

# TL2-C29

TLR2 signaling inhibitor

Catalog code: inh-c29

<https://www.invivogen.com/tlr2-in-c29>

For research use only

Version 22D11-NJ

## PRODUCT INFORMATION

### Contents

TL2-C29 is provided as a dried powder:

- 5 mg: inh-c29

### Storage and stability

- TL2-C29 is shipped at room temperature. Upon receipt, store at -20 °C.
- Upon resuspension, prepare aliquots and store at -20 °C. Resuspended product is stable for 2 months at -20 °C when properly stored. Avoid repeated freeze-thaw cycles.

### Quality control

- Purity: ≥95% (NMR)
- The inhibitory activity of human TLR2/1 and TLR2/6 signaling has been confirmed 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

TL2-C29 is a small-molecule inhibitor of Toll-like receptor 2 (TLR2)<sup>1</sup>. TLR2 plays an essential role in detecting a diverse range of microbial pathogen-associated molecular patterns (PAMPs) from bacteria, fungi, parasites, and viruses<sup>2</sup>. TLR1 and TLR6 co-receptors are crucial for TLR2 signaling and ligand specificity. Cell surface TLR2/TLR1 and TLR2/TLR6 heterodimers bind tri- and diacylated lipoproteins, respectively<sup>3,4</sup>. TLR2 signaling is initiated by ligand-induced dimerization of the essential cytoplasmic TIR (Toll/interleukin-1 receptor) domains of the TLR2 heterodimers and adapter proteins. Subsequent signaling cascades lead to a pro-inflammatory response.

TL2-C29 binds a pocket in the BB loop within the TLR2 TIR domain, thereby inhibiting TLR2 interaction with the MyD88 adapter molecule and downstream MAPK and NF-κB activation<sup>1</sup>. In human TLR2 signaling, TL2-C29 blocks both TLR2/1 and TLR2/6 pathways. In murine TLR2 signaling, TL2-C29 preferentially blocks the TLR2/1 pathway<sup>1</sup>.

1. Mistry, P. *et al.*, 2015. Inhibition of TLR2 signaling by small molecule inhibitors targeting a pocket within the TLR2 TIR domain. *PNAS*. 112(17):5455-5460. 2. Oliveira-Nascimento L. *et al.*, 2012. The role of TLR2 in infection and Immunity. *Front Immunol*. 3(79): doi:10.3389/fimmu.2012.00079. 3. Takeuchi O. *et al.*, 2001. Discrimination of bacterial lipoproteins by Toll-like receptor 6. *Int Immunol*. 13: 933-940. 4. Takeuchi O. *et al.*, 2002. Cutting edge: role of Toll-like receptor 1 in mediating immune response to microbial lipoproteins. *J Immunol*. 169: 10-14.

## CHEMICAL PROPERTIES

**Synonyms:** C29; TLR2-IN-C29; 3-[[[(2-hydroxy-3-methoxyphenyl)methylene]amino]-2-methyl-benzoic acid

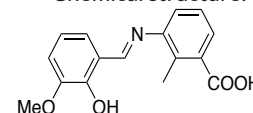
**CAS number:** 363600-92-4

**Formula:** C<sub>16</sub>H<sub>15</sub>NO<sub>4</sub>

**Molecular weight:** 285.3 g/mol

**Solubility:** 50 mM (14.3 mg/ml) in DMSO

**Chemical structure:**



## METHODS

### Preparation of stock solution (50 mM)

1. Add 350 µl DMSO to 5 mg TL2-C29.
2. Vortex until completely resuspended.
3. Prepare aliquots of TL2-C29 and store at -20 °C.
4. Once TL2-C29 is resuspended, further dilutions can be prepared using sterile saline buffers (e.g. PBS). *Note:* TL2-C29 is not soluble in water.

### Working concentration range in cellular assays:

- 25 µM - 200 µM for hTLR2/1 signaling inhibition
- 10 µM - 200 µM for hTLR2/6 signaling inhibition

## PROTOCOL

Below is a protocol using InvivoGen's HEK-Blue™ hTLR2 cells for studying the specific inhibition of human TLR2 signaling by TL2-C29. These cells express an inducible secreted embryonic alkaline phosphatase (SEAP) reporter to readily measure the activation of the NF-κB pathway. Changes in SEAP expression due to inhibition of TLR2 signaling can be assessed using QUANTI-Blue™ Solution, a SEAP detection reagent.

1. Add 20 µl TL2-C29 (10x conc) per well of a flat bottom 96-well plate.
2. Prepare a suspension of HEK-Blue™ hTLR2 cells (~310,000 cells per ml) in culture medium.
3. Add 160 µl of the cell suspension (~50,000 cells) to each well.
4. Incubate the plate at 37 °C in a CO<sub>2</sub> incubator for 3 hours.
5. Add 20 µl (10x conc) of an inducer of TLR2 signaling (e.g. Pam3CSK4 or FSL-1) and incubate the plate at 37 °C in a CO<sub>2</sub> incubator for 24 hours.
6. Prepare QUANTI-Blue™ Solution and carry out the measurements following the instructions on the data sheet.

## RELATED PRODUCTS

Product	Cat. Code
HEK-Blue™ hTLR2 cells	hkb-htlr2
QUANTI-Blue™ Solution	rep-qbs
Pam3CSK4	tlrl-pms
FSL-1	tlrl-fsl

## 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 3622-3480

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

 **InvivoGen**  
www.invivogen.com