

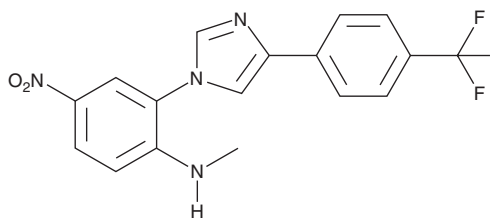
PRODUCT INFORMATION



CU-T12-9

Item No. 32867

CAS Registry No.: 1821387-73-8
Formal Name: N-methyl-4-nitro-2-[4-[4-(trifluoromethyl)phenyl]-1H-imidazol-1-yl]-benzenamine
MF: C₁₇H₁₃F₃N₄O₂
FW: 362.3
Purity: ≥98%
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

CU-T12-9 is supplied as a solid. A stock solution may be made by dissolving the CU-T12-9 in the solvent of choice, which should be purged with an inert gas. CU-T12-9 is soluble in the organic solvent DMSO. The solubility of CU-T12-9 in DMSO is approximately 30 mg/ml.

Description

CU-T12-9 is an agonist of the toll-like receptor 1/2 (TLR1/2) heterodimer.¹ It is selective for TLR1/2 over the TLR2/6 heterodimer in a secreted alkaline phosphatase (SEAP) assay at concentrations up to 10 µg/ml using HEK-Blue cells expressing human TLR2 and endogenously expressing both TLR1 and TLR6. CU-T12-9 (5 µM) induces NF-κB signaling in a reporter assay. It also increases nitric oxide (NO) production in RAW 264.7 cells and primary rat macrophages at 3.6 and 0.4 µM, respectively, and TNF-α production in RAW 264.7 cells (EC₅₀ = 0.06 µM).

Reference

1. Cheng, K., Gao, M., Godfroy, J.I., *et al.* Specific activation of the TLR1-TLR2 heterodimer by small-molecule agonists. *Sci. Adv.* **1(3)**, e1400139 (2015).

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, 11/29/2022

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