

Validation data for ODN 2006 (ODN 7909)

<https://www.invivogen.com/odn2006>

For research use only

Version 23K10-MM

ODN 2006 (ODN 7909) is a Class B CpG oligonucleotide (ODN). It is a ligand of choice for human Toll-like receptor 9 (TLR9). Activation of TLR9 with ODN 2006 triggers NF- κ B- and interferon regulatory factor (IRF)-mediated pro-inflammatory responses. The biological activity of ODN 2006 has been tested using InvivoGen's THP1-Dual™ hTLR9 cells (Figure 1). These cells express human TLR9 gene and two inducible reporter genes, allowing the concomitant study of the IRF and NF- κ B pathways, by monitoring the Lucia luciferase and SEAP (secreted embryonic alkaline phosphatase) activities, respectively.

Dose-dependent activation of human TLR9

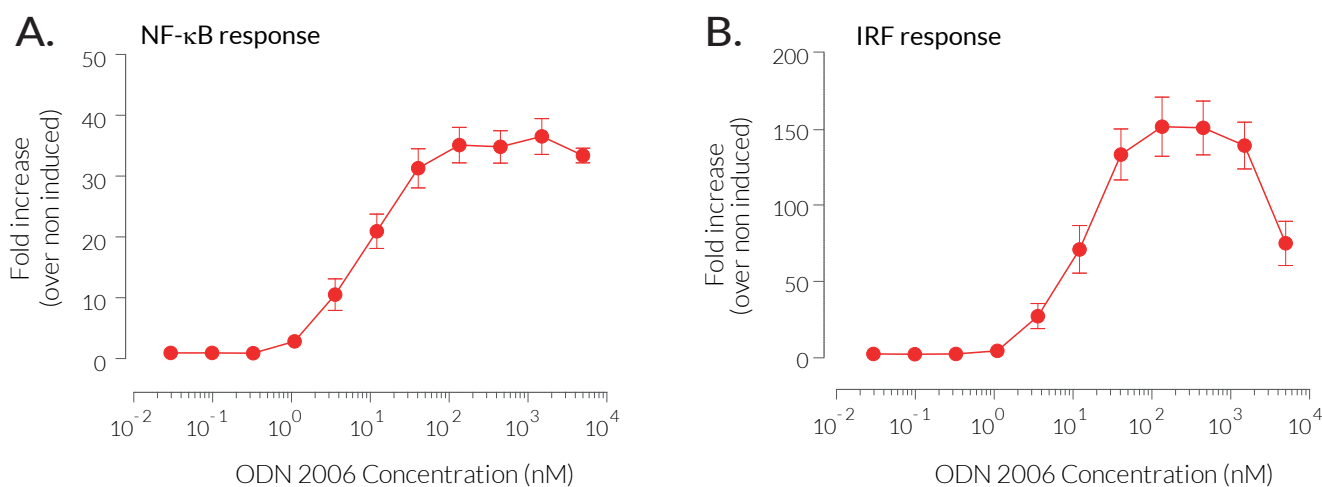


Figure 1. NF- κ B and IRF responses induced by ODN 2006. THP1-Dual™ hTLR9 cells were incubated with increasing concentrations of ODN 2006. After overnight incubation, the NF- κ B response was assessed by measuring the SEAP activity using QUANTI-Blue™ Solution (A) and the IRF response was assessed by measuring the Lucia luciferase activity using QUANTI-Luc™ 4 Lucia/Gaussia (B). Data are shown as a fold increase (mean \pm 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