

Validation data for ODN 2395

<https://www.invivogen.com/odn2395>

For research use only

Version 24C14-AK

ODN 2395 is a synthetic immunostimulatory oligonucleotide (ODN) that contains unmethylated CpG dinucleotides. ODN 2395 is a Class C CpG ODN specific for the human (h) and mouse (m) Toll-like receptor 9 (TLR9). In HEK-Blue™-derived reporter cells, ODN 2395 efficiently activates hTLR9 and mTLR9 (Figure 1). Moreover, ODN 2395 is able to activate the hTLR9-mediated NF- κ B and IRF pathways as verified using InvivoGen's THP1-Dual™ hTLR9 cells (Figure 2). This monocytic cell line expresses 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 NF- κ B response in HEK-Blue™-derived cells

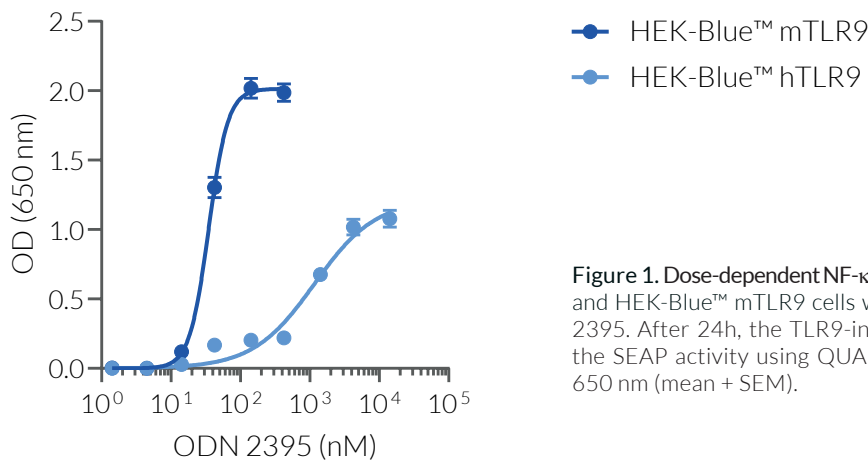


Figure 1. Dose-dependent NF- κ B response in HEK-Blue™-derived cells. HEK-Blue™ hTLR9 and HEK-Blue™ mTLR9 cells were incubated with increasing concentrations of ODN 2395. After 24h, the TLR9-induced NF- κ B responses were assessed by measuring the SEAP activity using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean + SEM).

Dose-dependent NF- κ B and IRF responses in THP1-Dual™ hTLR9 cells

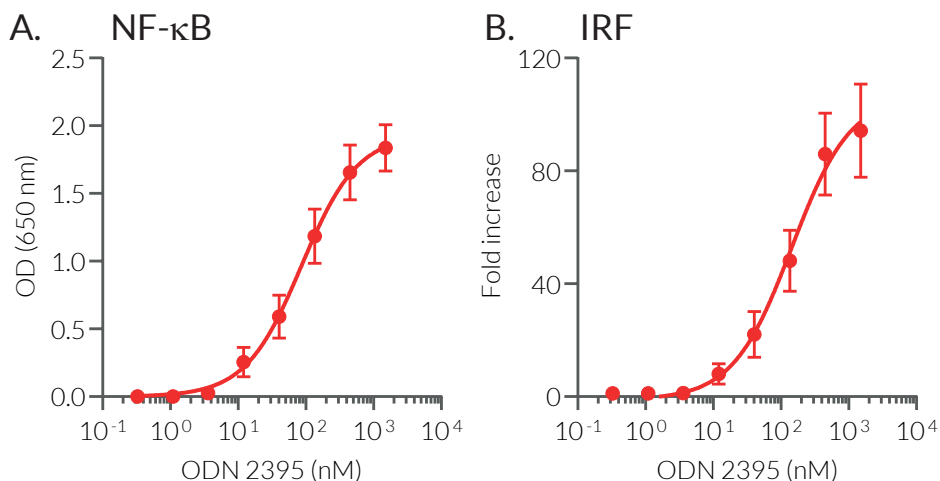


Figure 2. Dose-dependent NF- κ B and IRF responses in THP1-Dual™ hTLR9 cells. Cells were incubated with increasing concentrations of ODN 2395. After 24h, the hTLR9-induced (A) NF- κ B and (B) IRF responses were assessed by measuring SEAP and Lucia activity using QUANTI-Blue™ and QUANTI-Luc™, respectively. 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