

Validation data for RAW-Lucia™ ISG-KO-MDA5 cells

<https://www.invivogen.com/raw-lucia-isg-ko-md5>

For research use only

Version 19K14-MM

RAW-Lucia™ ISG-KO-MDA5 cells were generated from the RAW-Lucia™ ISG cell line through the stable knockout of the MDA-5 (melanoma-differentiation-associated gene 5, also known as Ifih1 or Helicard) gene. These cells derive from the murine RAW 264.7 macrophage cell line, which has been reported to express many pattern recognition receptors (PRRs), including the dsRNA sensors MDA-5 and RIG-I, along with their adaptor protein IPS-1 (also known as MAVS). The knockout of the MDA-5 gene in RAW-Lucia™ ISG-KO-MDA5 cells has been confirmed by PCR (see figure 1), sequencing and Western blot (figure 2).

PCR amplification

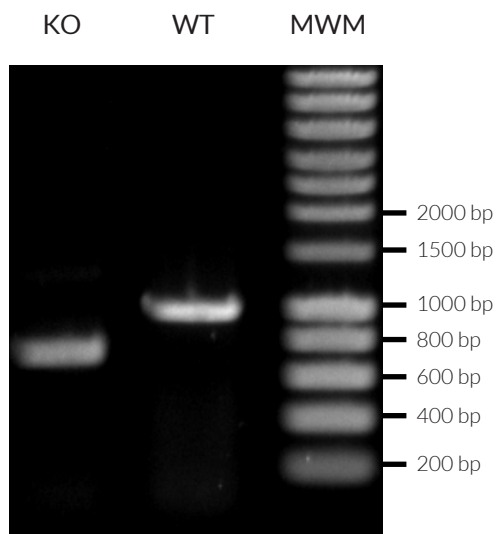


Figure 1: PCR amplification of the targeted region in the RAW-Lucia™ ISG-KO-MDA5 (KO) and RAW-Lucia™ ISG (WT) cells. MWM = molecular weight marker

Western blot

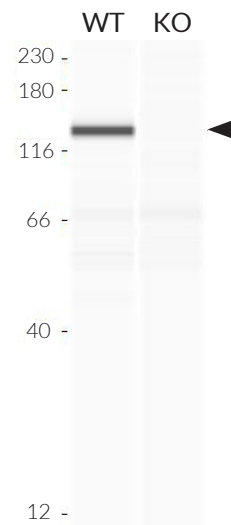


Figure 2: Validation of MDA-5 knockout by Western blot (Wes™). Analysis of lysates from the RAW-Lucia™ (WT) and RAW-Lucia™ ISG-KO-MDA5 (KO) cells using Anti-MDA-5, followed by an HRP-conjugated anti-rabbit secondary antibody. The arrow indicates the expected band for the murine MDA-5 protein (117 KDa).

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