

# Human & Mouse TLR2 Agonist Kit

Set of known agonists for human and mouse TLR2

Catalog code: tlr1-kit2hm

<https://www.invivogen.com/tlr2-agonist-kit>

For research use only

Version 18G26-MM

## PRODUCT INFORMATION

### Contents

- TLR6/2 agonist - **Pam2CSK4** (10 µg)
  - TLR1/2 agonist - **Pam3CSK4** (10 µg)
  - TLR6/2 agonist - **FSL-1** (10 µg)
  - TLR2 agonist - **HKLM** (10<sup>9</sup> cells)
  - TLR2 agonist - **LPS-PG** (100 µg)
  - TLR2 agonist - **LTA-SA standard** (100 µg)
  - TLR2 agonist - **PGN-SA** (100 µg)
- 2 x 1.5 ml endotoxin-free water

### Storage and stability

- Products are shipped at room temperature and should be stored according to the table below.

TLR Ligands	Lyophilized	Resuspended
<b>Pam2CSK4</b>	1 year @ 4°C	1 month @ 4°C, 6 months @ -20°C
<b>Pam3CSK4</b>	1 year @ 4°C	1 month @ 4°C, 6 months @ -20°C
<b>FSL-1</b>	1 year @ 4°C	6 months @ 4°C
<b>HKLM</b>	1 year @ 4°C	1 month @ 4°C, 6 months @ -20°C
<b>LPS-PG</b>	1 year @ 4°C	1 month @ 4°C, 6 months @ -20°C
<b>LTA-SA</b>	1 year @ -20°C	1 month @ 4°C, 6 months @ -20°C
<b>PGN-SA</b>	2 years @ -20°C	1 year @ -20°C

## DESCRIPTION

### • Pam2CSK4 - TLR6/2 agonist

Pam2CSK4 is a synthetic diacylated lipopeptide (LP). Bacterial lipoproteins are strong immune modulators that activate early innate host responses after infection. LP analogues of these lipoproteins signal either through TLR1/2 or TLR6/2 heterodimers. According to the current model, diacylated LPs induce signaling through TLR6/2. However, it was reported that diacylated LP, such as Pam2CSK4, induce signaling in a TLR6-independent manner<sup>1</sup>. This finding suggests that both the lipid and peptide part of lipoproteins take part in the specificity of recognition by TLR2 heterodimers.

### • Pam3CSK4 - TLR1/2 agonist

Pam3CSK4 is a synthetic tripalmitoylated lipopeptide that mimicks the acylated amino terminus of bacterial lipoproteins. Pam3CysSerLys4 (Pam3CSK4) is a potent activator of the proinflammatory transcription factor NF-κB<sup>2</sup>. Recognition of Pam3CSK4 is mediated by TLR2 which cooperates with TLR1 through their cytoplasmic domain to induce the signaling cascade leading to the activation of NF-κB<sup>3</sup>.

### • FSL1 - TLR6/2 agonist

FSL-1 (Pam2CGDPKHPKSF) is a synthetic lipoprotein that represents the N-terminal part of the 44-kDa lipoprotein LP44 of *Mycoplasma salivarium*<sup>4</sup>. The framework structure of FSL-1 is the same as that of MALP-2, a *Mycoplasma fermentans* derived lipopeptide (LP), but they differ in the amino acid sequence and length of the peptide portion<sup>5</sup>. FSL-1 is recognized by TLR2 and TLR6 inducing a MyD88-dependent signaling cascade that leads to the activation of NF-κB and the production of proinflammatory cytokines.

### • HKLM - TLR2 agonist

HKLM is a freeze-dried heat-killed preparation of *Listeria monocytogenes* (LM), a facultative intracellular Gram-positive bacterium. Infection with LM induces a strong nonspecific response characterized by the secretion of proinflammatory cytokines. This response is mediated by TLR2<sup>6</sup>. Stimulation with HKLM induces immediate activation of NF-κB and the production of proinflammatory cytokines<sup>7</sup>.

### • LPS-PG - TLR2 agonist

Recognition of LPS from *P. gingivalis* (LPS-PG), a Gram-negative bacteria, is unusual as it appears mediated by either TLR2 and TLR4<sup>8</sup>. Indeed, bone marrow cells obtained from TLR2-/- or TLR4-/- mice respond to LPS-PG while bone marrow cells obtained from TLR2 and TLR4 double-knockout do not. LPS-PG has also been reported to act as a TLR4 antagonist in some cell types<sup>9,10</sup>. This discrepancy may be explained by the ability of this bacterium to synthesize multiple, structurally different forms of lipid A. The TLR response to LPS-PG is dependent on the presence of key accessory molecules: CD14 is required for both TLR2 and TLR4 activation while MD-2 is only necessary for TLR4 activation<sup>8</sup>.

## 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 3622-3480

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

#### • LTA-SA - TLR2 agonist

LTA-SA is lipoteichoic acid (LTA) from *S.aureus* (SA). LTA is a major immunostimulatory component of Gram-positive bacteria. LTA is responsible for causing gram-positive sepsis. Like LPS, LTA is an amphiphile formed by a hydrophilic polyphosphate polymer linked to a neutral glycolipid. LTA stimulates immune cells through TLR2 to produce TNF- $\alpha$  and other inflammatory cytokines<sup>11</sup>. Recognition of LTA also involves LPS-binding protein (LBP) and CD14 but not TLR4 or MD2<sup>12</sup>. Activation of LTAs may require the involvement of TLR1<sup>13</sup>.

#### • PGN-SA - TLR2 agonist

PGN-SA is peptidoglycan (PGN) from *S.aureus* (SA). PGN is a major surface component of Gram-positive bacteria. It is embedded in a relatively thick cell wall and is usually covalently attached to other polymers, such as lipoproteins and LTAs. In Gram-negative bacteria, a thin layer of PGN is also found in the periplasmic space. PGN is known to be a potent activator of NF- $\kappa$ B and TNF- $\alpha$  through TLR2<sup>14</sup>. However, other pattern recognition proteins have been reported to mediate the immunostimulatory activity of PGN<sup>15, 16, 17</sup>. This discrepancy is correlated to the method of purification. PGN-SA which is purified by detergent lysis, enzymatic treatment, LiCl/EDTA and acetone cleaning is an activator of TLR2.

## METHODS

### Preparation of TLR agonist stock solutions

Product	Working concentration	Stock solution concentration	Volume of solvent
Pam2CSK4	1-100 ng/ml	100 $\mu$ g/ml	100 $\mu$ l H <sub>2</sub> O
Pam3CSK4	1-300 ng/ml	100 $\mu$ g/ml	100 $\mu$ l H <sub>2</sub> O
FSL-1	1-100 ng/ml	100 $\mu$ g/ml	100 $\mu$ l H <sub>2</sub> O
HKLM	10 <sup>7</sup> -10 <sup>8</sup> cells/ml	10 <sup>10</sup> cells/ml	100 $\mu$ l H <sub>2</sub> O
LPS-PG	100 ng-10 $\mu$ g/ml	1 mg/ml	100 $\mu$ l H <sub>2</sub> O
LTA-SA	100 ng-1 $\mu$ g/ml	200 $\mu$ g/ml	500 $\mu$ l H <sub>2</sub> O
PGN-SA	1-10 $\mu$ g/ml	200 $\mu$ g/ml	500 $\mu$ l H <sub>2</sub> O

### TLR stimulation

- Transfect your cell line with an NF- $\kappa$ B-inducible reporter plasmid, i.e. a plasmid carrying a reporter gene, such as SEAP or luciferase, under the control of an NF- $\kappa$ B-inducible ELAM-1 (E-selectin) promoter<sup>18</sup>.

*Note: InvivoGen provides pNiFty, a family of NF- $\kappa$ B-inducible reporter plasmids that can be transfected transiently (pNiFty) or stably (pNiFty2). pNiFty plasmids are available either with the SEAP or luciferase reporter genes.*

If your cell line does not naturally express TLRs, cotransfect with a plasmid expressing a given TLR gene, such as the pUNO plasmid family.

*Note: Alternatively, evaluate TLR2 activation using reporter cells, such as InvivoGen's HEK-Blue™ TLR2 cells which express the human and mouse TLR2 and SEAP reporter genes. NF- $\kappa$ B production in these cells can be easily quantified using a SEAP detection medium, such as QUANTI-Blue™ or HEK-Blue™ Detection.*

- Twenty-four to forty-eight hours after transfection, stimulate cells with the corresponding agonist for 6 hours to 24 hours.

- Determine TLR stimulation by assessing reporter gene expression using the appropriate detection system.

