

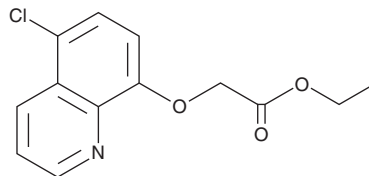
PRODUCT INFORMATION



A2793

Item No. 41081

CAS Registry No.: 88349-90-0
Formal Name: 2-[(5-chloro-8-quinolinyl)oxy]-acetic acid, ethyl ester
MF: C₁₃H₁₂ClNO₃
FW: 265.7
Purity: ≥98%
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

A2793 is supplied as a solid. A stock solution may be made by dissolving the A2793 in the solvent of choice, which should be purged with an inert gas. A2793 is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

A2793 is an inhibitor of the two-pore domain potassium channel K_{2P}18.1/TRESK (IC₅₀ = 6.8 μM in ionomycin-stimulated *Xenopus* oocytes expressing the mouse receptor).¹ It is selective for K_{2P}18.1/TRESK over K_{2P}2.1/TREK1, K_{2P}10.1/TREK2, K_{2P}4.1/TRAAK, K_{2P}3.2/TASK2, K_{2P}3.3/TASK3, K_{2P}16.1/TALK1, and K_{2P}13.1/THIK1 at 100 μM but does inhibit K_{2P}3.1/TASK1.

Reference

1. Lengyel, M., Erdélyi, F., Pergel, E., *et al.* Chemically modified derivatives of the activator compound cloxyquin exert inhibitory effect on TRESK (K_{2P}18.1) background potassium channel. *Mol. Pharmacol.* **95**(6), 652-660 (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, 05/31/2024

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