

# Validation data for CYT387

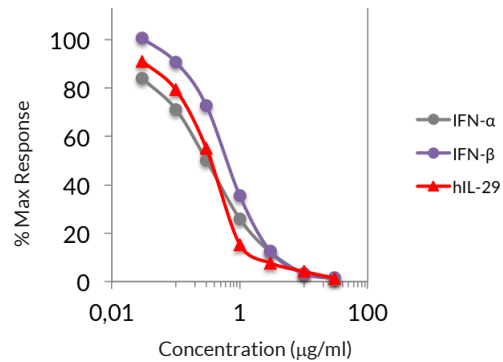
<https://www.invivogen.com/cyt387>

For research use only

Version 18125-MM

CYT387, also known as Momelotinib, is a potent ATP-competitive inhibitor of Janus kinases JAK1 and JAK2, thereby interrupting signaling via the JAK-STAT (signal transducers and activators of transcription) pathway. As a result, CYT387 prevents JAK-STAT activation which plays a pivotal role in cytokine receptor-mediated signal transduction, including interferon (IFN) signaling. Indeed, extracellular signals from IFNs are transduced by JAK-STAT signaling pathway, ultimately leading to the transcription of IFN stimulated genes (ISGs). The ability of CYT387 to inhibit IFN signaling pathway was tested in HEK-Blue™ IFN- $\alpha/\beta$  cells, which respond to type I and type III IFNs. Results obtained with this cell line show that CYT387 can effectively block the IFN- $\lambda$  (IL-29) and the IFN- $\alpha/\beta$  signaling pathway (figure 1).

## Evaluation of inhibitory activity of CYT387



**Figure 1. Effect of CYT387 on HEK-Blue™ IFN- $\alpha/\beta$  cell response to type I and type III IFNs:** HEK-Blue™ IFN- $\alpha/\beta$  cells were incubated with 3 U/ml hIFN- $\alpha$ 2b (grey), 1 U/ml hIFN- $\beta$ 1 (purple) or 10 ng/ml hIL-29 (hIFN- $\lambda$ 1) (red) and increasing concentrations of CYT387. After 24h incubation, IFN-induced ISG activation was assessed by measuring SEAP levels in the supernatant using QUANTI-Blue™. Percentages of maximal response (no inhibitor) for each cytokine are shown.

### 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 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)