

# CL097

## Imidazoquinoline Compound - TLR7/8 ligand

Catalog # tlr-c97, tlr-c97-5

For research use only

Version # 12F13-MM

### PRODUCT INFORMATION

#### Content:

- CL097 is provided lyophilized and is available in two quantities:
  - 500 µg: tlr-c97
  - 5 mg: tlr-c97-5
- sterile endotoxin-free water, 1.5 ml with #tlr-c97 and 10 ml with #tlr-c97-5

#### Storage:

- CL097 is provided as a pale yellow solid and shipped at room temperature. Store at -20°C. Lyophilized product is stable 1 year at -20°C when properly stored.
- Upon resuspension, prepare aliquots of CL097 and store at -20°C. Resuspended product is stable 6 months at -20°C. Avoid repeated freeze-thaw cycles.

### DESCRIPTION

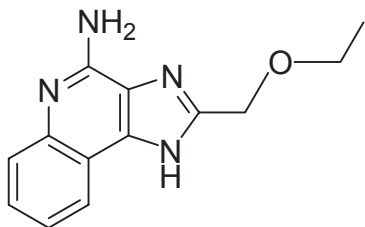
CL097 is a highly water-soluble derivative of the imidazoquinoline compound R848 ( $\geq 20$  mg/ml). Similarly to R848, CL097 is a TLR7 and TLR8 ligand<sup>1,2</sup>. It induces the activation of NF-κB at 0.4 µM (0.1 µg/ml) in TLR7-transfected HEK293 cells and at 4 µM (1 µg/ml) in TLR8-transfected HEK293 cells.

#### References

1. **Salio M. et al., 2007.** Modulation of human natural killer T cell ligands on TLR-mediated antigen-presenting cell activation. PNAS 104: 20490 - 20495. 2. **Butchi N.J. et al., 2008.** Analysis of the Neuroinflammatory Response to TLR7 Stimulation in the Brain: Comparison of Multiple TLR7 and/or TLR8 Agonists. J Immunol 180: 7604-7612. 3. **Schindler U. & Baichwal VR., 1994.** Three NF-κB binding sites in the human E-selectin gene required for maximal tumor necrosis factor alpha-induced expression. Mol Cell Biol, 14(9):5820-5831.

**Formula:** C<sub>13</sub>H<sub>14</sub>N<sub>4</sub>O

**Molecular weight:** 242.28



### METHODS

#### Preparation of CL097 stock solution (1 mg/ml)

- Stimulation of TLR7 can be achieved with 50 ng - 5 µg/ml CL097, and stimulation of TLR8 with 0.5 - 5 µg/ml CL097.
- Resuspend CL097 with sterile endotoxin-free water provided.
  - To obtain a 1 mg/ml stock solution:
    - Add 500 µl to 500 µg vial
    - Add 5 ml to 5 mg vial
  - Vortex until complete solubilization.
  - Prepare aliquots of CL097 and store at -20°C.

#### CL097 stimulation

Transfect your cell line with a pNiFty plasmid, an NF-κB reporter plasmid, i.e. a plasmid carrying a reporter gene such as SEAP or luciferase, under the control of an NF-κB-inducible ELAM-1 (E-selectin) promoter<sup>3</sup>.

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

- If your cell line does not naturally express TLR7 or TLR8, cotransfect with a TLR7 or TLR8 expressing plasmid, such as pUNO1-hTLR7 and pUNO1-hTLR8b.

*Note: Alternatively, evaluate TLR7 or TLR8 activation using reporter cells, such as InvivoGen's HEK-Blue™ hTLR7 or hTLR8 cells which express the human TLR7 or TLR8 gene and a SEAP reporter gene. NF-κB 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 50 ng - 5 µg/ml CL097 for 6 to 24 hours.
- Determine CL097 stimulation on TLR7 or TLR8 by assessing reporter gene expression using the appropriate detection system.

### RELATED PRODUCTS

Product	Catalog Code
HEK-Blue™ hTLR7 cells	hkb-htr7
HEK-Blue™ hTLR8 cells	hkb-htr8
pUNO1-hTLR7	puno1-htr7
pUNO1-hTLR8b	puno1-htr8b
pNiFty-Luc (Amp <sup>R</sup> )	pnifty-luc
pNiFty-SEAP (Amp <sup>R</sup> )	pnifty-seap
pNiFty2-Luc (Zeo <sup>R</sup> )	pnifty2-luc
pNiFty2-SEAP (Zeo <sup>R</sup> )	pnifty2-seap

#### 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

