

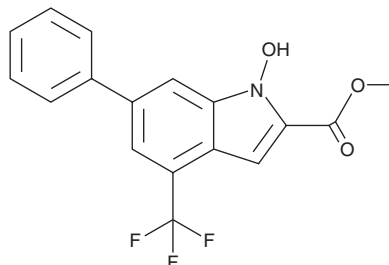
PRODUCT INFORMATION



NHI 2

Item No. 30201

CAS Registry No.: 1269802-97-2
Formal Name: 1-hydroxy-6-phenyl-4-(trifluoromethyl)-1H-indole-2-carboxylic acid, methyl ester
MF: C₁₇H₁₂F₃NO₃
FW: 335.3
Purity: ≥98%
UV/Vis.: λ_{max}: 256, 322 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

NHI 2 is supplied as a crystalline solid. A stock solution may be made by dissolving the NHI 2 in the solvent of choice, which should be purged with an inert gas. NHI 2 is soluble in organic solvents such as ethanol and DMSO. The solubility of NHI 2 in these solvents is approximately 25 mg/ml.

Description

NHI 2 is an inhibitor of lactate dehydrogenase A (LDHA; IC₅₀ = 14.7 μM).¹ It is selective for LDHA over LDHB (IC₅₀ = 55.8 μM). It inhibits lactate production in HeLa cells when used at concentrations of 100 and 200 μM. NHI 2 is cytotoxic to HeLa cells (IC₅₀ = 33.4 μM).

Reference

1. Granchi, C., Calvaresi, E.C., Tuccinardi, T., *et al.* Assessing the differential action on cancer cells of LDH-A inhibitors based on the *N*-hydroxyindole-2-carboxylate (NHI) and malonic (Mal) scaffolds. *Org. Biomol. Chem.* **11(38)**, 6588-6596 (2013).

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, 05/04/2020

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