

ODN 1668 FITC

FITC labeled CpG oligonucleotide - Mouse TLR9 ligand

Catalog # tlr1-1668f (formerly tlr1-fmodnb)

For research use only

Version # 16J28-MM

PRODUCT INFORMATION

Content:

- 50 µg (7.25 nmol) lyophilized ODN1668 labeled with FITC at the 3' terminus.
- 1.5 ml endotoxin-free water

ODN 1668 sequence

5'-tccatgacgttctctgatgct-3' (20 mer)

Note: Bases are phosphorothioate (nuclease resistant).

Molecular weight: 6947 g/mol

Storage :

- ODN 1668 FITC is shipped at room temperature. Store at -20°C. Protect from light.
- Resuspended product should be stored at -20°C and protected from light. Product is stable for 6 months. 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)¹. 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 effects². Three classes of stimulatory CpG ODNs have been identified, classes A, B and C, which differ in their immune-stimulatory activities³⁻⁴. 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 CpG ODN solution (500 µM)

- TLR9 activation can be achieved with 1-5 µM ODN 1668 FITC.
- Resuspend lyophilized ODN 1668 FITC with 14 µl of endotoxin-free water (provided).
 - Vortex until complete solubilization.
 - Store at -20°C.

TLR9 stimulation using ODN 1668 FITC

ODN 1668 FITC 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

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 x10⁴) to each ODN-containing well.
- Incubate for 6-24 h at 37°C, 5% CO₂.
- Determine TLR9 stimulation by assessing cytokine expression using ELISA, or SEAP expression using QUANTI-Blue™, a SEAP detection medium.
- Evaluate CpG ODN cellular uptake and localization by confocal laser-scanning microscopy (excitation 495 nm, emission 520 nm) or flow cytometry.

RELATED PRODUCTS

Product	Catalog Code
ODN 1668	tlr1-1668
pUNO-mTLR9 (mouse TLR9 gene)	puno-mtlr9
HEK-Blue™ mTLR9 Cells	hkb-mtlr9
QUANTI-Blue™	rep-qb1
293/mTLR9 Cells	293-mtlr9
pNiFty2-Luc (Zeo [®])	pnifty2-luc
pNiFty2-SEAP (Zeo [®])	pnifty2-seap

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

