

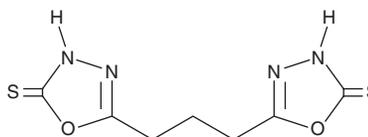
# PRODUCT INFORMATION



## CAY10761

Item No. 29865

**CAS Registry No.:** 333409-31-7  
**Formal Name:** 5,5'-(1,3-propanediyl)bis-1,3,4-oxadiazole-2(3H)-thione  
**MF:** C<sub>7</sub>H<sub>8</sub>N<sub>4</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 244.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 266 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

CAY10761 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10761 in the solvent of choice, which should be purged with an inert gas. CAY10761 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10761 in these solvents is approximately 30 mg/ml.

CAY10761 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10761 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10761 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

CAY10761 is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase 1 (ENPP1; IC<sub>50</sub>s = 467 and 429 μM for the human and snake venom enzymes, respectively).<sup>1,2</sup> It also inhibits mushroom tyrosinase (K<sub>i</sub> = 1.9 μM) and urease from jack bean, *P. mirabilis*, and *B. pasteurii* (IC<sub>50</sub>s = 0.093, <0.125, and 0.089 mM, respectively, at pH 8.2).<sup>3,4</sup>

### References

1. Khan, K.M., Fatima, N., Rasheed, M., *et al.* 1,3,4-Oxadiazole-2(3H)-thione and its analogues: A new class of non-competitive nucleotide pyrophosphatases/phosphodiesterases 1 inhibitors. *Bioorg. Med. Chem.* **17(22)**, 7816-7822 (2009).
2. Onyedibe, K.I., Wang, M., and Sintim, H.O. ENPP1, an old enzyme with new functions, and small molecule inhibitors - A STING in the tale of ENPP1. *Molecules* **24(22)**, E4192 (2019).
3. Ghani, U., and Ullah, N. New potent inhibitors of tyrosinase: Novel clues to binding of 1,3,4-thiadiazole-2(3H)-thiones, 1,3,4-oxadiazole-2(3H)-thiones, 4-amino-1,2,4-triazole-5(4H)-thiones, and substituted hydrazides to the dicopper active site. *Bioorg. Med. Chem.* **18(11)**, 4042-4048 (2010).
4. Amtul, Z., Rasheed, M., Choudhary, M.I., *et al.* Kinetics of novel competitive inhibitors of urease enzymes by a focused library of oxadiazoles/thiadiazoles and triazoles. *Biochem. Biophys. Res. Commun.* **319(3)**, 1053-1063 (2004).

#### 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, 03/21/2024

#### 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