

PRODUCT INFORMATION



AL 34662 (formate)

Item No. 40489

Formal Name: 1-[(2S)-2-aminopropyl]-1H-indazol-6-ol, monoformic acid

MF: C₁₀H₁₃N₃O • CH₂O₂

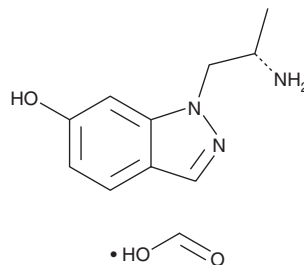
FW: 237.3

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

AL 34662 (formate) is supplied as a solid. A stock solution may be made by dissolving the AL 34662 (formate) in the solvent of choice, which should be purged with an inert gas. AL 34662 (formate) is soluble in the organic solvent DMSO.

Description

AL 34662 is an agonist of the serotonin (5-HT) receptor subtype 5-HT₂.¹ It selectively binds to 5-HT₂ receptors (IC₅₀ = 1.5 and 0.77 nM in isolated and washed human or rat cerebral cortex homogenates, respectively) over a panel of 19 neurotransmitter and steroid receptors at 10 μM. AL 34662 induces calcium mobilization in a reporter assay using primary human ciliary muscle cells (EC₅₀ = 140 nM). It reduces intraocular pressure (IOP) in a cynomolgus monkey model of trabeculoplasty-induced ocular hypertension when administered at doses of 100 or 300 μg/eye.

Reference

1. Sharif, N.A., McLaughlin, M.A., and Kelly, C.R. AL-34662: A potent, selective, and efficacious ocular hypotensive serotonin-2 receptor agonist. *J. Ocul. Pharmacol. Ther.* **23**(1), 1-13 (2007).

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, 02/13/2025

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