

Validation data for B16-Blue™ IFN- α/β cells

<https://www.invivogen.com/b16-blue-ifnab>

For research use only

Version 24G19-NJ

B16-Blue™ IFN- α/β cells allow the detection of bioactive murine type I interferons (i.e. mIFN- α and mIFN- β) by monitoring the activation of the JAK/STAT/ISGF3 pathway. These cells were generated by stable transfection of the murine melanoma B16 cell line with the secreted embryonic alkaline phosphatase (SEAP) reporter under the control of the IFN-inducible ISG54 promoter. B16-Blue™ IFN- α/β cells respond specifically to mIFN- α/β in a dose-dependent manner (Figure 1) and do not respond to human IFN- α/β (Figure 2). Of note, they do not respond to mIFN- γ (Figure 2). Stimulation of these cells with type I IFN inducers, such as poly(dA:dT) delivered intracellularly, triggers the production of SEAP by the activation of the IRF-inducible promoter (Figure 2).

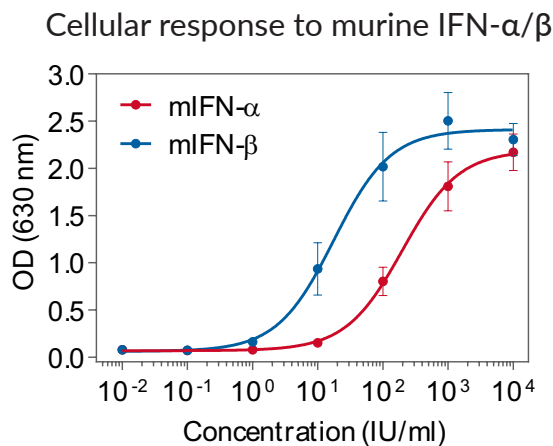


Figure 1. Dose-response of B16-Blue™ IFN- α/β cells to recombinant murine IFN- α/β . Cells were stimulated with increasing concentrations of recombinant murine IFN- α (also known as mIFN- α 3) and mIFN- β . After overnight incubation, the ISGF3 response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent, and reading the optical density (OD) at 630 nm. The OD at 630 nm is shown as mean \pm SEM.

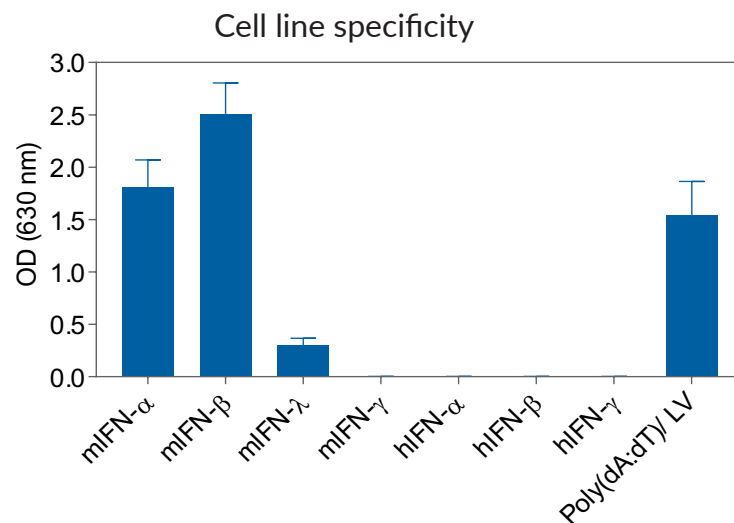


Figure 2. Response of B16-Blue™ IFN- α/β cells to a panel of cytokines. Cells were stimulated with various human and murine recombinant cytokines: 1000 IU/ml mIFN- α , mIFN- β , human IFN- α 2a (hIFN- α 2a), hIFN- β , 100 ng/ml of mIFN- λ , mIFN- γ , hIFN- γ , or 1 μ g/ml of the type I IFN inducer poly(dA:dT) complexed extemporaneously with the transfection reagent LyoVec™ (Poly(dA:dT)/LV). After overnight incubation, SEAP activity was assessed using QUANTI-Blue™ Solution. The OD at 630 nm is shown as mean \pm SEM.

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 Asia: +852 3622-3480
E-mail: info@invivogen.com