003) 19:225] demonstrated the inhibition of IFN-alpha signaling in a co-culture of pDCs with CD40 stimulated B-cells reduces IgG secretion, while inhibition of IFN-beta signaling does not occur. Ku *et al.* [*J. Exp. Med.* (2004) 200:917] showed epithelial cells surrounding Varicella Zoster lesions express IFN-alpha by immunohistochemistry (IHC) in a SCID-human skin xenograft model.

**Shipping Conditions:** Dry Ice

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

**Storage Conditions/Comments:** After receipt, the product may be stored at -20°C for short-term use ( $\leq 6$  months). For long-term storage, we recommend storing the product at -70°C or below for retention of full activity. When thawing, 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.pblassaysci.com](http://www.pblassaysci.com).

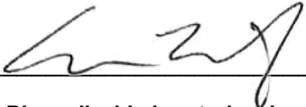


# pbl assay science

131 Ethel Road West, Suite 6 | Piscataway, NJ 08854 USA | T:+1.732.777.9123 | F:+1.732.777.9141 | E: info@pbl assaysci.com | W: pbl assaysci.com

---


## Authorization

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

Date: March 26, 2024

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.

Rev. 05

A decorative footer consisting of a wavy, multi-colored gradient bar transitioning from light green on the left to bright blue on the right.