

# Validation data for LumiKine™ Xpress mIFN-β 2.0

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

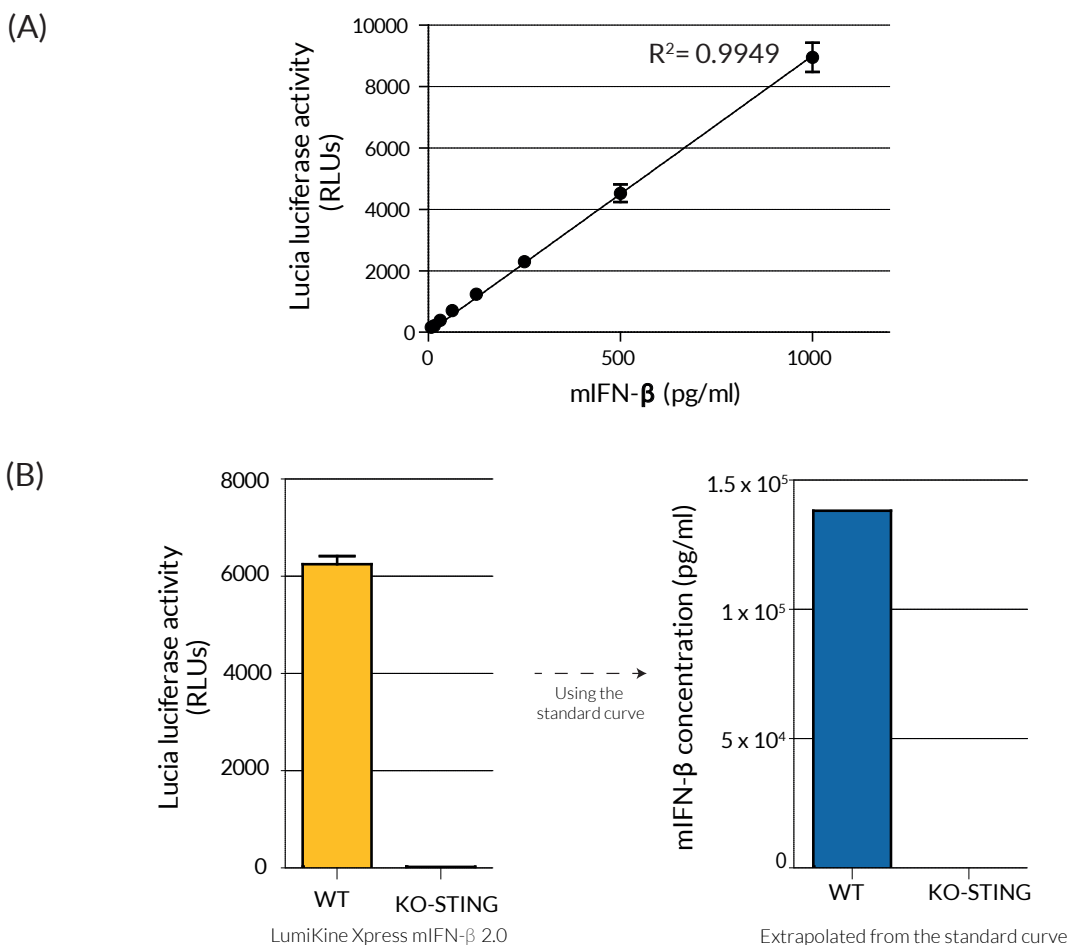
For research use only

Version 19E14-ED

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

## Determining unknown mIFN-β concentrations

A 7-point standard curve was generated using the standard mIFN-β provided in the LumiKine™ Xpress mIFN-β 2.0 kit (Figure 1a). From this, 'unknown' mIFN-β concentrations in sera were determined for wild type (WT) and STING-knock out (KO-STING) mice injected with CL592 (cat code #tlrl-nacai), a synthetic STING agonist. mIFN-β was successfully detected (yellow) and quantified (blue) in the supernatant of WT samples upon activation of the STING signaling pathway, whereas mIFN-β levels were 'not detectable' in the KO-STING samples (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 mIFN-β. (B) Wild type (WT) and STING knock out (KO-STING) mice were injected intraperitoneally with CL592 (2 x 100 mg/kg), a STING agonist. After 4 hours, serum was isolated and diluted 1/200 using DMEM and 10% heat inactivated (HI)-FBS. The concentration of mIFN-β in both samples was quantified using LumiKine™ Xpress mIFN-β 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)