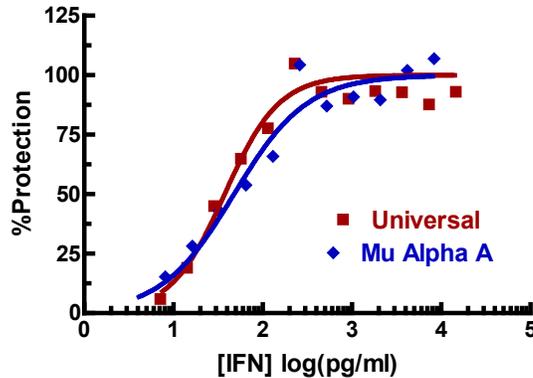




incubation in cold tap water until just thawed – the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein 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).

### Comparison of Universal IFN and Mouse Alpha A Antiviral Activity on Mouse Cells



**Figure 1:** The activity of Mouse IFN Alpha-A and Universal Type-I IFN was compared in an L929/EMCV CPE assay. The EC<sub>50</sub> for Universal IFN in this experiment was 38 pg/ml while the EC<sub>50</sub> for Mu Alpha-A was 48 pg/ml. Similar results were obtained for several batches of Mouse Alpha-A. Results are representative and may vary depending upon experimental conditions.

#### 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.