

Validation data for G3-YSD

<http://www.invivogen.com/g3-ysd>

For research use only

Version 18E18-MM

G3-YSD (G3-ended Y-form Short DNA) is a potent agonist for the cytosolic DNA sensor cGAS. This 26-mer DNA palindromic sequence derived from HIV-1 contains unpaired guanosines trimers (G3) in Y-form dsDNA-ssDNA junctions. These guanosine-overhangs have been identified as minimal recognition motifs for cGAS. The activity of this ligand has been tested using the 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- κ B-inducible SEAP).

Stimulation of these cells with the CDS ligands, G3-YSD and VACV-70, complexed to LyoVec™ leads to a significant IRF response (Figure 1). Of note, this response is stronger when using G3-YSD compared to VACV-70. As expected, intracellular delivery of G3-YSD Control does not induce an IRF response in these cells (Figure 1). Importantly, the IRF response to intracellular delivery of G3-YSD is strictly cGAS-dependent (Figure 2).

Evaluation of IRF induction

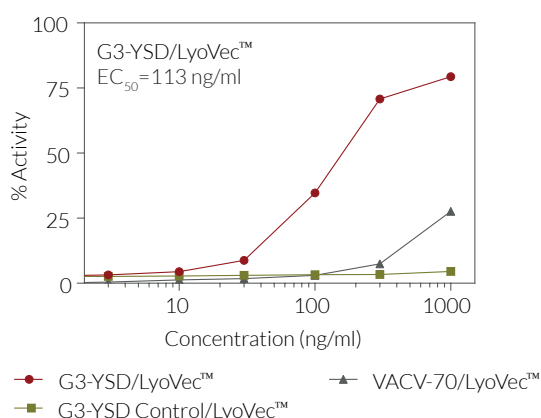


Figure 1: Intracellular delivery of G3-YSD induces a potent IRF response in a dose-dependent manner.

THP1-Dual™ cells were stimulated with increasing concentrations of G3-YSD, G3-YSD Control or VACV-70 complexed with LyoVec™, or 1×10^4 U/ml hIFN- β . After overnight incubation, the IRF response was assessed by determining Lucia luciferase activity in the supernatant using QUANTI-Luc™. For each ligand, the response is expressed relative to that of hIFN- β (taken as 100%).

Evaluation of signaling pathway

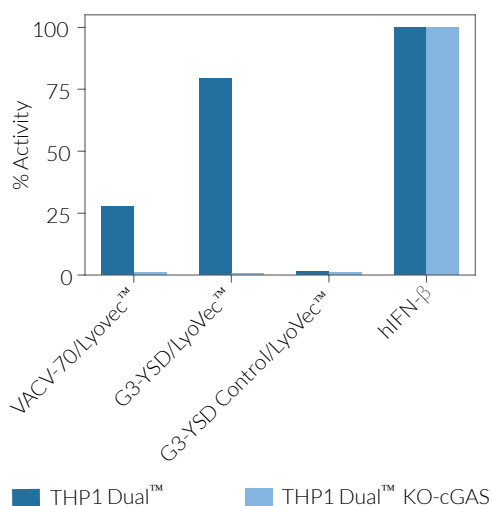


Figure 2: IRF response upon intracellular delivery of G3-YSD is strictly dependent on cGAS.

THP1-Dual™ and THP1-Dual™ KO-cGAS cells were stimulated with $1 \mu\text{g/ml}$ G3-YSD, G3-YSD Control or VACV-70 complexed with LyoVec™, or 1×10^4 U/ml hIFN- β . After overnight incubation, the IRF response was assessed using QUANTI-Luc™ and expressed relative to that of hIFN- β (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