

LumiKine™ Xpress mIFN-β 2.0

Second generation murine IFN-beta bioluminescent ELISA

Catalog code: luex-mifnbv2

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

For research use only

Version 19E14-ED

PRODUCT INFORMATION

Contents (for five plates)

- 150 µg lyophilized mIFN-β capture antibody
- 5 µg lyophilized Lucia-conjugated detection antibody
- 100 ng lyophilized mIFN-β standard
- 2 pouches of QUANTI-Luc™ Plus
- 5 white flat-bottom MaxiSorp® 96-well plates and plate sealers

Storage and stability

- Products are shipped at room temperature.
- Store antibodies, cytokine and QUANTI-Luc™ Plus pouches at -20°C. Lyophilized products are stable for 12 months when properly stored.
- Resuspended antibodies are stable for 1 month when stored at 4°C or 12 months when aliquoted and stored at -20°C. Avoid repeated freeze-thaw cycles.
- Resuspended cytokine is stable for 1 month at 4°C or 12 months when aliquoted and stored at -80°C. Avoid repeated freeze-thaw cycles.
- Reconstituted QUANTI-Luc™ Plus is stable for 1 week at 4°C and for at least 1 month at -20°C. Prepare aliquots to avoid repeated freeze-thaw cycles. Protect from light.

Quality control

- The sensitivity of LumiKine™ Xpress mIFN-β 2.0 has been validated.

PRODUCT DESCRIPTION

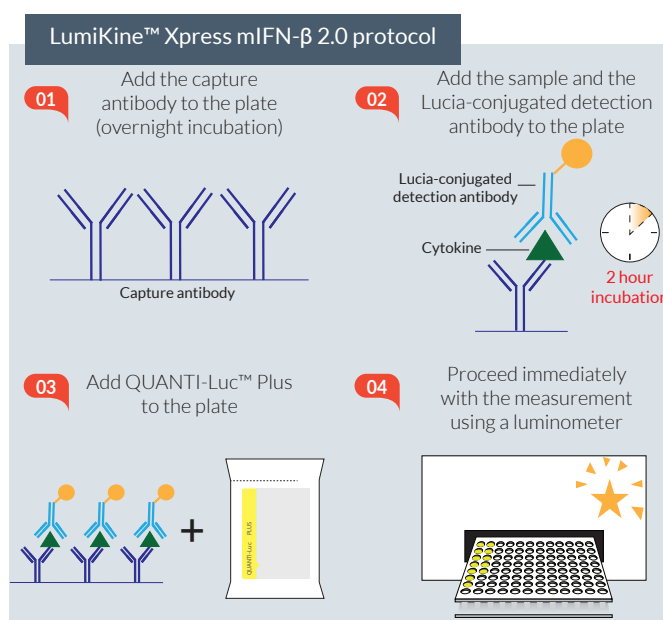
LumiKine™ Xpress mIFN-β 2.0 is a bioluminescent ELISA kit designed to rapidly quantify the levels of murine interferon-β (mIFN-β) in cell culture supernatant, serum, and plasma samples. IFN-β is a type I IFN secreted in response to viral and bacterial infections. Expression of IFN-β is induced by the TBK1-IRF3 signaling axis involving a number of different pattern recognition receptors (PRRs) such as cyclic GMP-AMP synthase (cGAS) and Toll-like receptors (TLRs)¹. The binding of IFN-β to its receptor initiates a signaling cascade, ultimately resulting in the activation of IFN stimulated genes (ISGs)².

LumiKine™ Xpress mIFN-β 2.0 has been enhanced for both faster results and increased sensitivity. This kit uses an optimized pair of capture and detection antibodies. During a single 2-hour incubation, the mIFN-β antigen will bind to **1**) an immobilized mIFN-β capture antibody, as well as **2**) a detection antibody transcriptionally fused to Lucia luciferase; eliminating a step and saving time. Levels of mIFN-β in the sample are determined using QUANTI-Luc™ Plus, a Lucia luciferase detection reagent. The bioluminescent signal is emitted instantaneously and is assessed by luminometry. The intensity of this signal is directly proportional to the concentration of mIFN-β present in the samples.

