

# ODN 1826 Biotin

Biotin labeled CpG oligonucleotide, type B; Mouse TLR9 ligand

Catalog # tlr1-1826b

<http://www.invivogen.com/odn1826-biotin>

For research use only

Version # 17A02-MM

## PRODUCT INFORMATION

### Content

- 2 x 50 µg (7.32 nmol) lyophilized ODN 1826 labeled with biotin at the 3' terminus.
- 1.5 ml endotoxin-free water

### ODN 1826 sequence

5'-tccatgacgttcctgacgtt-3' (20 mer)

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

**Molecular weight:** 6824 g/mol

### Storage

- ODN 1826 Biotin is shipped at room temperature. Store at -20°C. Lyophilized product is stable for 12 months.
- Resuspended product should be stored at -20°C. Resuspended 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-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 ODN 1826 Biotin solution (500 µM)

- TLR9 activation can be achieved with 1-5 µM of ODN 1826 Biotin.
1. Resuspend 50 µg of lyophilized ODN 1826 Biotin with 15 µl of endotoxin-free water (provided).
  2. Vortex gently until completely dissolved.
  3. Store at -20°C.

### TLR9 stimulation using ODN 1826 Biotin

ODN 1826 Biotin 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/hek-blue-tlr9](http://www.invivogen.com/hek-blue-tlr9)

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

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

## RELATED PRODUCTS

Product	Catalog Code
ODN 1826	tlrl-1826
ODN 1826 FITC	tlrl-1826f
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)