

# ODN 1668 Control

Negative control oligonucleotide for murine TLR9 ligand ODN 1668

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

For research use only

Version # 16E24-MM

## PRODUCT INFORMATION

### Content

- ODN 1668 Control is provided lyophilized and is available in 3 quantities:
  - 200 µg (**31.42 nmol**): tlr1-1668c (formerly tlr1-modnbc)
  - 1 mg (**157.10 nmol**): tlr1-1668c-1 (formerly tlr1-modnbc-1)
  - 5 x 1 mg (5 mg; **785.50 nmol**): tlr1-1668c-5 (formerly tlr1-modnbc-5)

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

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

### ODN1668 Control sequence

5'-tccatgagcttctgatgct-3' (20 mer)

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

**Molecular weight:** 6364 g/mol

### Storage and stability

- ODN 1668 Control is shipped at room temperature. Upon receipt, store at -20 °C.

- Upon resuspension, prepare aliquots of ODN 1668 Control 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

- 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 Control contains GpC dinucleotides instead of CpGs and can be used as a negative control for ODN 1668 (class B CpG ODN).

**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)

- Resuspend ODN 1668 Control with endotoxin-free water (provided).
  - Add 63 µl to 200 µg of ODN 1668 Control
  - Add 315 µl to 1 mg of ODN 1668 Control
- 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

ODN 1668 Control can be used as a control ODN to study the stimulatory effect of ODN 1668 on TLR9 in HEK-Blue™ TLR9 cells. HEK-Blue™ TLR9 cells stably overexpress the TLR9 gene and an NF- $\kappa$ B-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.

*Note: Use the ODN 1668 Control at the same concentration as the CpG-containing ODN 1668.*

- 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                      | tlr1-1668    |
| 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)