

# ODN 1826 FITC

## FITC labeled CpG oligonucleotide - Mouse TLR9 ligand

Catalog # tlr1-1826f (formerly tlr1-fmodn)

For research use only

Version # 16J28-MM

### PRODUCT INFORMATION

#### Content

- 50 µg (7.31 nmol) lyophilized ODN 1826 labeled with FITC at the 3' terminus
- 1.5 ml endotoxin-free water

#### ODN 1826 sequence

5'-tccatgacgttctctgacgtt-3' (20 mer)

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

**Molecular weight:** 6901 g/mol

#### Storage

- ODN 1826 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)<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 1826 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- $\alpha$  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 1826 FITC.
- Resuspend lyophilized ODN 1826 FITC with 14 µl of endotoxin-free water (provided).
  - Vortex until completely dissolved.
  - Store at -20 °C.

#### TLR9 stimulation using ODN 1826 FITC

ODN 1826 FITC can be used to stimulate 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.

- 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.
- 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 1826	tlr1-1826
pUNO1-mTLR9 (mouse TLR9 gene)	puno1-mtlr9
HEK-Blue™ mTLR9 Cells	hkb-mtlr9
QUANTI-Blue™	rep-qb1
293/mTLR9 Cells	293-mtlr9
pNiFty2-Luc (Zeo <sup>®</sup> )	pnifty2-luc
pNiFty2-SEAP (Zeo <sup>®</sup> )	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](mailto:info@invivogen.com)

