

FITC ODN 2006

FITC labeled CpG oligonucleotide - Human TLR9 ligand

Catalog # tlr1-2006f (formerly tlr1-fhodnb)

For research use only

Version # 16J20-MM

PRODUCT INFORMATION

Content

- 50 µg (6.2 nmol) lyophilized ODN 2006 labeled with FITC at the 3' terminus.
- 1.5 ml endotoxin-free water

ODN 2006 sequence

5'-tcgtcgtttgtcgtttgtcgtt-3' (24 mer)

Note: Bases are phosphorothioate (nuclease resistant).

Molecular weight: 8171 g/mol

Storage and stability

- FITC ODN 2006 is shipped at room temperature and should be stored at -20 °C for up to 1 year. Protect from light.
- Resuspended product should be stored at -20 °C and protected from light. Product is stable for 6 months. Avoid repeated cycles of freeze-thaw.

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 immunostimulatory activities^{3,4}.

ODN 2006 (also known as ODN 7909 or PF-3512676) is a CpG ODN class B with a preference for human TLR9. B-class CpG ODNs contain a full phosphorothioate backbone with one or more CpG dinucleotides. They strongly activate B cells but weakly stimulate 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- α /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 2006 FITC.
- Resuspend lyophilized labeled ODN 2006 with 12 µl of sterile endotoxin-free water provided.
 - Vortex until complete solubilization.
 - Aliquote and store at -20 °C.

TLR9 stimulation using ODN 2006 FITC

ODN 2006 FITC can be used to stimulate TLR9 in HEK-Blue™ TLR9 cells. HEK-Blue™ TLR9 cells stably overexpress the TLR9 gene and an NF- κ B-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 2006 (ODN 7909)	tlr1-2006
pUNO1-hTLR9a (human TLR9 gene)	puno1-htlr9a
HEK-Blue™ hTLR9 cells	hkb-htlr9
QUANTI-Blue™	rep-qb1
293XL/hTLR9	293xl-htlr9
pNiFty2-Luc (Zeo ^R)	pnifty2-luc
pNiFty2-SEAP (Zeo ^R)	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