

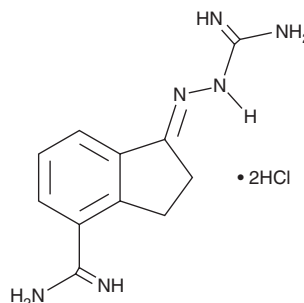
PRODUCT INFORMATION



Sardomozide (hydrochloride)

Item No. 34901

CAS Registry No.: 138794-73-7
Formal Name: 2-[4-(aminoiminomethyl)-2,3-dihydro-1H-inden-1-ylidene]-hydrazinecarboximidamide, dihydrochloride
Synonyms: CGP 48664, SAM486A
MF: C₁₁H₁₄N₆ • 2HCl
FW: 303.2
Purity: ≥95%
UV/Vis.: λ_{max}: 232, 272, 281, 309 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Sardomozide (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the sardomozide (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Sardomozide (hydrochloride) is soluble in ethanol.

Description

Sardomozide is an inhibitor of S-adenosylmethionine decarboxylase (SAMDC), an enzyme involved in polyamine biosynthesis (IC₅₀ = 0.005 μM for the rat liver enzyme).¹ It is selective for SAMDC over diamine oxidase (DAO; IC₅₀ = 18 μM for the rat small intestine enzyme). Sardomozide inhibits the proliferation of T24 bladder cancer cells (IC₅₀ = 0.71 μM). It decreases intracellular levels of spermidine and spermine and increases intracellular putrescine levels in L1210 murine leukemia cells when used at a concentration of 3 μM.² Sardomozide (0.2 and 0.4 μM) inhibits HIV-1 replication in PM1 cells.³ It reduces tumor growth in a SK-MEL-24 melanoma mouse xenograft model when administered at doses of 0.5 and 5 mg/kg.²

References

1. Stanek, J., Caravatti, G., Frei, J., et al. 4-Amidinoindan-1-one 2'-amidinohydrazone: A new potent and selective inhibitor of S-Adenosylmethionine decarboxylase. *J. Med. Chem.* **36(15)**, 2168-2171 (1993).
2. Regenass, U., Mett, H., Stanek, J., et al. CGP 48664, a new S-adenosylmethionine decarboxylase inhibitor with broad spectrum antiproliferative and antitumor activity. *Cancer Res.* **54(12)**, 3210-3217 (1994).
3. Schäfer, B., Hauber, I., Bunk, A., et al. Inhibition of multidrug-resistant HIV-1 by interference with cellular S-adenosylmethionine decarboxylase activity. *J. Infect. Dis.* **194(6)**, 740-750 (2006).

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, 09/10/2021

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