

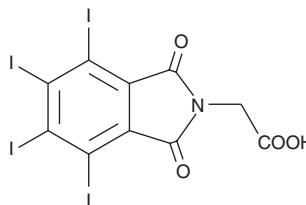
PRODUCT INFORMATION



CAY10578

Item No. 10011264

CAS Registry No.: 19231-60-8
Formal Name: 4,5,6,7-tetraiodo-1,3-dioxo-2-isoindolineacetic acid
MF: C₁₀H₃I₄NO₄
FW: 708.8
Purity: ≥95%
UV/Vis.: λ_{max}: 263 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

CAY10578 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10578 in the solvent of choice, which should be purged with an inert gas. CAY10578 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CAY10578 in these solvents is approximately 5 and 10 mg/ml, respectively.

CAY10578 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10578 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10578 has a solubility of approximately 0.15 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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, which lends promise as an anticancer drug target. CAY10578 is a potent inhibitor of CK2 with an IC₅₀ value of 0.3 μM.¹ This compound is ATP competitive, displaying a K_i value of 0.2 μM and only minimally inhibits the activities of other protein kinases (e.g., GSK3, Cdk5, and MSK1).¹

Reference

1. Golub, A.G., Yakovenko, O.Y., Prykhod'ko, A.O., *et al.* Evaluation of 4,5,6,7-tetrahalogeno-1*H*-isoindole-1,3(2*H*)-diones as inhibitors of human protein kinase CK2. *Biochem. Biophys. Acta* **1784**(1), 143-149 (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, 11/14/2022

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