

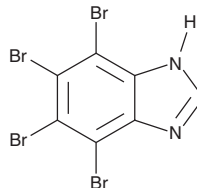
PRODUCT INFORMATION



4,5,6,7-Tetrabromobenzimidazole

Item No. 18886

CAS Registry No.: 577779-57-8
Formal Name: 4,5,6,7-tetrabromo-1H-benzimidazole
Synonyms: Casein Kinase II Inhibitor XII, TBBz
MF: C₇H₂Br₄N₂
FW: 433.7
Purity: ≥98%
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

4,5,6,7-Tetrabromobenzimidazole (TBBz) is supplied as a crystalline solid. A stock solution may be made by dissolving the TBBz in the solvent of choice, which should be purged with an inert gas. TBBz is soluble in organic solvents such as ethanol and DMSO.

Description

Casein kinase 2 (CK2) is a constitutive protein kinase involved in many signal transduction pathways. CK2 is known to negatively regulate apoptosis, and its activity is increased in many proliferating tissues and tumors. TBBz is a selective, ATP-competitive inhibitor of CK2 with K_i values ranging from 0.5-1 μ M in rat liver, *S. cerevisiae*, *C. tropicalis*, and *N. crassa*.¹ It is inactive against PKA and PKC and a relatively weak inhibitor of CK1 ($K_i = 54.4$ -64.2 μ M).¹

Reference

1. Zien, P., Bretner, M., Zastapilo, K., *et al.* Selectivity of 4,5,6,7-tetrabromobenzimidazole as an ATP-competitive potent inhibitor of protein kinase CK2 from various sources. *Biochem. Biophys. Res. Commun.* **306**(1), 129-133 (2003).

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, 11/04/2022

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