

Validation data for diABZI

<https://www.invivogen.com/diabzi>

For research use only
Version 21E28-ED

diABZI also known as diABZI (compound 3) trihydrochloride, is a potent, non-cyclic dinucleotide STING agonist. The activation of STING leads to IFN regulatory factor (IRF3)-dependent type I IFN production and NF- κ B-dependent inflammatory cytokine production. The activity diABZI has been validated using InvivoGen's THP1-Dual™ cells, a cell line derived from the human THP-1 monocyte cell line by stable integration of two inducible reporter constructs. As a result, these cells allow the simultaneous study of the IRF pathway by assessing the activity of a secreted Lucia luciferase, and the NF- κ B pathway by monitoring the activity of SEAP. Stimulation of THP1-Dual™ cells with diABZI results in a dose-dependent activation of STING. Furthermore, diABZI is more potent when compared to the canonical STING ligand, 2'3'-cGAMP (Figure 1).

Evaluation of STING activation by diABZI

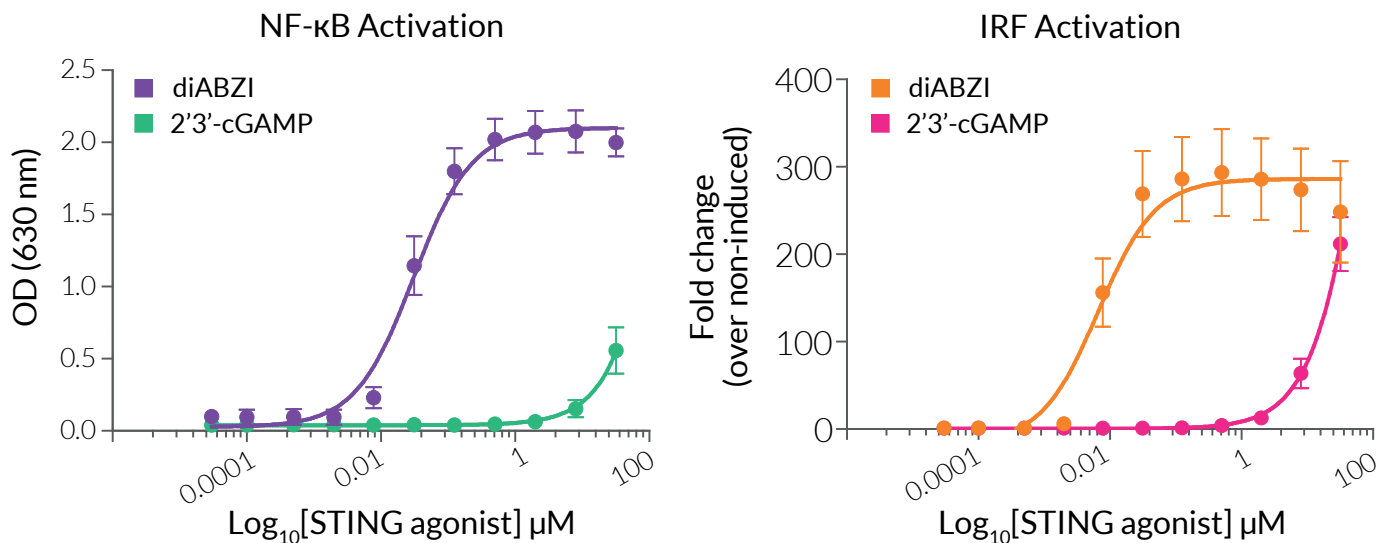


Figure 1: diABZI induces a dose-dependent response in THP1-Dual™ cells. The cells were incubated with increasing concentrations of diABZI or 2'3'-cGAMP. After overnight incubation (A) the NF- κ B activity was assessed by measuring the SEAP activity in the supernatant using QUANTI-Blue™ Solution. Optical density (OD) at 630 nm is shown. (B) The IRF response was assessed by measuring the activity of Lucia luciferase in the supernatant using QUANTI-Luc™. Fold change over non-induced cells 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