1. Trinchieri G., 2010. Type I interferon: friend or foe? JEM 207(10):2053 -2063.
2. Schreiber G. 2017. The molecular basis for differential type I interferon signaling. J. Biol. Chem. 292:7285-94.

KEY FEATURES

- **Incubation Time:** 2 hours (after plate preparation)
- **Limit of detection:** 5 pg/ml
- **Specificity:** No cross-reactivity with human (h)IFN-α2, hIFN-β, or mIFN-α2
- **Target:** Natural and recombinant mIFN-β
- **Standard cytokine:** HEK293-expressed mIFN-β
- **Measurement:** Bioluminescent ELISA - relative light units (RLUs)
- **Sample size and type:** 50 µl of cell culture supernatant, serum, or plasma



METHODS

Solutions Required

Prepare the following solutions:

Note: All solutions should be filtered through a 0.2 µm filter before use.

- **Coating buffer:** 0.2M carbonate/bicarbonate buffer (pH 9.4),
Note: Alternatively, you can use sterile PBS (phosphate buffered saline).
- **Blocking buffer:** PBS containing 3% BSA and 0.05% Tween 20
- **Wash buffer:** PBS containing 0.05% Tween 20
- **Reagent diluent:** DMEM, 10% heat inactivated (HI)-FBS
Note: The reagent diluent selected for use can alter the performance of the immunoassay. Optimization of the reagent diluent for samples with complex matrices, such as serum and plasma, may improve the performance of the assay.

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

Reagent Preparation

Note: Bring all reagents to room temperature before use. Allow all components to sit for a minimum of 15 minutes with gentle agitation after initial reconstitution. Working dilutions should be prepared and used immediately.

a) mIFN- β capture antibody stock solution (300 $\mu\text{g/ml}$)

- Add 500 μl of sterile PBS to the vial and mix by pipetting until completely dissolved.
- Use immediately or prepare aliquots and store at -20°C . Avoid repeated freeze-thaw cycles.
- Dilute with coating buffer (1/60) to a working concentration of 5 $\mu\text{g/ml}$.

b) Lucia-conjugated detection antibody stock solution (10 $\mu\text{g/ml}$)

- Add 500 μl of sterile PBS to the vial and mix by pipetting until completely dissolved.
- Use immediately or prepare aliquots and store at -20°C . Avoid repeated freeze-thaw cycles.
- Dilute with reagent diluent to a working concentration of 30 ng/ml.

c) mIFN- β standard stock solution (100 ng/ml)

- Add 1 ml of PBS containing 1% BSA and 0.05% Tween 20 (0.2 μm filtered) to the vial, and mix by pipetting until completely dissolved.
- Use immediately (see 'General ELISA protocol' section (b)) or prepare aliquots and store at -80°C . Avoid repeated freeze-thaw cycles.

d) QUANTI-Luc™ Plus solution

- Pour the pouch contents into a 50 ml screw cap tube.
- Add 25 ml of sterile water.
- Gently swirl product until powder is completely dissolved.
- Use QUANTI-Luc™ Plus solution immediately or aliquot and store at -20°C until required. Avoid repeated freeze-thaw cycles.

Note: This product is photosensitive and must be protected from light.

