

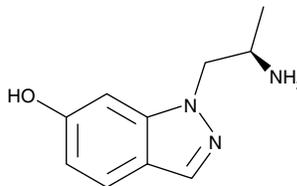
Product Information



AL 34662

Item No. 10011546

CAS Registry No.: 210580-75-9
Formal Name: 1-[(2S)-2-aminopropyl]-1H-indazol-6-ol
Synonym: AL 34497
MF: C₁₀H₁₃N₃O
FW: 191.2
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 214, 289 nm



Laboratory Procedures

For long term storage, we suggest that AL 34662 be stored as supplied at -20°C. It should be stable for at least two years.

AL 34662 is supplied as a crystalline solid. A stock solution may be made by dissolving the AL 34662 in an organic solvent purged with an inert gas. AL 34662 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AL 34662 in ethanol is approximately 1 mg/ml and approximately 10 mg/ml in DMSO and DMF.

AL 34662 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AL 34662 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AL 34662 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The transduction of neurobehavioral effects by serotonin (5-hydroxy tryptamine; 5-HT) is mediated *via* at least six major 5-HT receptor subtypes, including 5-HT₂. AL 34662 is a potent 5-HT₂ receptor agonist with ocular hypotensive activity. It binds to the human and rat 5-HT₂ receptors in cerebral cortex homogenates with IC₅₀ values of 1.5 and 0.77 nM, respectively.¹ AL 34662 binds to the recombinant human 5-HT_{2A}, 5-HT_{2B}, and 5-HT_{2C} receptors with IC₅₀ values of 14.5, 8.1, and 3 nM, respectively. In ocular hypertensive cynomolgus monkey eyes, AL 34662 lowered intraocular pressure (IOP) 25% at a dose of 100 µg and 33% at 300 µg (six hours post dose) with minimal side effects.¹

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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011546

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog**.

Copyright Cayman Chemical Company, 06/24/2014

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com