



Certificate of Analysis

Anti-Pig Interferon Alpha, Clone F17 (MAb)

Catalog No: 27105-1

Lot No: 7579

Size: 500 µg/vial

Description: Mouse Monoclonal Antibody against Pig Interferon Alpha, neutralizing

Clone: F17

Volume: 0.258 ml

Concentration: 1.94 mg/ml

Buffer: 31 mM Citric Acid, 139 mM Sodium Phosphate Dibasic, 50% Glycerol pH 6.4

Antigen: Pig Interferon Alpha

Isotype: Mouse IgG₁

Endotoxin: < 1 EU/µg

Purity: > 95%

Purification Method: Affinity chromatography via Protein G

Bioactivity: Neutralizes pig interferon alpha

Assay Used to Measure Bioactivity: Neutralization of Pig Interferon Alpha on MDBK cells in cytopathic effect inhibition assay.

Suggested Applications: Western blot; immunoprecipitation; immunohistochemistry; ELISA

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

References:

1. L'Haridon *et al.* (1991) *Hybridoma*, 10: 35-47.
2. Splichal *et al.* (1994) *Immunology Letters*, 43: 203-208. (ELISA)

Shipping Conditions: Wet Ice

Physical State of Product During Shipping: Liquid

Storage Conditions/Comments: After receipt, this product should be kept at -20°C for retention of full activity. Do not store below -20°C. For more information on protein handling, visit our Resource Library at www.pblassaysci.com.

Authorization

Released by: _____

Date: January 12, 2023

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.