

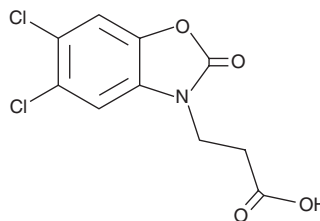
PRODUCT INFORMATION



GSK180

Item No. 19217

CAS Registry No.: 1799725-26-0
Formal Name: 5,6-dichloro-2-oxo-3(2H)-benzoxazolepropanoic acid
MF: C₁₀H₇Cl₂NO₄
FW: 276.1
Purity: ≥95%
Supplied as: A 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

GSK180 is supplied as a solid. A stock solution may be made by dissolving the GSK180 in the solvent of choice, which should be purged with an inert gas. GSK180 is soluble (≥10 mg/ml) in DMSO.

Description

GSK180 is an inhibitor of kynurenine 3-monooxygenase (IC₅₀ = 6.3 nM).¹ It is selective for human over *P. fluorescens* kynurenine 3-monooxygenase (IC₅₀ = 500 nM). GSK180 (27 mg/kg) increases circulating levels of kynurenine and kynurenic acid in mice.² It protects against secondary acute lung injury and renal injury in a rat model of experimentally induced acute pancreatitis multiple organ dysfunction syndrome (AP-MODS).

References

1. Liddle, J., Beaufils, B., Binnie, M., *et al.* The discovery of potent and selective kynurenine 3-monooxygenase inhibitors for the treatment of acute pancreatitis. *Bioorg. Med. Chem. Lett.* **27(9)**, 2023-2028 (2017).
2. Mole, D.J., Webster, S.P., Uings, I., *et al.* Kynurenine-3-monooxygenase inhibition prevents multiple organ failure in rodent models of acute pancreatitis. *Nat. Med.* **22(2)**, 202-209 (2016).

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, 10/15/2024

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