

CL264

Adenine analog; TLR7 ligand

Catalog code: tlrl-c264e, tlrl-c264e-5

<https://www.invivogen.com/cl264>

For research use only

Version 19K29-MM

PRODUCT INFORMATION

Contents

• CL264 is provided as a lyophilized powder containing Tris-acetate-EDTA buffer salts. CL264 is available in two quantities:

- 500 µg CL264: tlrl-c264e

- 5 mg CL264: tlrl-c264e-5

• Sterile endotoxin-free water, 1.5 ml with #tlrl-c264e and 10 ml with #tlrl-c264e-5.

Storage and stability

- CL264 is shipped at room temperature. Store lyophilized product at -20°C. Lyophilized product is stable for 1 year at -20°C.

- Upon resuspension, prepare aliquots of CL264 and store at -20°C. Resuspended product is stable for 6 months at -20°C. Avoid repeated freeze-thaw cycles.

Quality control

- The biological activity has been validated 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

CL264 is a 9-benzyl-8-hydroxyadenine derivative containing a glycine on the benzyl group (in para). Similarly to SM360320, CL264 induces the activation of NF-κB and the secretion of IFN-α in TLR7-expressing cells¹. CL264 is a TLR7-specific ligand, it does not stimulate TLR8 even at high concentrations (> 10 µg/ml). In TLR7-transfected HEK293 cells, CL264 triggers NF-κB activation at a concentration of 0.1 µM which is 5-10 times less than Imiquimod.

1. Lee J. et al., 2006. Activation of anti-hepatitis C virus responses via Toll-like receptor 7. PNAS 103(6):1828-33.

CHEMICAL PROPERTIES

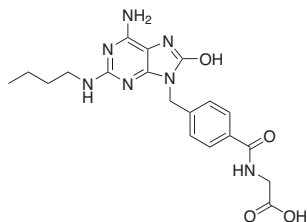
Synonym: 2-(4-((6-amino-2-(butylamino)-8-hydroxy-9H_purin_9-yl)methyl)benzamido)acetic acid

Formula: C₁₉H₂₃N₇O₄

Molecular weight: 413.43 g/mol

Solubility: 1 mg/ml in water

Structure:



METHODS

Preparation of CL264 stock solution (1 mg/ml)

Stimulation of TLR7 can be achieved with 50 ng - 10 µg/ml CL264.

To obtain a 1 mg/ml stock solution:

- Add 500 µl sterile water (provided) to 500 µg of CL264.

- Add 5 ml sterile water (provided) to 5 mg of CL264.

This generates a stock solution at 1 mg/ml with 40 mM Tris, 20 mM acetic acid and 0.25 mM EDTA. Vortex until complete solubilization. Once CL264 is solubilized, prepare aliquots and store at -20°C.

An example of TLR7 stimulation using CL264

CL264 can be used to stimulate TLR7 in HEK-Blue™ TLR7 cells. These cells stably express an NF-κB-inducible secreted embryonic alkaline phosphatase (SEAP) and overexpress the TLR7 gene. For more information visit: <https://www.invivogen.com/hek-blue-trl7>.

1. Add a range of concentrations of CL264 (50 ng - 10 µg/ml final concentration) to HEK-Blue™ TLR7 cells (prepare a cell suspension according to data sheet).
2. Incubate cells and CL264 for 6-24 h at 37°C, 5% CO₂.
3. Determine TLR7 stimulation with CL264 by assessing SEAP expression using a SEAP detection reagent, such as QUANTI-Blue™ Solution or HEK-Blue™ Detection.

RELATED PRODUCTS

Product	Description	Cat. Code
Gardiquimod™	TLR7 ligand	tlrl-gdq
HEK-Blue™ hTLR7 cells	Human TLR7 reporter cells	hkb-htrl7
HEK-Blue™ mTLR7 cells	Murine TLR7 reporter cells	hkb-mtrl7
HEK-Blue™ Detection	SEAP detection reagent	hb-det2
Imiquimod	TLR7 ligand	tlrl-imq
Loxoribine	TLR7 ligand	tlrl-lox
QUANTI-Blue™ Solution	SEAP detection reagent	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 Hong Kong: +852 3622-3480

E-mail: info@invivogen.com