

# ODN 2007

## Class B CpG oligonucleotide; Bovine / porcine TLR9 ligand

Catalog code: tlr1-2007-1, tlr1-2007-5

<https://www.invivogen.com/odn2007>

For research use only

Version 23K22-MM

## PRODUCT INFORMATION

### Contents

- ODN 2007 is provided lyophilized and is available in two quantities:
  - 1 mg (**141.65 nmol**); tlr1-2007-1
  - 5 x 1 mg (5 mg; **708.25 nmol**); tlr1-2007-5

*Note:* ODN 2007 is sterile filtered prior to lyophilization.

- endotoxin-free water; 1.5 ml with #tlr1-2007-1, and 10 ml with #tlr1-2007-5.

### ODN 2007 sequence

5'- tcg tcg ttg tcg ttt tgt cgt t -3' (22 mer)

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

**Molecular weight:** 7058 g/mol

### Storage and stability

- ODN 2007 is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension, prepare aliquots of ODN 2007 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

- TLR9 activity has been tested using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

ODN 2007 is a Class B CpG oligonucleotide (ODN). It is a short synthetic single-stranded DNA molecule containing unmethylated CpG dinucleotides (CpG motifs). These unmethylated CpG motifs mimic microbial DNA and act as immunostimulants. ODN 2007 is a ligand of choice for bovine and porcine Toll-like receptor 9 (TLR9). Activation of TLR9 triggers NF- $\kappa$ B- and interferon regulatory factor (IRF)-mediated pro-inflammatory responses upon the recognition of unmethylated cytosine-phosphorothioate-guanosine (CpG) forms of DNA<sup>1,2</sup>. Unmethylated CpG dinucleotides are a hallmark of microbial (bacterial, viral, fungal, and parasite) DNA, as well as mitochondrial self-DNA<sup>2,3</sup>. Class B (also called Type K) CpG ODNs, such as ODN 2007, contain a full phosphorothioate backbone with one or more CpG dinucleotides. They strongly activate B cells but weakly stimulate IFN- $\alpha$  secretion in plasmacytoid dendritic cells<sup>4</sup>.

1. Kumagai Y. *et al.*, 2008. TLR9 as a key receptor of the recognition of DNA. *Adv. Drug. Deliv. Rev.* 60(7):795-804. 2. Kayraklioglu N. *et al.*, 2021. CpG oligonucleotides as vaccine adjuvants. *DNA Vaccines: Methods and Protocols*. *Methods in Molecular Biology*. Vol. 2197, p51-77. 3. Kumar V., 2021. The trinity of cGAS, TLR9, and ALRs: guardians of the cellular galaxy against host-derived self-DNA. *Front. Immunol.* 11:624597. 4. Krieg A.M. *et al.*, 1995. CpG motifs in bacterial DNA trigger direct B-cell activation. *Nature.* 374(6522):546-9.

## METHODS

### Preparation of CpG ODN solution (500 $\mu$ M)

TLR9 activation can be achieved with 1-5  $\mu$ M ODN 2007.

- Add 285  $\mu$ l of endotoxin-free water (provided) to 1 mg vial of ODN 2007.
- 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 using ODN 2007

ODN 2007 can be used to stimulate TLR9 in cellular assays. If your cell line does not express TLR9, transfect with a TLR9 plasmid, such as pUNO1-bTLR9 or pUNO1-pTLR9 plasmids expressing the bovine and pig TLR9 genes, respectively. InvivoGen also provides a collection of inducible reporter plasmids for monitoring PRR (pattern recognition receptor) activation and cytokine signaling. Two families of pNiFty plasmids are available pNiFty2-N family (NF- $\kappa$ B-inducible reporter plasmids) and pNiFty3-I family (IRF-inducible reporter plasmids). For more information, visit: [www.invivogen.com/innate-immunity-pnifty](http://www.invivogen.com/innate-immunity-pnifty).

In addition, ODN 2007 can be used to stimulate TLR9 in HEK-Blue™ TLR9 cells. HEK-Blue™ TLR9 cells stably overexpress an NF- $\kappa$ B-inducible secreted embryonic alkaline phosphatase (SEAP) reporter gene and the human TLR9 (hTLR9) or murine TLR9 (mTLR9) gene. For more information, visit: <https://www.invivogen.com/hek-blue-tlr9>.

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

- Dispense 20  $\mu$ l 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™ Solution, a SEAP detection medium.

## RELATED PRODUCTS

Product	Description	Cat. Code
HEK-Blue™ hTLR9 cells	Reporter cells	hkb-htr9
HEK-Blue™ mTLR9 cells	Reporter cells	hkb-mtr9
pNiFty2-N-SEAP-Zeo	Reporter plasmid	prf2-sp
pUNO1-bTLR9	Bovine TLR9 gene	puno1-btr9
pUNO1-pTLR9	Pig TLR9 gene	puno1-ptlr9
QUANTI-Blue™ Solution	SEAP detection medium	rep-qbs

## 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 Asia: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)