General ELISA Protocol

a) Plate Preparation

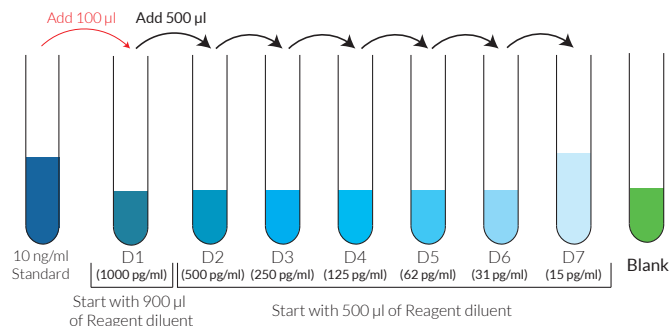
1. Add 50 μl of mIFN- β capture antibody (diluted to 5 $\mu\text{g/ml}$ in coating buffer) to each well of a white flat-bottom MaxiSorp® 96-well plate.
2. Cover the plate with an adhesive seal and incubate overnight at room temperature.
3. Remove excess capture antibody by flicking the plate over a sink and patting it against clean paper towels to remove any remaining drops.
4. Add 200 μl of blocking buffer to each well and incubate for 2 hours at 37°C .
5. Remove blocking buffer by flicking the plate over a sink and patting it against clean paper towels. The plate is now ready for sample addition. Alternatively, the plate can be covered with an adhesive seal and stored at -20°C for 3 months.

b) General Standard Curve Setup

Below is an example protocol for setting up a 7-point standard curve using a 2-fold serial dilution with the highest standard at 1000 pg/ml.

1. To prepare a 10 ng/ml mIFN- β standard solution, add 30 μl of the standard stock solution (100 ng/ml) to 270 μl of reagent diluent
2. Add 100 μl of the 10 ng/ml standard solution to 900 μl of reagent diluent. This is the first dilution (D1; 1000 pg/ml)
3. Add 500 μl of D1 to 500 μl of reagent diluent. This is the second dilution (D2)
4. Continue to perform a 2-fold serial dilution until D7.
5. Ensure you have a negative control (reagent diluent only)

How to set up the 7-point standard curve:



c) Assay Procedure

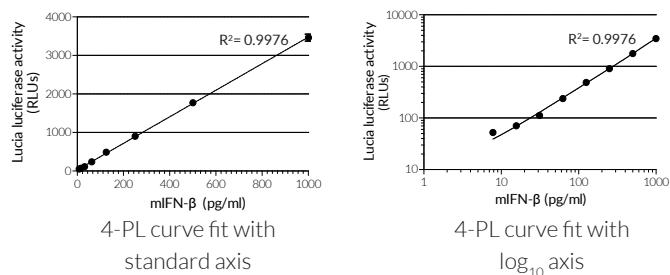
Note: We recommend to test all standards and samples in triplicate.

1. Add 50 μl of each sample (diluted in reagent diluent or an appropriate diluent) per well of the pre-coated plate.
2. Add 50 μl of each prepared standard (including a Blank) to additional wells.
3. Straight away add 50 μl of Lucia-conjugated detection antibody (diluted to 30 ng/ml in reagent diluent) to all wells.
4. Cover with an adhesive seal and incubate for 2 hours at 37°C .
5. Prepare the QUANTI-Luc™ Plus solution (see 'Reagent preparation section (d)'). If frozen bring to room temperature.
6. Remove the liquid by flicking the plate over a sink. Fill each well with 200 μl of wash buffer. Repeat the washing process twice for a total of three washes. After the last wash, remove any remaining Wash buffer by patting the plate against clean paper towels.
7. Set luminometer reading time to the minimum value (0.1-0.5 second).
8. Add 50 μl of reconstituted QUANTI-Luc™ Plus to each well and proceed IMMEDIATELY with the measurement.

STANDARD CURVE DETERMINATION

Create a standard curve using computer software capable of generating a four parameter logistic (4-PL) curve fit. As an alternative, construct a standard curve by plotting the relative light units (RLUs) (triplicate data averaged) on the y-axis, and the different standard concentrations on the x-axis. Draw a best fit curve through as many points as possible on the graph. The data may be linearized by plotting the log of the standard concentrations versus the log of the relative light units (RLUs) and the best fit line can be determined by a regression analysis. This procedure will produce an adequate but less precise fit of the data.

Typical Data



These standard curves are only for demonstration purposes. A standard curve should be generated for each set of samples assayed.

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