



Human Cell-Expressed Human Interferon Lambda Proteins

IFN Lambda 2/IL-28A; IFN Lambda 1/IL-29; IFN Lambda 3/IL-28B

Catalog No. 11720-1, 11725-1, 11730-1

- Activity quantified versus the lambda 1 international standard
- Authentic human cell origin
- Provided in units of activity (100,000 units/vial)

Figure 1. IFN Lambda 2/IL-28A (11720-1)

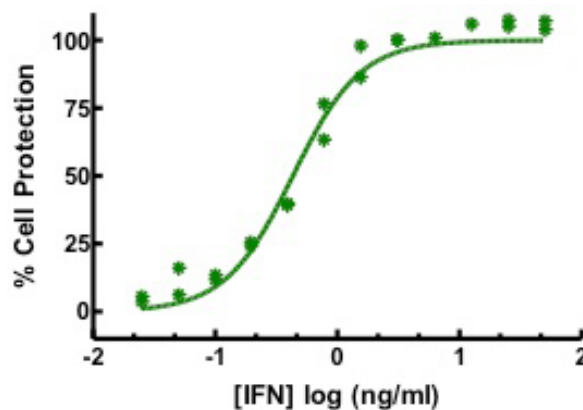


Figure 2. IFN Lambda 1/IL-29 (11725-1)

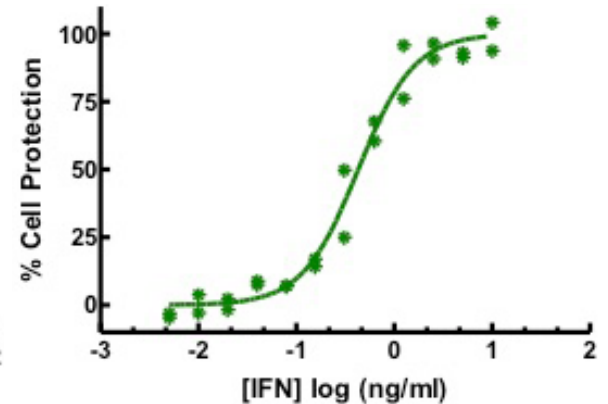
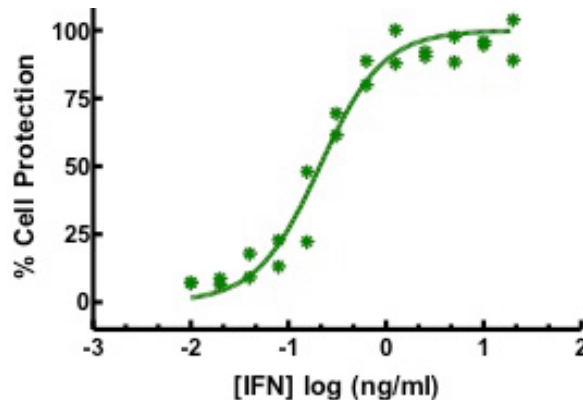


Figure 3. IFN Lambda 3/IL-28B (11730-1)



Activities of Lambda interferons were quantified using the cytopathic effect inhibition assay (CPE; A549/EMCV). The EC_{50} for interferon in this assay is ~ 1 U/ml. The units are determined versus a Human IFN Lambda 1 (Hu-IFN- $\lambda 1$) laboratory standard calibrated to the international reference standard for Human Interferon Lambda 1 (Hu-IFN- $\lambda 1$) 10/176 provided by the National Institute for Biological Standards and Control (NIBSC).