

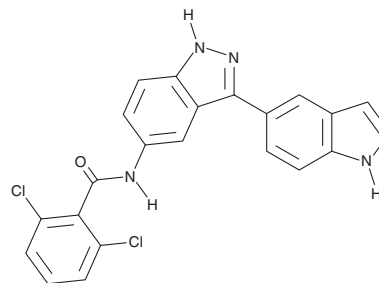
PRODUCT INFORMATION



MD2-TLR4-I

Item No. 38328

CAS Registry No.: 2249801-12-3
Formal Name: 2,6-dichloro-N-[3-(1H-indol-5-yl)-1H-indazol-5-yl]-benzamide
Synonym: Myeloid Differentiation Factor 2-Toll-like Receptor 4-I
MF: C₂₂H₁₄Cl₂N₄O
FW: 421.3
Purity: ≥98%
UV/Vis.: λ_{max}: 242 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

MD2-TLR4-I is supplied as a solid. A stock solution may be made by dissolving the MD2-TLR4-I in the solvent of choice, which should be purged with an inert gas. MD2-TLR4-I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of MD2-TLR4-I in these solvents is approximately 10 mg/ml.

Description

MD2-TLR4-I is a toll-like receptor 4 (TLR4) antagonist.¹ It inhibits LPS-induced production of TNF-α and IL-6 in primary mouse macrophages (IC₅₀s = 0.89 and 0.53 μM, respectively). MD2-TLR4-I inhibits LPS-induced, TLR4-mediated NF-κB activation in a reporter assay using RAW 264.7 cells. *In vivo*, MD2-TLR4-I (10 mg/kg) reduces lung macrophage infiltration and prevents lung edema and alveolar thickening in a mouse model of LPS-induced acute lung injury.

Reference

1. Liu, Z., Chen, L., Yu, P., *et al.* Discovery of 3-(indol-5-yl)-indazole derivatives as novel myeloid differentiation protein 2/Toll-like receptor 4 antagonists for treatment of acute lung injury. *J. Med. Chem.* **62**(11), 5453-5469 (2019).

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, 05/31/2023

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