

# Validation data for ODN 1826

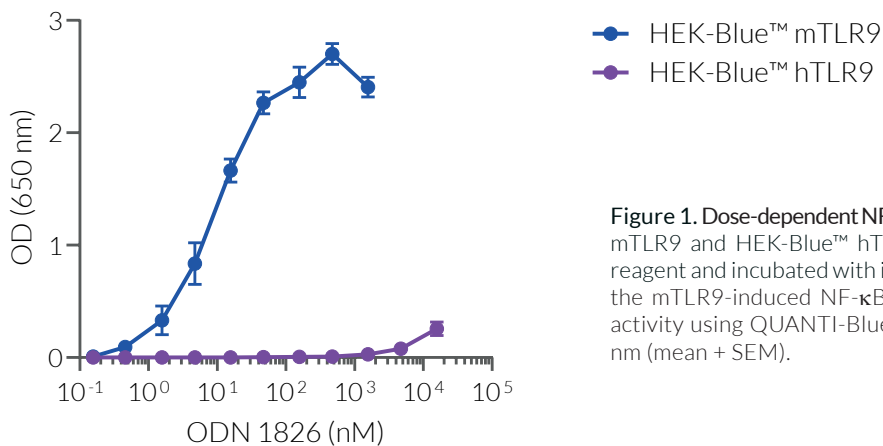
<https://www.invivogen.com/odn1826>

For research use only

Version 24B22-AK

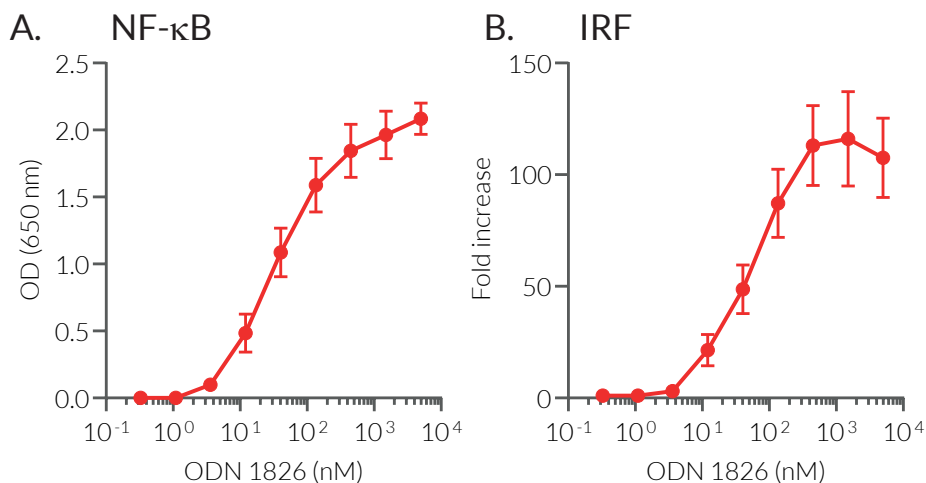
ODN 1826 is a synthetic immunostimulatory oligonucleotide (ODN) that contains unmethylated CpG dinucleotides. ODN 1826 is a Class A CpG ODN with a preference for the mouse Toll-like receptor 9 (mTLR9). In HEK-Blue™-derived reporter cells, ODN 1826 efficiently activates mTLR9, but not human (h)TLR9 (Figure 1). Interestingly, ODN 1826 is able to activate the hTLR9-mediated NF-κB and IRF pathways as verified using InvivoGen's THP1-Dual™ hTLR9 cells (Figure 2). This monocytic cell line overexpresses the human *TLR9* gene as well as two inducible reporter genes for the NF-κB-inducible SEAP (secreted embryonic alkaline phosphatase) and IRF-inducible Lucia luciferase.

## Dose-dependent NF-κB response in HEK-Blue™-derived cells



**Figure 1.** Dose-dependent NF-κB response in HEK-Blue™-derived cells. HEK-Blue™ mTLR9 and HEK-Blue™ hTLR9 cells were cultured in HEK-Blue™ Detection reagent and incubated with increasing concentrations of ODN 1826. After 24h, the mTLR9-induced NF-κB response was assessed by measuring the SEAP activity using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (mean + SEM).

## Dose-dependent NF-κB and IRF responses in THP1-Dual™ hTLR9 cells



**Figure 2.** Dose-dependent NF-κB and IRF responses in THP1-Dual™ hTLR9 cells. Cells were incubated with increasing concentrations of ODN 1826. After 24h, the hTLR9-induced (A) NF-κB 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](mailto:info@invivogen.com)