

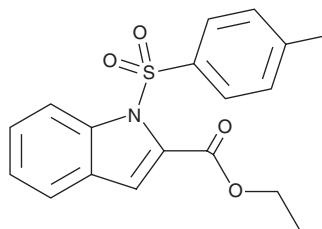
PRODUCT INFORMATION



NOD-IN-1

Item No. 28370

CAS Registry No.: 132819-92-2
Formal Name: 1-[(4-methylphenyl)sulfonyl]-1H-indole-2-carboxylic acid, ethyl ester
MF: C₁₈H₁₇NO₄S
FW: 343.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 271 nm
Supplied as: A crystalline 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

NOD-IN-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the NOD-IN-1 in the solvent of choice, which should be purged with an inert gas. NOD-IN-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NOD-IN-1 in these solvents is approximately 30 mg/ml.

NOD-IN-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NOD-IN-1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NOD-IN-1 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NOD-IN-1 is an inhibitor of nucleotide-binding oligomerization domain (NOD) receptors and a derivative of nodinitib-1 (Item No. 11040).¹ NOD-IN-1 inhibits C12-iE-DAP-induced NF-κB activation in HEK-Blue cells expressing human recombinant NOD1 or NOD2 (IC₅₀s = 5.74 and 6.45 μM, respectively).

Reference

1. Keček Plešec, K., Urbančič, D., Gobec, M., *et al.* Identification of indole scaffold-based dual inhibitors of NOD1 and NOD2. *Bioorg. Med. Chem.* **24**(21), 5221-5234 (2016).

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, 10/25/2022

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