

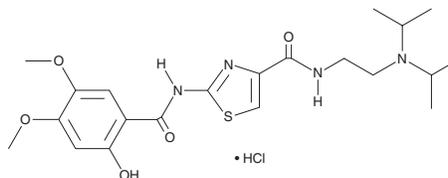
PRODUCT INFORMATION



Acotiamide (hydrochloride)

Item No. 43731

CAS Registry No.: 185104-11-4
Formal Name: N-[2-[bis(1-methylethyl)amino]ethyl]-2-[(2-hydroxy-4,5-dimethoxybenzoyl)amino]-4-thiazolecarboxamide, monohydrochloride
Synonym: YM-443
MF: C₂₁H₃₀N₄O₅S • HCl
FW: 487.0
Purity: ≥98%
Supplied as: A crystalline 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 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the acotiamide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Acotiamide (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. Acotiamide (hydrochloride) is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in 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, 07/14/2025

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