

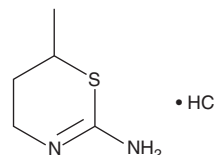
PRODUCT INFORMATION



AMT (hydrochloride)

Item No. 81010

CAS Registry No.: 21463-31-0
Formal Name: 5,6-dihydro-6-methyl-4H-1,3-thiazin-2-amine, hydrochloride
MF: $C_5H_{10}N_2S \cdot HCl$
FW: 166.7
Purity: $\geq 98\%$
Supplied as: A crystalline 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

AMT (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the AMT (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AMT (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). It is also soluble in water. The solubility of AMT (hydrochloride) in ethanol and water is approximately 25 mg/ml and approximately 16 and 8 mg/ml in DMSO and DMF, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

AMT is a potent, relatively selective inhibitor of iNOS with a K_i of 4.2 nM in mouse macrophages. It inhibits other isoforms of NOS, but at higher concentrations; the IC_{50} values for rat nNOS and bovine eNOS are 34 and 150 nM, respectively.¹

Reference

1. Nakane, M., Klinghofer, V., Kuk, J.E., *et al.* Novel potent and selective inhibitors of inducible nitric oxide synthase. *Mol. Pharmacol.* **47**(4), 831-834 (1995).

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, 04/04/2023

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