

Validation data for recombinant B18R

<https://www.invivogen.com/b18r>

For research use only

Version 23J26-AK

B18R, also called the soluble IFN- α/β receptor B18, is a viral protein naturally encoded by the western reserve (WR) strain of the vaccinia virus. It acts as a decoy receptor of type I interferons (IFNs). B18R has also been reported to increase cell viability and mRNA transfection efficacy, especially during pluripotent (iPSCs) cellular reprogramming. InvivoGen's recombinant B18R successfully blocks type I IFN-mediated signaling, as verified in HEK-Blue™ IFN- α/β cells stimulated for 0 min, 15 min, or 1 hour with recombinant human IFN- α (Figure 1). Also, it increases RNA-transfection efficacy, as verified by transfecting THP1-Dual™ cells with GFP mRNA \pm B18R (Figure 2).

Dose-dependent neutralization of hIFN- α signaling

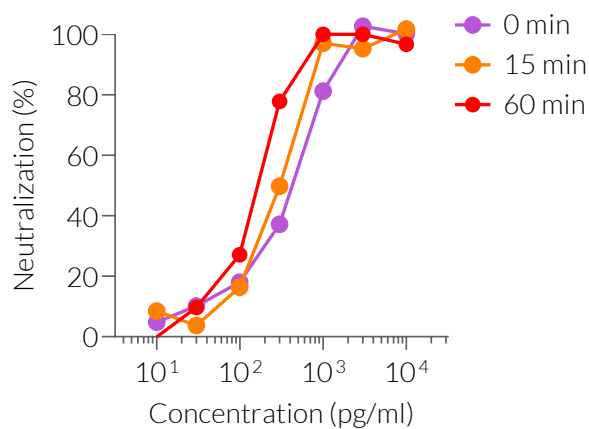


Figure 1. Recombinant B18R is a potent inhibitor of human IFN- α . Increasing concentrations of recombinant B18R were incubated for 0 min, 15 min, or 1 hour with the recombinant human IFN- α 2b (0.1 ng/ml) prior to the addition of HEK-Blue™ IFN- α/β cells. After overnight incubation, SEAP activity in the cell culture supernatant was assessed using QUANTI-Blue™ Solution. Data are shown in percentage (%) of neutralization.

Enhanced mRNA transfection in THP-1 cells

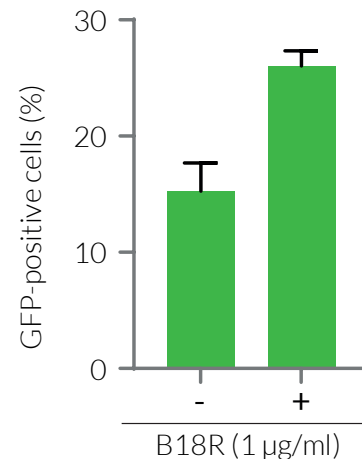


Figure 2. Enhanced mRNA transfection in THP-1 cells. The transfection efficacy of B18R (1 µg/ml) transfecting THP1-Dual™ cells with 100 ng of GFP mRNA using lipofection technology was assessed using flow cytometry. The percentage (%) of GFP-expressing and viable cells is shown.

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 Asia: +852 3622-3480

E-mail: info@invivogen.com