

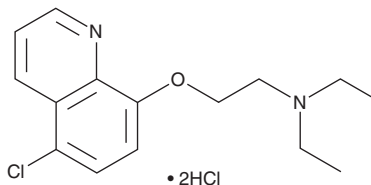
PRODUCT INFORMATION



A2764 (hydrochloride)

Item No. 36943

CAS Registry No.: 861038-72-4
Formal Name: 2-[(5-chloro-8-quinolinyl)oxy]-N,N-diethyl-ethanamine, dihydrochloride
MF: C₁₅H₁₉ClN₂O • 2HCl
FW: 351.7
Purity: ≥98%
UV/Vis.: λ_{max}: 244 nm
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

A2764 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the A2764 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. A2764 (hydrochloride) is slightly soluble in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of A2764 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of A2764 (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

A2764 is an inhibitor of the two-pore domain potassium channel K_{2p}18.1/TRESK (IC₅₀ = 11.8 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}3.1/TASK1, K_{2p}3.2/TASK2, K_{2p}3.3/TASK3, K_{2p}16.1/TALK1, and K_{2p}13.1/THIK1 at 100 μM. A2764 (100 μM) inhibits basal potassium currents in isolated mouse dorsal root ganglion neurons.

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, 12/21/2022

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