

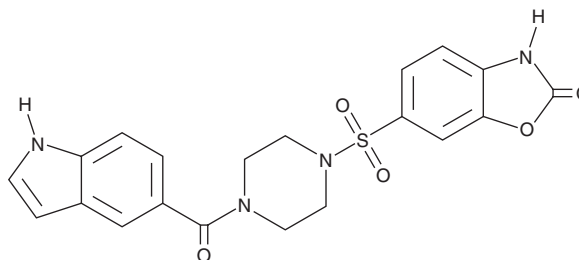
# PRODUCT INFORMATION



## TH-1760

Item No. 40469

**CAS Registry No.:** 2567914-01-4  
**Formal Name:** 6-[[4-(1H-indol-5-ylcarbonyl)-1-piperazinyl]sulfonyl]-2(3H)-benzoxazolone  
**MF:** C<sub>20</sub>H<sub>18</sub>N<sub>4</sub>O<sub>5</sub>S  
**FW:** 426.4  
**Purity:** ≥95%  
**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

TH-1760 is supplied as a solid. A stock solution may be made by dissolving the TH-1760 in the solvent of choice, which should be purged with an inert gas. TH-1760 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

### Description

TH-1760 is an inhibitor of nucleotide triphosphate diphosphatase NUDT15 (IC<sub>50</sub> = 25 nM).<sup>1</sup> It is selective for NUDT15 over a panel of additional Nudix family hydrolases and nucleotide pyrophosphatases at 100 μM, as well as a panel of 44 kinases at 10 μM. TH-1760 enhances thiopurine-induced cell death in NB4 and HL-60 cancer cells in a NUDT15-dependent manner.

### Reference

1. Zhang, S.M., Desroses, M., Hagenkort, A., *et al.* Development of a chemical probe against NUDT15. *Nat. Chem. Biol.* **16**(10), 1120-1128 (2020).

#### 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, 04/08/2024

#### 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](http://WWW.CAYMANCHEM.COM)