

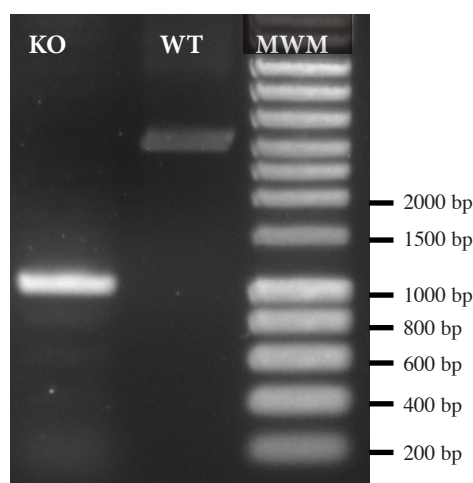
# Validation data for RAW-Lucia™ ISG-KO-IRF3 cells

For research use only

Version # 16C15-MM

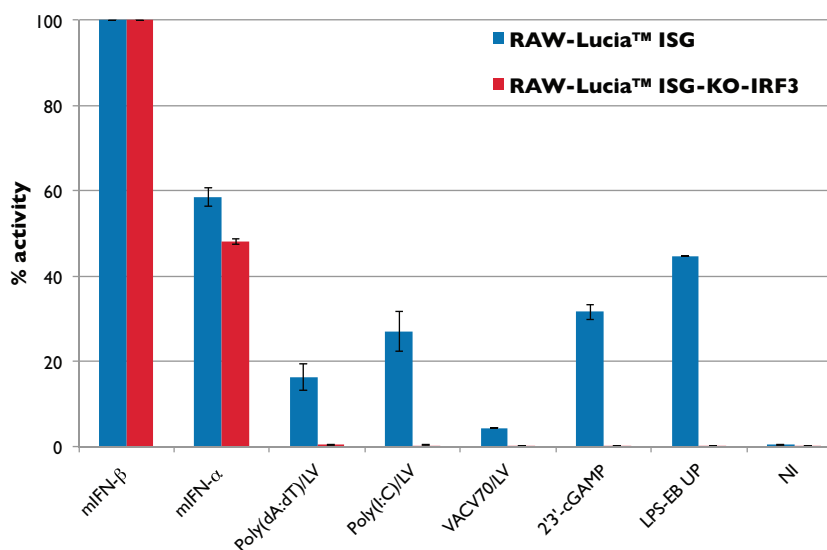
RAW-Lucia™ ISG-KO-IRF3 cells were generated from the RAW-Lucia™ ISG cell line through the stable knockout of the IRF3 gene. These cells derive from the murine RAW 264.7 macrophage cell line. The knockout of the IRF3 gene in RAW-Lucia™ ISG-KO-IRF3 cells has been confirmed by PCR (see figure 1) and sequencing. Biological activity has been assessed using the Lucia luciferase reporter assay to monitor IRF induction (see figure 2).

## PCR AMPLIFICATION



**Figure 1:** PCR amplification of the targeted region in the RAW-Lucia™ ISG-KO-IRF3 (KO) and RAW-Lucia™ ISG (WT) cells. MWM = molecular weight marker

## IRF INDUCTION (Lucia luciferase reporter)



**Figure 2:** Stimulation of RAW-Lucia™ ISG-KO-IRF3 and RAW-Lucia™ ISG cells (parental cell line) with poly(dA:dT)/LyoVec™ (1 µg/ml), VACV70/LyoVec™ (1 µg/ml), poly(I:C)/LyoVec™ (1 µg/ml), and 2'3'-cGAMP (3 µg/ml). Mouse IFN-α (1x10<sup>4</sup> U/ml) and IFN-β (1x10<sup>4</sup> U/ml) serve as positive controls. Non-induced cells (NI) have been included as a negative control. After a 24h incubation, IRF activation was determined by measuring the relative light units (RLUs) in a luminometer using QUANTI-Luc™, a Lucia luciferase detection reagent. The IRF induction of each ligand is expressed relative to that of mIFN-β at 1x10<sup>4</sup> U/ml (taken as 100%).

### 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