

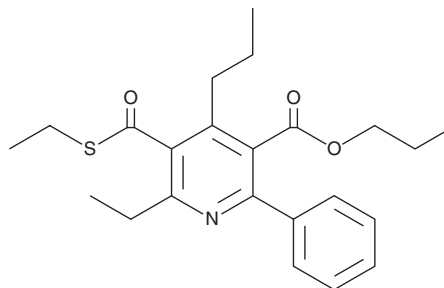
PRODUCT INFORMATION



MRS1523

Item No. 17128

CAS Registry No.: 212329-37-8
Formal Name: 6-ethyl-5-[(ethylthio)carbonyl]-2-phenyl-4-propyl-3-pyridinecarboxylic acid propyl ester
MF: C₂₃H₂₉NO₃S
FW: 399.5
Purity: ≥98%
UV/Vis.: λ_{max}: 247 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

MRS1523 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of MRS1523 in these solvents is approximately 50 mg/ml.

Description

MRS1523 is an adenosine A₃ receptor antagonist.¹ It binds to the human adenosine A₃ receptor (K_i = 0.0189 μM), as well as selectively binds rat adenosine A₃ over A₁ and A_{2A} receptors (K_is = 0.113, 15.6, and 2.05 μM, respectively) in radioligand binding assays. It inhibits the proliferation of B16-BL6 mouse melanoma cells when used at a concentration of 1 μM.²

Reference

1. Li, A.H., Moro, S., Melman, N., *et al.* Structure-activity relationships and molecular modeling of 3, 5-diacyl-2,4-dialkylpyridine derivatives as selective A₃ adenosine receptor antagonists. *J. Med. Chem.* **41(17)**, 3186-3201 (1998).
2. Yoshiyama, N., Yamada, S., Takeuchi, C., *et al.* Cordycepin (3'-deoxyadenosine) inhibits the growth of B16-BL6 mouse melanoma cells through the stimulation of adenosine A₃ receptor followed by glycogen synthase kinase-3β activation and cyclin D₁ suppression. *Naunyn Schmiedebergs Arch Pharmacol.* **377(4-6)**, 591-595 (2007).

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, 08/24/2022

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