Certificate of Analysis

Anti-Pig Interferon Alpha, Clone F17 (MAb)

Catalog No: 27105-1

Lot No:

Size: 500 µg/vial

Description: Mouse Monoclonal Antibody against Pig Interferon Alpha, neutralizing

Clone: F17

Volume: ml

Concentration: mg/ml

Buffer:

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.

below -20 C. For more information on protein harding, visit our resource Library at www.pbiassaysci.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.