

# Validation data for HEK-Lucia™ RIG-I Cells

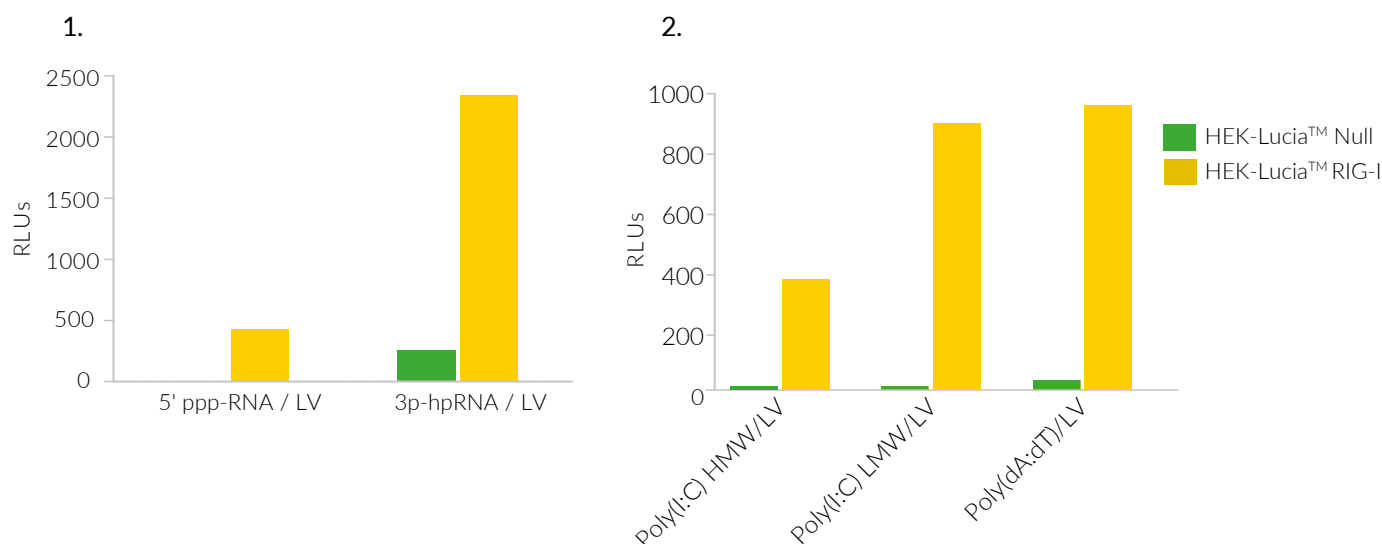
<http://www.invivogen.com/hek-lucia-rigi>

For research use only

Version # 17L08-NJ

HEK-Lucia™ RIG-I cells were generated from HEK-Lucia™ Null cells, HEK293-derived cells that stably express the secreted Lucia luciferase reporter gene. This reporter gene is under the control of an IFN-inducible ISG54 promoter enhanced by a multimeric IFN-stimulated response element (ISRE). HEK-Lucia™ RIG-I cells stably express high levels of human RIG-I, and together with their parental cell line, they can be used to study the role of RIG-I by monitoring IRF-induced Lucia luciferase activity with QUANTI-Luc™. HEK-Lucia™ RIG-I cells respond strongly to cytosolic RIG-I ligands such as 3p-hpRNA and 5'ppp-dsRNA.

## IRF INDUCTION (Lucia luciferase reporter)



HEK-Lucia™ Null and HEK-Lucia™ RIG-I cells were stimulated with 1 µg/ml 5'ppp-dsRNA or 3p-hpRNA (panel 1), or with 100 ng/ml of Poly(I:C) HMW/Lyovec, 100 ng/ml of Poly(I:C) LMW/Lyovec, or 100 ng/ml of Poly(dA:dT)/Lyovec (panel 2). After overnight incubation, the ISG response was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™ and expressed as relative light units (RLUs).

### 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](mailto:info@invivogen.com)