

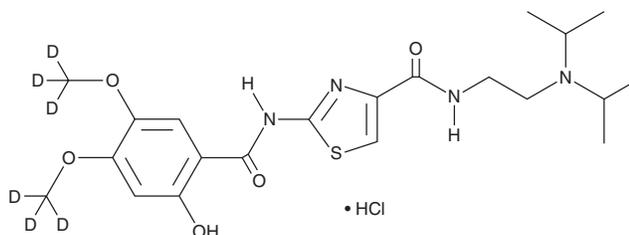
PRODUCT INFORMATION



Acotiamide-d₆ (hydrochloride)

Item No. 43489

Formal Name: N-(2-(diisopropylamino)ethyl)-2-(2-hydroxy-4,5-bis(methoxy-d₃)benzamido)thiazole-4-carboxamide, monohydrochloride
MF: C₂₁H₂₄D₆N₄O₅S • HCl
FW: 493.0
Chemical Purity: ≥95% (Acotiamide)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
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

Acotiamide-d₆ is intended for use as an internal standard for the quantification of acotiamide by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Acotiamide-d₆ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the acotiamide-d₆ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Acotiamide-d₆ (hydrochloride) is soluble (≥10 mg/ml) in methanol and DMSO.

Description

Acotiamide is an inhibitor of acetylcholinesterase (AChE; IC₅₀ = 0.384 μM). It potentiates ACh-induced contraction of isolated guinea pig gastric and antrum strips when used at a concentration of 3 μM.¹ Formulations containing acotiamide have been used in the treatment of functional dyspepsia.

Reference

1. Ito, K., Kawachi, M., Matsunaga, Y., *et al.* Acotiamide hydrochloride, a therapeutic agent for functional dyspepsia, enhances acetylcholine-induced contraction via inhibition of acetylcholinesterase activity in circular muscle strips of guinea pig stomach. *Drug Res. (Stuttg.)* **66(4)**, 196-202 (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, 05/08/2025

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