

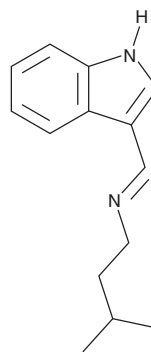
PRODUCT INFORMATION



Indolimine-214

Item No. 37953

Formal Name: (E)-1-(1H-indol-3-yl)-N-isopentylmethanimine
Synonym: Isoamylindolimine
MF: C₁₄H₁₈N₂
FW: 214.3
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 260, 272, 293 nm
Supplied as: A neat oil
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Indolimine-214 is supplied as a neat oil. A stock solution may be made by dissolving the indolimine-214 in the solvent of choice, which should be purged with an inert gas. Indolimine-214 is soluble in the organic solvent methanol.

Description

Indolimine-214 is a metabolite of *M. morganii*, a bacterium that is increased in the gut microbiome of patients with inflammatory bowel disease (IBD) and colorectal cancer (CRC).¹ Indolimine-214 (25 and 100 µg/ml) increases γH2AX levels in HeLa cells and increases the tail DNA percentage in a comet assay using HeLa cells when used at a concentration of 100 µg/ml, both markers of DNA damage, but does not induce DNA damage when used alone in a circular pUC19 plasmid-based genotoxicity assay.

Reference

1. Cao, Y., Oh, J., Xue, M., *et al.* Commensal microbiota from patients with inflammatory bowel disease produce genotoxic metabolites. *Science* **378**(6618), eabm3233 (2022).

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, 01/05/2023

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