

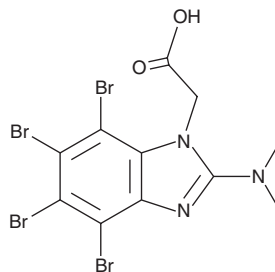
PRODUCT INFORMATION



TMCB

Item No. 27652

CAS Registry No.: 1085822-09-8
Formal Name: 4,5,6,7-tetrabromo-2-(dimethylamino)-1H-benzimidazole-1-acetic acid
MF: C₁₁H₉Br₄N₃O₂
FW: 534.8
Purity: ≥98%
UV/Vis.: λ_{max}: 240, 274 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

TMCB is supplied as a crystalline solid. A stock solution may be made by dissolving the TMCB in the solvent of choice, which should be purged with an inert gas. TMCB is soluble in the organic solvent DMSO at a concentration of approximately 25 mM.

Description

TMCB is a dual inhibitor of casein kinase 2 (CK2) and ERK8 (IC₅₀ = 0.5 μM for both).¹ It also binds to Pim-1, HIPK2, and DYRK1A (K_s = 8.65, 15.25, and 11.9 μM, respectively) and inhibits greater than 50% of the activity of five kinases, including CK2, ERK8, and DYRK2, in a panel of 78 kinases at 10 μM. TMCB induces apoptosis in Jurkat cells when used at concentrations ranging from 10 to 50 μM.

References

1. Pagano, M.A., Bain, J., Kazimierczuk, Z., *et al.* The selectivity of inhibitors of protein kinase CK2: An update. *Biochem. J.* **415**(3), 353-365 (2008).

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, 12/08/2022

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