

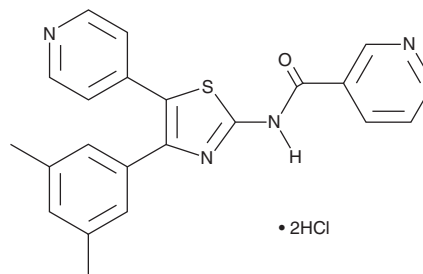
PRODUCT INFORMATION



DPTN (hydrochloride)

Item No. 41540

CAS Registry No.: 325767-87-1
Formal Name: N-[4-(3,5-dimethylphenyl)-5-(4-pyridinyl)-2-thiazolyl]-3-pyridinecarboxamide, dihydrochloride
MF: $C_{22}H_{18}N_4OS \cdot 2HCl$
FW: 459.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Description

DPTN is an adenosine A_3 receptor antagonist ($K_i = 0.36$ nM for the human receptor).¹ It is selective for the adenosine A_3 receptor over the adenosine A_1 and A_{2A} receptors ($K_{is} = >666$ and >826 nM, respectively). DPTN (10 mg/kg) prevents dye leakage in a rat model of plasma extravasation induced by the adenosine A_3 receptor agonist IB-MECA (piclidenoson; Item No. 28445). It, alone and in combination with the synthetic glucocorticoid dexamethasone (Item No. 11015), inhibits airway hypersensitivity to acetylcholine in a rat model of allergen-induced asthma stimulated by ovalbumin inhalation when administered at a dose of 10 mg/kg.

Reference

1. Miwatashi, S., Arikawa, Y., Matsumoto, T., *et al.* Synthesis and biological activities of 4-phenyl-5-pyridyl-1,3-thiazole derivatives as selective adenosine A_3 antagonists. *Chem. Pharm. Bull.* **56**(8), 1126-1137 (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, 08/23/2024

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