

# ODN 1668

## Class B CpG oligonucleotide; a murine TLR9 ligand

Catalog # tlr1-1668, tlr1-1668-1, tlr1-1668-5

For research use only

Version # 16E17-MM

### PRODUCT INFORMATION

#### Content:

- ODN 1668 is provided lyophilized and is available in three quantities:
  - 200 µg (**31.42 nmol**): tlr1-1668 (formerly tlr1-modnb)
  - 1 mg (**157.1 nmol**): tlr1-1668-1 (formerly tlr1-modnb-1)
  - 5 x 1 mg (5 mg; **785.5 nmol**): tlr1-1668-5 (formerly tlr1-modnb-5)

*Note: ODN 1668 is sterile filtered prior to lyophilization.*

- endotoxin-free water; 1.5 ml with #tlr1-1668 and tlr1-1668-1, and 10 ml with #tlr1-1668-5.

#### ODN 1668 sequence

5'-tccatgacgttcctgatgct-3' (20 mer)

*Note: Bases are phosphorothioate-linked (nuclease resistant).*

**Molecular weight:** 6364 g/mol

#### Storage and stability

- ODN 1668 is shipped at room temperature. Upon receipt, store at -20 °C.
- Upon resuspension, prepare aliquots of ODN 1668 and store at -20 °C. Resuspended product is stable for 6 months at -20 °C when properly stored. Avoid repeated freeze-thaw cycles.

#### Quality control

- TLR9 activity has been tested using HEK-Blue™ TLR9 cells.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

### DESCRIPTION

CpG ODNs are synthetic oligonucleotides that contain unmethylated CpG dinucleotides in particular sequence contexts (CpG motifs)<sup>1</sup>. These CpG motifs are present at a 20-fold greater frequency in bacterial DNA compared to mammalian DNA. CpG ODNs are recognized by Toll-like receptor 9 (TLR9) leading to strong immunostimulatory effects<sup>2</sup>. Three classes of stimulatory CpG ODNs have been identified, classes A, B and C, which differ in their immune-stimulatory activities<sup>3,4</sup>.

ODN 1668 is a class B CpG ODN with a preference for mouse TLR9. Class B CpG ODNs contain a full phosphorothioate backbone with one or more CpG dinucleotides. They strongly activate B cells but stimulate weakly IFN-α secretion.

**1. Krieg, A. et al., 1995.** CpG motifs in bacterial DNA trigger direct B-cell activation. *Nature*, 374:546-9. **2. Bauer, S. et al., 2001.** Human TLR9 confers responsiveness to bacterial DNA via species-specific CpG motif recognition. *PNAS*, 98:9237-42. **3. Krug A. et al., 2001.** Identification of CpG oligonucleotide sequences with high induction of IFN-alpha/beta in plasmacytoid dendritic cells. *Eur J Immunol*, 31:2154-63. **4. Marshall J. et al., 2005.** Superior activity of the type C class of ISS in vitro and in vivo across multiple species. *DNA Cell Biol.* 24(2):63-72.

### METHODS

#### Preparation of stock solution (500 µM)

TLR9 activation can be achieved with 1-5 µM ODN 1668.

- Resuspend ODN 1668 with endotoxin-free water (provided).
  - Add 63 µl to 200 µg of ODN 1668
  - Add 315 µl to 1 mg of ODN 1668
- Vortex until completely dissolved. Prepare aliquots and store at -20 °C.
- Prepare serial dilutions using endotoxin-free water.

*Note: The working concentration may vary depending on the levels of TLR9 gene expression and the species from which the gene was obtained.*

#### TLR9 stimulation using ODN 1668

ODN 1668 can be used to stimulate TLR9 in HEK-Blue™ TLR9 cells. HEK-Blue™ TLR9 cells stably overexpress the TLR9 gene and an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene. For more information, visit: [www.invivogen.com](http://www.invivogen.com)

Below is a protocol to study TLR9 stimulation using HEK-Blue™ TLR9 cells in a 96-well plate.

- Dispense 20 µl of stimulatory or control ODN per well of a 96-well plate.
- Prepare cell suspension of HEK-Blue™ TLR9 cells according to the data sheet.
- Add HEK-Blue™ TLR9 cells (4-8 x10<sup>4</sup>) to each ODN-containing well.
- Incubate for 6-24 h at 37 °C, 5% CO<sub>2</sub>.
- Determine TLR9 stimulation by assessing cytokine expression using ELISA, or SEAP expression using QUANTI-Blue™, a SEAP detection medium.

### RELATED PRODUCT

Product	Catalog Code
ODN 1668 Control	tlr1-1668c
pUNO1-mTLR9 (mouse TLR9 gene)	puno1-mtlr9
HEK-Blue™ mTLR9 Cells	hkb-mtlr9
QUANTI-Blue™	rep-qb1

#### 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 Hong Kong: +852 3-622-34-80  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)