

# Validation data for G3-YSD Control

<https://www.invivogen.com/g3-ysdc>

For research use only

Version 24F17-MM

G3-YSD Control is a negative control for G3-YSD, a potent cGAS (cyclic GMP-AMP synthase, cGAMP synthase) agonist. This negative control differs from the G3-YSD agonist in the hairpin-flanking nucleoside trimers. While G3-YSD is flanked with guanosine trimers (G3) conferring its agonist activity, G3-YSD Control is flanked with cytidine trimers (C3) abrogating cGAS activation. The activity of this ligand has been tested using THP1-Dual™ cells which express multiple cytosolic DNA sensors (CDSs) and two inducible reporter genes (interferon regulatory factor (IRF)-inducible secreted Lucia luciferase and NF- $\kappa$ B-inducible SEAP). Stimulation of these cells with G3-YSD complexed to LyoVec™ leads to a significant IRF response. As expected, intracellular delivery of G3-YSD Control does not induce an IRF response in these cells (Figure 1).

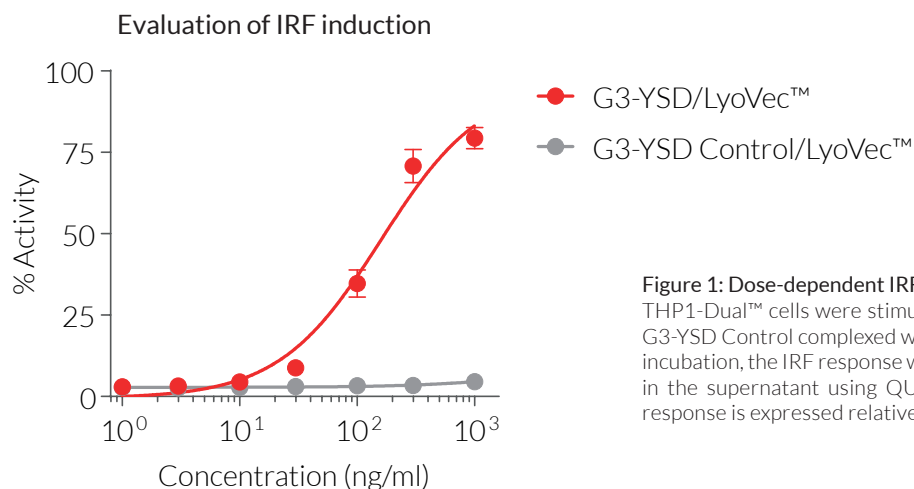


Figure 1: Dose-dependent IRF response in THP1-Dual™ cells.

THP1-Dual™ cells were stimulated with increasing concentrations of G3-YSD or G3-YSD Control complexed with LyoVec™, or  $1 \times 10^4$  U/ml hIFN- $\beta$ . After overnight incubation, the IRF response was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™ 4 Lucia/Gaussia. For each ligand, the response is expressed relative to that of hIFN- $\beta$  (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 Asia: +852 3-622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)