

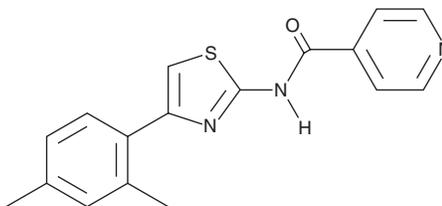
# PRODUCT INFORMATION



## N-[4-(2,4-Dimethylphenyl)-2-thiazolyl]-4-pyridinecarboxamide

Item No. 36050

**CAS Registry No.:** 560103-80-2  
**Formal Name:** N-(4-(2,4-dimethylphenyl)thiazol-2-yl)isonicotinamide  
**Synonym:** Nec2/Hec1 Inhibitor  
**MF:** C<sub>17</sub>H<sub>15</sub>N<sub>3</sub>OS  
**FW:** 309.4  
**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

N-[4-(2,4-Dimethylphenyl)-2-thiazolyl]-4-pyridinecarboxamide is supplied as a solid. A stock solution may be made by dissolving the N-[4-(2,4-dimethylphenyl)-2-thiazolyl]-4-pyridinecarboxamide in the solvent of choice, which should be purged with an inert gas. N-[4-(2,4-Dimethylphenyl)-2-thiazolyl]-4-pyridinecarboxamide is soluble in DMSO.

### Description

N-[4-(2,4-Dimethylphenyl)-2-thiazolyl]-4-pyridinecarboxamide is a derivative of INH1 (Item No. 29015) with anticancer activity.<sup>1</sup> It is cytotoxic to MDA-MB-231, MDA-MB-488, HeLa, and K562 cells (IC<sub>50</sub>s = 1.1, 3.4, 1.6, and 1.2 μM, respectively).

### Reference

1. Qiu, X.-L., Li, G., Wu, G., *et al.* Synthesis and biological evaluation of a series of novel inhibitor of Nek2/Hec1 analogues. *J. Med. Chem.* **52**(6), 1757-1767 (2009).

#### 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, 12/19/2022

#### 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