

PRODUCT INFORMATION



Fenethylamine (hydrochloride)

Item No. 14146

CAS Registry No.: 1892-80-4
Formal Name: 3,7-dihydro-1,3-dimethyl-7-[2-[(1-methyl-2-phenylethyl)amino]ethyl]-1H-purine-2,6-dione, monohydrochloride

Synonyms: Amphetaminoethyltheophylline, Amfetyline, NSC 169872

MF: C₁₈H₂₃N₅O₂ • HCl
FW: 377.9

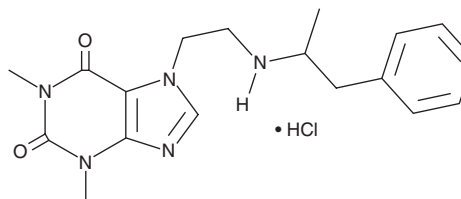
Purity: ≥98%

UV/Vis.: λ_{max}: 275 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

Fenethylamine (hydrochloride) (Item No. 14146) is an analytical reference standard categorized as a stimulant.^{1,2} Fenethylamine is a theophylline derivative of amphetamine. It has been found in tablets seized by authorities.¹ Formulations containing fenethylamine have been used in the treatment of attention-deficit/hyperactivity disorder (ADHD) and narcolepsy. Fenethylamine is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. Kristen, G., Schaefer, A., and von Schlichtegroll, A. Fenethylamine: Therapeutic use, misuse and/or abuse. *Drug Alcohol Depend.* **17(2-3)**, 259-271 (1986).
2. Katselou, M., Papoutsis, I., Nikolaou, P., et al. Fenethylamine (captagon) abuse - local problems from an old drug become universal. *Basic Clin. Pharmacol. Toxicol.* **119(2)**, 133-140 (2016).

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, 03/11/2024

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