



Mammalian Cell-Expressed Rat Interferon Alpha Proteins

Catalog No. 13101-1 (*IFN Alpha 1*), 13114-1 (*IFN Alpha 14*)

- Bioactivity measured on rat epithelial-like cells (RATEC)
 - Recommend adding the reducing agent glutathione to culture medium for full activity
- Typical Activity (EC_{50}) is 20-100 and 2-10 $\mu\text{g/ml}$ for IFN Alpha 1 and 14, respectively
- Greater than 95% purity

Figure 1. Mammalian cell-expressed Rat IFN Alpha 1 CPE Assay

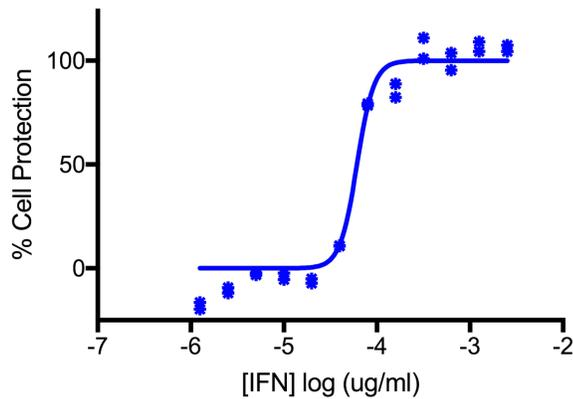


Figure 2. Mammalian cell-expressed Rat IFN Alpha 14 CPE Assay

