

# Validation data for ODN 2007

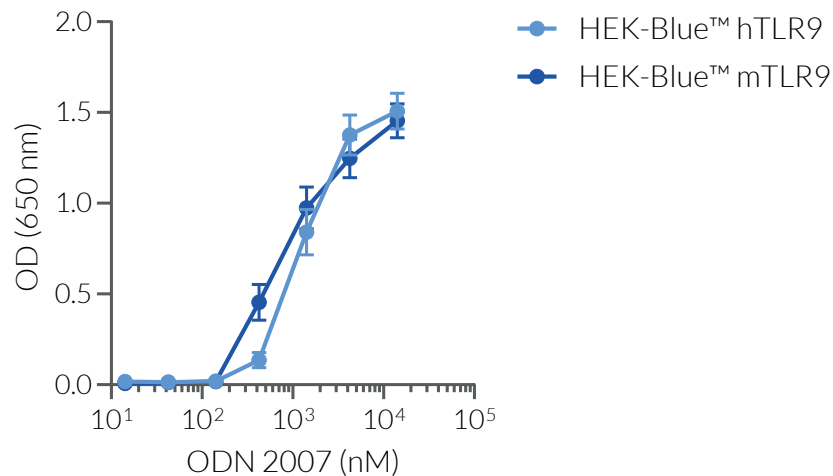
<https://www.invivogen.com/odn2007>

For research use only

Version 24C14-AK

ODN 2007 is a synthetic immunostimulatory oligonucleotide (ODN) that contains unmethylated CpG dinucleotides. ODN 2007 is a Class B CpG ODN with a preference for the Toll-like receptor 9 (TLR9) of cows, pigs, and chickens. It is also able to activate the NF- $\kappa$ B pathway mediated by the human (h) and mouse (m) TLR9, as verified using InvivoGen's HEK-Blue™-derived reporter cells (Figure 1). These cell lines express the human or murine *TLR9* gene as well as an inducible reporter gene for the NF- $\kappa$ B-inducible SEAP (secreted embryonic alkaline phosphatase).

## Dose-dependent NF- $\kappa$ B response in HEK-Blue™-derived cells



**Figure 1.** Dose-dependent NF- $\kappa$ B response in HEK-Blue™-derived cells. HEK-Blue™ hTLR9 and HEK-Blue™ mTLR9 cells were incubated with increasing concentrations of ODN 2007. After 24h, the TLR9-induced NF- $\kappa$ B responses were assessed by measuring the SEAP activity using QUANTI-Blue™. Data are shown as optical density (OD) at 650 nm (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)