

Validation data for ODN 2336

<https://www.invivogen.com/odn2336>

For research use only

Version 24B20-AK

ODN 2336 is a synthetic immunostimulatory oligonucleotide (ODN) that contains unmethylated CpG dinucleotides. ODN 2336 is a Class A CpG ODN with a preference for the human Toll-like receptor 9 (hTLR9). ODN 2336 is able to activate the hTLR9-mediated NF- κ B and IRF pathways as verified using InvivoGen's HEK-Dual™ hTLR9 cells (Figure 1). These cells express the human *TLR9* gene as well as two inducible reporter genes for the NF- κ B-inducible SEAP (secreted embryonic alkaline phosphatase) and IRF-inducible Lucia luciferase.

Dose-dependent activation of human TLR9

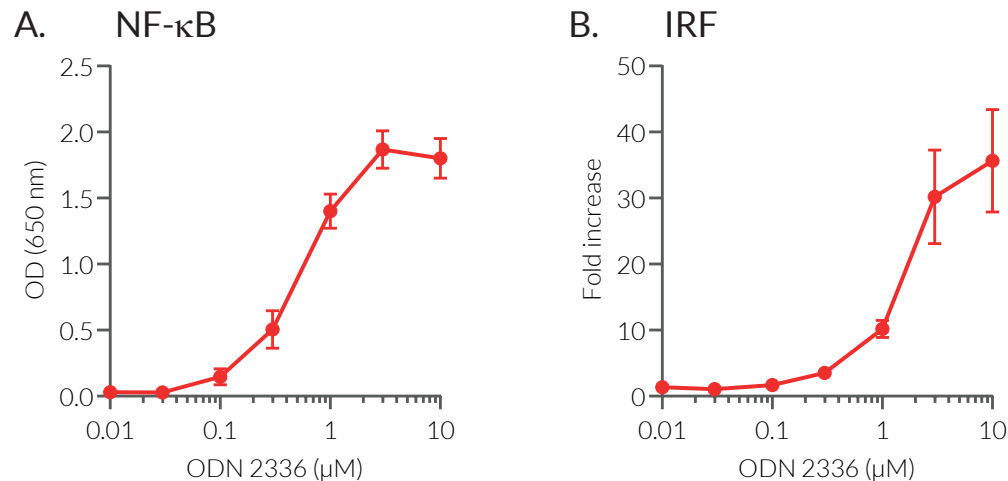


Figure 1. ODN 2336 is a potent activator of human TLR9. HEK-Dual™ hTLR9 cells were incubated with increasing concentrations of ODN 2336. After 24h, the hTLR9-induced (A) NF- κ B and (B) IRF responses were assessed by measuring SEAP and Lucia activity using QUANTI-Blue™ and QUANTI-Luc™, respectively. Data are shown as optical density (OD) at 650 nm or in fold increase over non-induced cells (mean + SEM).

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