

# CL572

## Dual TLR2 (human only) & TLR7 ligand

Catalog # tlrl-c572

For research use only

Version # 15B04-MM

### PRODUCT INFORMATION

#### Content:

- 500 µg CL572 provided as a lyophilized powder

#### Storage:

- CL572 is shipped at room temperature. Store lyophilized product at -20 °C. Lyophilized product is stable for 1 year at -20 °C.
- Upon resuspension, store at -20 °C. Resuspended product is stable for 6 months at 4 °C. Do not store resuspended product in plastic tubes.

### DESCRIPTION

CL572 is a 8-hydroxy-adenine compound conjugated to a monoacyl-ethyl-cystein group via a glutamic acid derivative. Monoacyl-ethyl-cystein-containing dipeptides have been recently shown to specifically activate human TLR2<sup>1</sup>. Indeed, CL572 is a robust inducer of human TLR2 but is unable to stimulate mouse TLR2. CL572 is also a potent inducer of TLR7.

1. Agnihotri G. *et al.*, 2011. Structure-activity relationships in toll-like receptor 2-agonists leading to simplified monoacyl lipopeptides. *J Med Chem.* 54(23):8148-60.

### BACKGROUND

InvivoGen has developed a series of novel molecules designed to induce potent immune responses through the combined activation of several pattern recognition receptors (PRRs) that trigger different innate immune signaling pathways. These molecules are agonists for TLR2, TLR7 or both. Agonists that activate TLR2 are derived from the well-established TLR2 ligand, Pam2CSK4, and those recognized by TLR7 are derived from the 8-hydroxyadenine derivative CL264, a TLR7 agonist recently developed by InvivoGen (see Related Products overleaf).

TLR2 and TLR7 are two PRRs with distinct characteristics. TLR2 is a cell surface receptor expressed by many cell types, while TLR7 is an endosomal receptor expressed predominantly in plasmacytoid dendritic cells (pDC) and to a lesser extent in B cells. TLR2 signaling triggers the NF-κB pathway and the production of pro-inflammatory cytokines, such as TNF-α, whereas TLR7 signaling induces mainly the IRF pathway and the production of IFN-α. Combined activation of these different pathways results in robust immune responses with potential therapeutic effects. InvivoGen's multi-PRR agonists are promising candidates for antitumor and vaccine applications.

### APPLICATIONS

CL572 can be used to stimulate both TLR2 and TLR7.

### CHEMICAL PROPERTIES

**Synonym:** S-(2-myristoyloxy ethyl)-(R)-cysteinyl 4-((6-amino-2-(butylamino)-8-hydroxy-9H-purin-9-yl)methyl) aniline

**Formula:** C<sub>41</sub>H<sub>65</sub>N<sub>9</sub>O<sub>7</sub>S

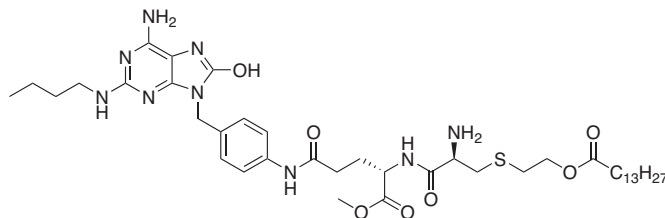
**Molecular weight:** 828 g/mol

**Solubility:** DMSO (1 mg/ml)

**Working concentration:** 0.5 ng - 1 µg/ml (~1 nM - 10 µM)

**Endotoxin level:** <0.001 EU/µg

#### Structure:



### METHODS

#### Preparation of CL572 stock solution (1 mg/ml)

- Add 500 µl DMSO to 500 µg CL572. Vortex until complete solubilization.
- Further dilutions can be prepared using aqueous solutions.

#### TLR stimulation with CL572 using HEK-Blue cells

CL572 can be used to stimulate TLR2 in HEK-Blue™ TLR2 cells and TLR7 in HEK-Blue™ TLR7 cells. These cells stably express an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP) and overexpress the appropriate TLR gene. For more information visit: [www.invivogen.com/hek-blue-cells](http://www.invivogen.com/hek-blue-cells)

1. Stimulate HEK-Blue™ TLR2 cells with 0.5 ng - 1 µg/ml CL572 and HEK-Blue™ TLR7 cells with 100 ng - 10 µg/ml CL572.
2. Incubate for 6 - 24 h at 37 °C, 5% CO<sub>2</sub>.
3. Determine TLR stimulation using a SEAP detection medium, such as QUANTI-Blue™ or HEK-Blue™ Detection (see Related Products, overleaf) or by assessing cytokine expression using an ELISA.

#### 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)

## RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ hTLR2 Cells	hkb-htlr2
HEK-Blue™ mTLR2 Cells	hkb-mtlr2
HEK-Blue™ hTLR7 Cells	hkb-htlr7
HEK-Blue™ mTLR7 Cells	hkb-mtlr7
RAW-Blue™ Cells (Mouse macrophage reporter cells)	raw-sp
HEK-Blue™ Detection (SEAP detection medium)	hb-det2
QUANTI-Blue™ (SEAP detection medium)	rep-qb1
<b>Dual TLR2 &amp; TLR7 ligands:</b>	
CL401 (TLR2 & TLR7 ligand)	tlrl-c401
Adilipoline™ (CL413; TLR2 & TLR7 ligand)	tlrl-c413
CL531 (TLR2 & TLR7 ligand)	tlrl-c531
<b>TLR ligands &amp; nucleic acid carriers:</b>	
AdiFectin™ (CL347; TLR7 ligand & nucleic acid carrier)	tlrl-c347
CL419 (TLR2 ligand & nucleic acid carrier)	tlrl-c419
PamadiFectin™ (CL553; TLR7 & TLR2 ligand & nucleic acid carrier)	tlrl-c553
<b>TLR2 ligands:</b>	
HKLM (Heat killed <i>Listeria monocytogenes</i> )	tlrl-hklm
FSL-1 (Synthetic diacylated lipoprotein)	tlrl-fsl
Pam2CSK4 (Synthetic diacylated lipoprotein)	tlrl-pm2s
<b>TLR7 ligands:</b>	
CL264 (8-hydroxyadenine derivative)	tlrl-c264e
CL307 (Spermine covalently linked to CL264 )	tlrl-c307
Gardiquimod™ (Imidazoquinoline compound)	tlrl-gdqs

---

### 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)