



Universal Type I Interferon

Recombinant Human IFN Alpha hybrid protein expressed in *E. coli*

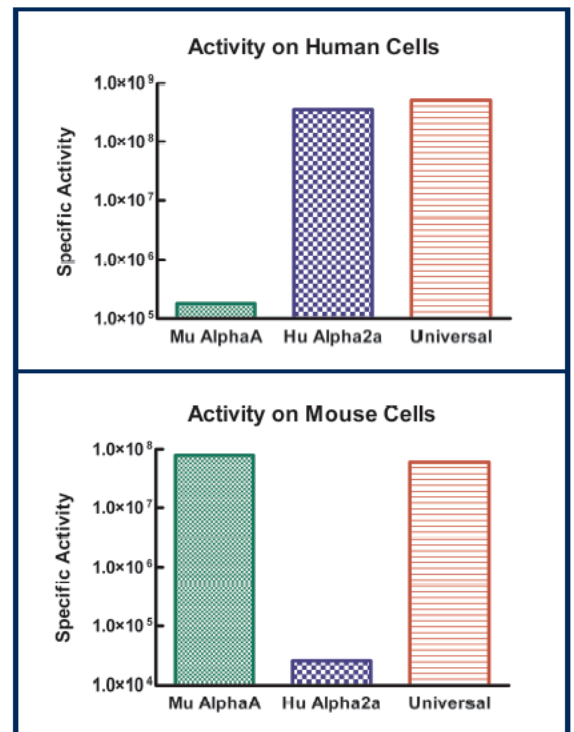
Catalog No. 11200

- Best human IFN protein for broad crossover to multiple animal models, *in vivo* and *in vitro*
- Suitable for animal studies where species-specific IFN alpha is not available
- High specific activity on multiple animal cell lines
- Useful from cross-species comparative studies of Type I IFN effects

Figure 1. Comparison of the Activity of Human IFN Alpha-2a, Mouse IFN Alpha-A, and 'Universal' IFN, on human and mouse cells.

Specific activity of the three IFNs was determined in an A549/EMCV CPE assay (top panel) and a L929/EMCV CPE assay (bottom panel).

Both Universal and Human IFN Alpha-2a are potent on human cells while the Mouse IFN Alpha-A has little activity on these cells. Likewise the Mouse IFN Alpha-A and Universal IFN are potent on mouse cells while the Human IFN Alpha-2a has little activity.



Product Information

Catalog Number	11200-1, -2
Description	Universal Type I Interferon
Size	1 x 10 ⁵ units, 1 x 10 ⁶ units

Product Citations

1. Parainfluenza virus 5 genome are located in viral cytoplasmic bodies whilst the virus dismantles the interferon-induced antiviral state of cells, Carlos *et al.*, (2009), *Journal of Genetic Virology*, 90:2147.
2. Abortive versus Productive Viral Infection of Dendritic Cells with a Paramyxovirus Results in Differential Upregulating of Select Costimulatory Molecules, Alexander-Miller *et al.*, (2005), *Journal of Virology*, 79(12):7544.



For more information on this product, visit our website.