## References

1. Buwitt-Beckmann U. *et al.*, 2005. Toll-like receptor 6-independent signaling by diacylated lipopeptides. *Eur J Immunol.* 35(1):282-9.
2. Aliprantis AO. *et al.*, 1999. Cell activation and apoptosis by bacterial lipoproteins through toll-like receptor-2. *Science.* 285(5428):736-9.
3. Ozinsky A. *et al.*, 2000. The repertoire for pattern recognition of pathogens by the innate immune system is defined by cooperation between toll-like receptors. *PNAS.* 97(25):13766-71.
4. Shibata KI. *et al.*, 2000. The N-terminal lipopeptide of a 44-kDa membrane-bound lipoprotein of *Mycoplasma salivarium* is responsible for the expression of intercellular adhesion molecule-1 on the cell surface of normal human gingival fibroblasts. *J. Immunol.* 165:6538-6544.
5. Okusawa T. *et al.* 2004. Relationship between Structures and Biological Activities of Mycoplasmal Diacylated Lipopeptides and Their Recognition by Toll-Like Receptors 2 and 6. *Infect Immun.* 72(3): 1657-1665.
6. Flo TH. *et al.*, 2000. Human toll-like receptor 2 mediates monocyte activation by *Listeria monocytogenes*, but not by group B streptococci or lipopolysaccharide. *J Immunol.* 164(4):2064-9
7. Hauf N. *et al.*, 1997. *Listeria monocytogenes* infection of P388D1 macrophages results in a biphasic NF- $\kappa$ B (RelA/p50) activation induced by lipoteichoic acid and phospholipases and mediated by I $\kappa$ B $\alpha$  and I $\kappa$ B $\beta$  degradation. *Proc. Natl. Acad. Sci. USA* 94(17):9394-9
8. Darveau RP. *et al.*, 2004. Porphyromonas gingivalis lipopolysaccharide contains multiple lipid A species that functionally interact with both toll-like receptors 2 and 4. *Infect Immun.* 72(9):5041-51.
9. Yoshimura A. *et al.*, 2002. Lipopolysaccharides from periodontopathic bacteria Porphyromonas gingivalis and Capnocytophaga ochracea are antagonists for human toll-like receptor 4. *Infect Immun.* 70(1):218-25.
10. Coats SR. *et al.*, 2003. Porphyromonas gingivalis lipopolysaccharide antagonizes Escherichia coli lipopolysaccharide at toll-like receptor 4 in human endothelial cells. *Infect Immun.* 71(12):6799-807.
11. Schwandner R. *et al.*, 1999. Peptidoglycan- and lipoteichoic acid-induced cell activation is mediated by toll-like receptor 2. *J Biol Chem.* 274(25):17406-9.
12. Schroder NW. *et al.*, 2003. Lipoteichoic acid (LTA) of *Streptococcus pneumoniae* and *Staphylococcus aureus* activates immune cells via Toll-like receptor (TLR)-2, lipopolysaccharide-binding protein (LBP), and CD14, whereas TLR-4 and MD-2 are not involved. *J Biol Chem.* 278(18):15587-94.
13. Han SH. *et al.*, 2003. Pneumococcal lipoteichoic acid (LTA) is not as potent as staphylococcal LTA in stimulating Toll-like receptor 2. *Infect Immun.* 71(10):5541-8.
14. Takeuchi O. *et al.*, 1999. Differential roles of TLR2 and TLR4 in recognition of gram-negative and gram-positive bacterial cell wall components. *Immunity.* 11(4):443-51.
15. Travassos LH. *et al.*, 2004. Toll-like receptor 2-dependent bacterial sensing does not occur via peptidoglycan recognition. *EMBO Rep.* 5(10):1000-1006.
16. Girardin SE *et al.*, 2003. Peptidoglycan molecular requirements allowing detection by Nod1 and Nod2. *J Biol Chem.* 278(43):41702-8.
17. Dziarski R., 2004. Peptidoglycan recognition proteins (PGRPs). *Mol Immunol.* 40(12):877-86.
18. Schindler U. & Baichwal VR., 1994. Three NF- $\kappa$ B binding sites in the human E-selectin gene required for maximal tumor necrosis factor alpha-induced expression. *Mol Cell Biol.* 14(9):5820-5831.

## 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 3622-3480  
E-mail: info@invivogen.com

