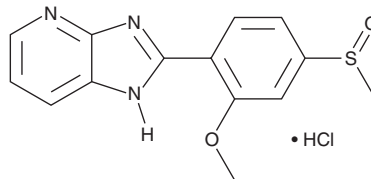


PRODUCT INFORMATION

Sulmazole (hydrochloride)

Item No. 45128

CAS Registry No.: 53930-42-0
Formal Name: 2-[2-methoxy-4-(methylsulfinyl)phenyl]-3H-imidazo[4,5-b]pyridine, monohydrochloride
Synonyms: AR-L 115BS, Sumazol
MF: C₁₄H₁₃N₃O₂S • HCl
FW: 323.8
Purity: ≥98%
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

Sulmazole is supplied as a solid. A stock solution may be made by dissolving the sulmazole in the solvent of choice, which should be purged with an inert gas. Sulmazole is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of sulmazole can be prepared by directly dissolving the solid in aqueous buffers. Sulmazole is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Sulmazole is a cardiotonic agent.^{1,2} It increases cAMP accumulation *via* antagonism of the adenosine A₁ receptor (IC₅₀ = 78.1 μM), inhibition of phosphodiesterase (PDE), and reduction of G_i-mediated inhibition of adenylate cyclase in rat adipocyte membranes.¹ Sulmazole also selectively inhibits PDE5 over PDE3 (IC₅₀s = 21 and 150 μM, respectively) in a cell-free assay.² It restores cardiac output, the force of right ventricular contraction, and heart rate, and decreases right atrial pressure in a dog heart-lung preparation model of pentobarbital-induced depressed cardiac function.³

References

1. Parsons, W.J., Ramkumar, V., and Stiles, G.L. The new cardiotonic agent sulmazole is an A₁ adenosine receptor antagonist and functionally blocks the inhibitory regulator, G_i. *Mol. Pharmacol.* **33**(4), 441-448 (1988).
2. Coates, W.J., Connolly, B., Dhanak, D., *et al.* Cyclic nucleotide phosphodiesterase inhibition by imidazopyridines: Analogues of sulmazole and isomazole as inhibitors of the cGMP specific phosphodiesterase. *J. Med. Chem.* **36**(10), 1387-1392 (1993).
3. Morita, T., Satoh, K., and Taira, N. Improvement of cardiac performance by AR-L 115 BS, a new cardiotonic drug, in the experimental failing dog heart. *Jpn. Heart J.* **25**(3), 439-445 (1984).

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, 03/11/2026

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