

Certificate of Analysis

Rat Interferon Alpha, mammalian

Catalog No: 13101-9

Lot No: 7812

Size: 0.1 mg/vial

Description: Recombinant Rat Interferon Alpha (Rat IFN- α)

Concentration: 0.505 mg/ml

Volume: 0.198 ml/vial

Typical Activity (EC₅₀): 20–100 pg/ml

Buffer: 50 mM Tris/Citrate, pH 5.0; 600 mM NaCl; 6% Glycerol

Molecular Weight: 19.0 kDa

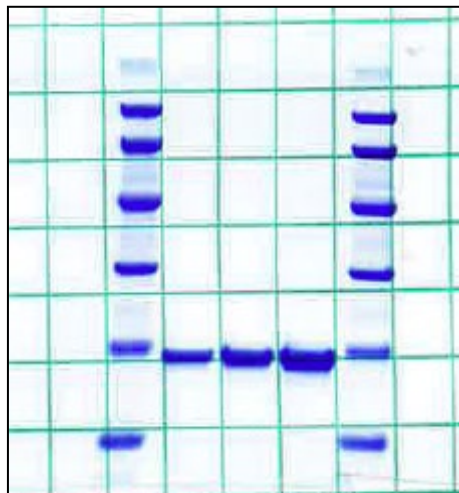
Purity: > 95%

Source: Rat Alpha 1 gene expressed in mammalian cells

Accession #: X00336

Assay Used to Measure Bioactivity: Activity was measured on rat epithelial-like (RATEC) cells challenged with VSV in the presence of 2 mM Glutathione. For full bioactivity, it is recommended that the reducing agent glutathione is added to the culture medium. Interferon was titrated with the use of the cytopathic effect inhibition assay as described [Rubinstein, S., Familletti, P.C., and Pestka, S. (1981) "Convenient Assay for Interferons," *J. Virol.* **37**, 755-758; Familletti, P.C., Rubinstein, S., and Pestka, S. (1981) "A Convenient and Rapid Cytopathic Effect Inhibition Assay for Interferon," in *Methods in Enzymology*, Vol. 78 (S. Pestka, ed.), Academic Press, New York, 387-394].

Image of the Purity Gel:



Lane	Sample
1	Marker
2	13101 - 1 μ g
3	13101 - 2 μ g
4	13101 - 3 μ g
5	Marker

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) or Tissue Culture Media with serum. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____



Date: September 22, 2025

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. 02