

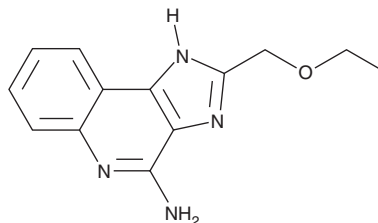
PRODUCT INFORMATION



CL097

Item No. 39513

CAS Registry No.: 1026249-18-2
Formal Name: 2-(ethoxymethyl)-3H-imidazo[4,5-c]quinolin-4-amine
MF: C₁₃H₁₄N₄O
FW: 242.3
Purity: ≥98%
UV/Vis.: λ_{max}: 248, 322 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CL097 is supplied as a solid. A stock solution may be made by dissolving the CL097 in the solvent of choice, which should be purged with an inert gas. CL097 is soluble in DMSO.

Description

CL097 is a dual agonist of toll-like receptor 7 (TLR7) and TLR8.¹ It induces activation of NF-κB in HEK293 cells expressing human TLR7 or TLR8 when used at concentrations of 0.1 and 1 μg/ml, respectively. CL097 induces TNF-α production and enhances LPS- or hyaluronic acid-induced TNF-α, IL-10, and IL-12 production in isolated human monocyte-derived macrophages.² *In vivo*, CL097 (5 mg/kg) increases pancreatic lymph node IFN-α levels in pre-diabetic non-obese diabetic (NOD) mice.³

References

- Schön, M.P. and Schön, M. TLR7 and TLR8 as targets in cancer therapy. *Oncogene* **27(2)**, 190-199 (2008).
- Petricevic, B., Wessner, B., Sachet, M., *et al.* CL097, a TLR7/8 ligand, inhibits TLR-4-dependent activation of IRAK-M and BCL-3 expression. *Shock* **32(5)**, 484-490 (2009).
- Lee, A.S., Ghoreishi, M., Cheng, W.K., *et al.* Toll-like receptor 7 stimulation promotes autoimmune diabetes in the NOD mouse. *Diabetologia* **54(6)**, 1407-1416 (2011).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/03/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM