

# Validation data for LumiKine™ Xpress hIFN-α 2.0

<https://www.invivogen.com/lumikine-xpress-hifna>

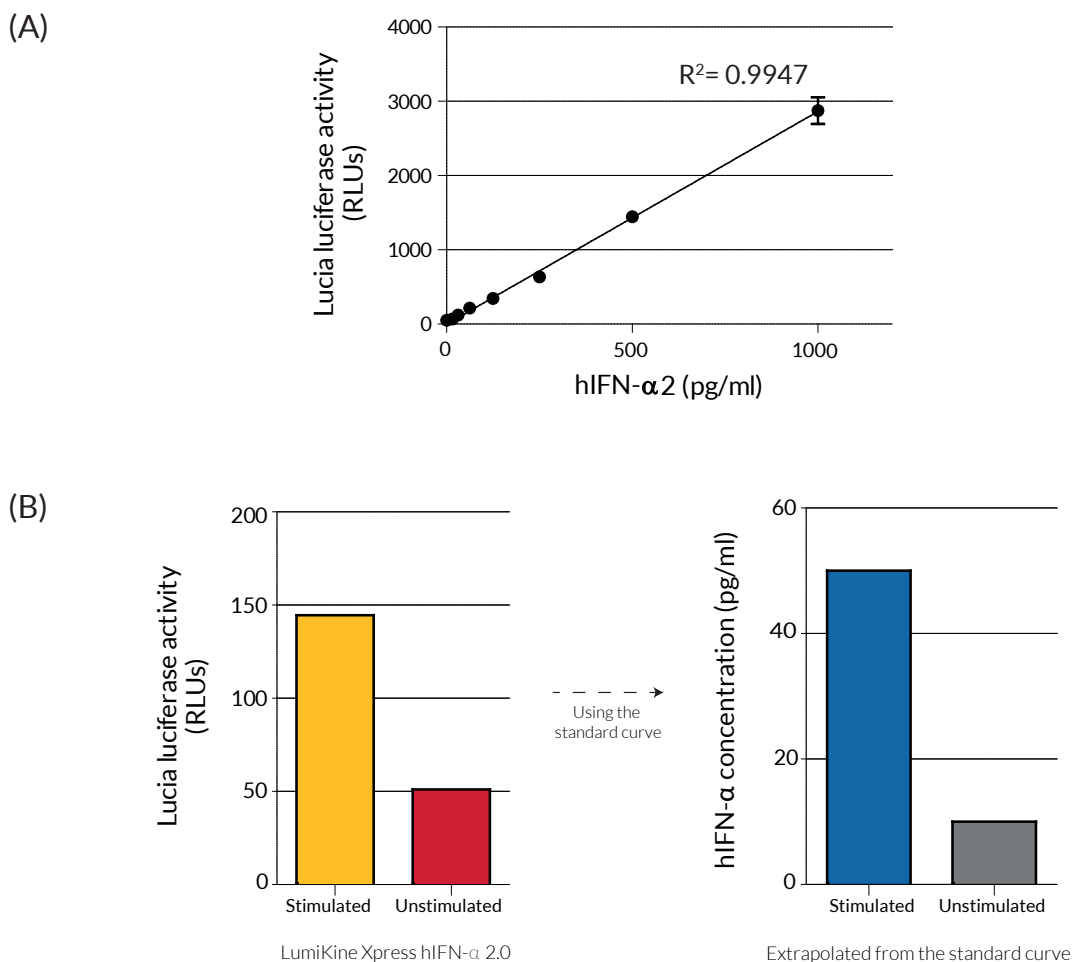
For research use only

Version 19E14-ED

LumiKine™ Xpress hIFN-α 2.0 is a bioluminescent ELISA kit designed to rapidly quantify the levels of human interferon-α (hIFN-α) in cell culture supernatant, serum, and plasma samples. Expression of IFN-α is induced by a number of innate immune pathways including the cGAS-STING signaling pathway upon detection of cytosolic DNA. Unknown hIFN-α levels have been successfully quantified using LumiKine™ Xpress hIFN-α 2.0.

## Determining unknown hIFN-α concentrations

A 7-point standard curve was generated using the standard hIFN-α2 provided in the LumiKine™ Xpress hIFN-α 2.0 kit (Figure 1a). From this, 'unknown' hIFN-α concentrations were determined upon stimulation of human peripheral blood mononuclear cells (PBMCs) with 2'3' cGAMP, a natural STING agonist. hIFN-α was successfully detected (yellow) and quantified (blue) for stimulated cells upon activation of the STING signaling pathway (Figure 1b).



**Figure 1:** (A) A 7-point standard curve (beginning at 1000 pg/ml) was generated using a two-fold serial dilution of standard hIFN-α2. (B) Human peripheral blood mononuclear cells (PBMCs) were either stimulated with 2'3' cGAMP, a STING agonist, or left unstimulated. After 16 hours, the supernatant was isolated and the concentration of hIFN-α in both samples was quantified using LumiKine™ Xpress hIFN-α 2.0.

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)