

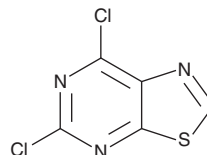
# PRODUCT INFORMATION



## 5,7-Dichlorothiazolo[5,4-d]pyrimidine

Item No. 36489

CAS Registry No.: 13479-88-4  
Formal Name: 5,7-dichloro-thiazolo[5,4-d]pyrimidine  
MF:  $C_5HCl_2N_3S$   
FW: 206.1  
Purity:  $\geq 95\%$   
UV/Vis.:  $\lambda_{max}$ : 222 nm  
Supplied as: A solid  
Storage:  $-20^\circ C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

5,7-Dichlorothiazolo[5,4-d]pyrimidine is supplied as a solid. A stock solution may be made by dissolving the 5,7-dichlorothiazolo[5,4-d]pyrimidine in the solvent of choice, which should be purged with an inert gas. 5,7-Dichlorothiazolo[5,4-d]pyrimidine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5,7-dichlorothiazolo[5,4-d]pyrimidine in these solvents is approximately 1 and 2 mg/ml, respectively.

### Description

5,7-Dichlorothiazolo[5,4-d]pyrimidine is a building block.<sup>1-4</sup> It has been used in the synthesis of a variety of compounds, including spleen tyrosine kinase (Syk), mammalian target of rapamycin (mTOR), and PI3K $\delta$  inhibitors, as well as HIV-1 non-nucleoside reverse transcriptase inhibitors (NNRTIs) with antiviral activity.

### References

1. Lucas, M.C., Goldstein, D.M., Hermann, J.C., *et al.* Rational design of highly selective spleen tyrosine kinase inhibitors. *J. Med. Chem.* **55**(23), 10414-10423 (2012).
2. Bonazzi, S., Goold, C.P., Gray, A., *et al.* Discovery of a brain-penetrant ATP-competitive inhibitor of the mechanistic target of rapamycin (mTOR) for CNS disorders. *J. Med. Chem.* **63**(3), 1068-1083 (2020).
3. Murray, J.M., Sweeney, Z.K., Chan, B.K., *et al.* Potent and highly selective benzimidazole inhibitors of PI3-kinase delta. *J. Med. Chem.* **55**(17), 7686-7695 (2012).
4. Kang, D., Zhao, T., Wang, Z., *et al.* Discovery of piperidine-substituted thiazolo[5,4-d]pyrimidine derivatives as potent and orally bioavailable HIV-1 non-nucleoside reverse transcriptase inhibitors. *Commun. Chem.* **2**, 74 (2019).

#### 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