

Validation data for STG-968

<https://www.invivogen.com/sting-conjugatable-ligands>

For research use only

Version 23J12-NJ

STG-968 is a “click-chemistry compatible” conjugatable STING ligand featuring an azido group added to the N-terminus of an analog of CL656, a well-known STING agonist. STG-968 efficiently triggers IRF- and NF- κ B-mediated cellular responses (Figure 1). STG-968 can be used to generate immunostimulatory antibody-drug conjugates (ADCs) as conjugation to a Anti-HER2-hlgG1 and subsequent activation of STING has been validated using cellular assays. The Anti-HER2/STG-968 ADC is more potent at inducing a STING-mediated response than STG-968 only in cells expressing HER2 (Figure 2A). Of note, at high concentrations, STG-968, Anti-HER2/STG-968 ADC and Anti- β -Gal/STG-968 ADC, induce a STING-mediated cellular response, independently of HER2 expression (Figure 2A and B). This observation could be explained by cellular uptake through endocytosis/pinocytosis.

Biological activity of STG-968

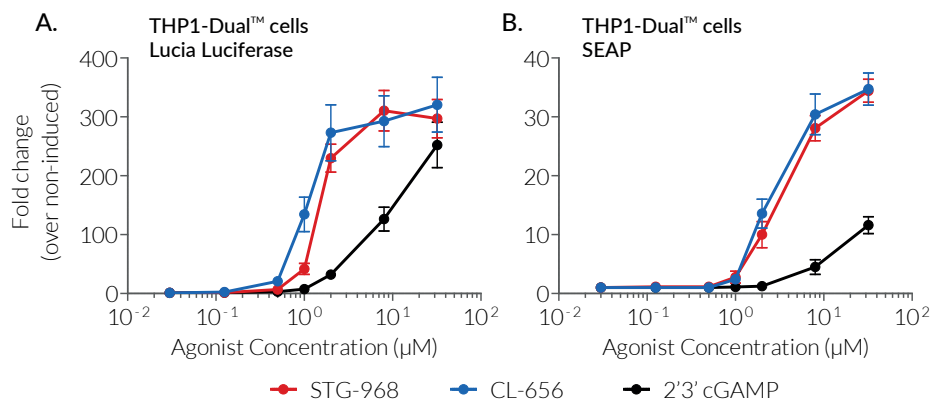


Figure 1: IRF and NF- κ B responses induced by STING conjugatable ligand STG-968.

THP1-Dual™ cells were stimulated with increasing concentrations of STG-968, CL-656, or 2'3'-cGAMP. After overnight incubation, the IRF and NF- κ B responses were determined by measuring Lucia luciferase and SEAP activity in the supernatant using QUANTI-Luc™ (A), or QUANTI-Blue™ Solution (B), respectively. Data are shown as a fold increase (mean \pm SEM) over non-induced cells.

Biological activity of STG-968 conjugated to Anti-HER2-hlgG1

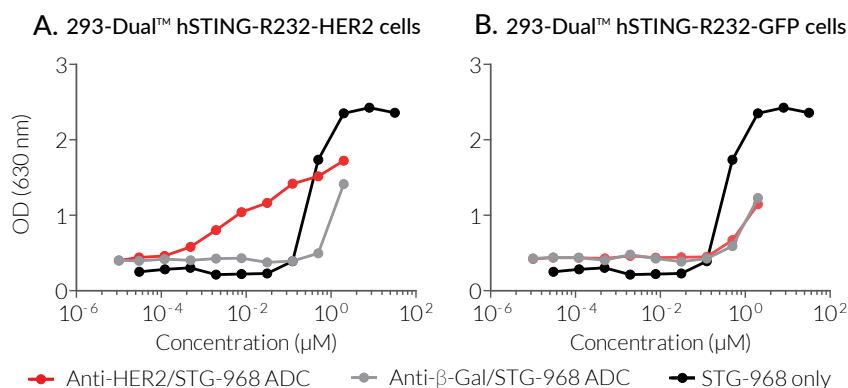


Figure 2: Dose-response of HER2-expressing STING reporter cells to Anti-HER2/STG-968 ISAC.

~ 5x10⁵ 293-Dual™ hSTING-R232-HER2 cells (A) or 293-Dual™ hSTING-R232-GFP control cells (B) were stimulated with increasing concentrations of Anti-HER2/STG-968 ADC (Ratio 1:5), Anti- β Gal/STG-968 ADC (Ratio 1:5), or STG-968 only. After overnight incubation, the STING response was determined using QUANTI-Blue™ Solution, a SEAP detection reagent. The optical density (OD) at 630 nm is shown.

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 3622-3480
E-mail: info@invivogen.com