

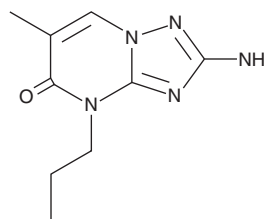
PRODUCT INFORMATION



ICI 63197

Item No. 35538

CAS Registry No.: 27277-00-5
Formal Name: 2-amino-6-methyl-4-propyl-[1,2,4]triazolo[1,5-a]pyrimidin-5(4H)-one
MF: C₉H₁₃N₅O
FW: 207.2
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 307 nm
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

ICI 63197 is supplied as a solid. A stock solution may be made by dissolving the ICI 63197 in the solvent of choice, which should be purged with an inert gas. ICI 63197 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ICI 63197 in these solvents is approximately 10 mg/ml. It is also slightly soluble 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 ICI 63197 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of ICI 63197 in PBS (pH 7.2) is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ICI 63197 is an inhibitor of phosphodiesterase 3 (PDE3) and PDE4 (K_is = 9 and 10 μM, respectively).¹ It is selective for PDE3 and PDE4 over PDE1 and PDE2 (K_is = 487 and >1,000 μM, respectively). ICI 63197 inhibits acetylcholine-induced bronchospasm in isolated guinea pig lungs and histamine-induced bronchospasm in conscious guinea pigs.² It decreases motor activity and rectal temperature and induces forepaw shaking, head twitches, and grooming in rats in a dose-dependent manner.³

References

1. Hoey, M. and Houslay, M.D. Identification and selective inhibition of four distinct soluble forms of cyclic nucleotide phosphodiesterase activity from kidney. *Biochem. Pharmacol.* **40**(2), 193-202 (1990).
2. Davies, G.E. Antibronchoconstrictor activity of two new phosphodiesterase inhibitors, a triazolopyrazine (ICI 58 301) and a triazolopyrimidine (ICI 63 197). *J. Pharm. Pharmacol.* **25**(9), 681-689 (1973).
3. Wachtel, H. Characteristic behavioural alterations in rats induced by rolipram and other selective adenosine cyclic 3',5'-monophosphate phosphodiesterase inhibitors. *Psychopharmacology (Berl.)* **77**(4), 309-316 (1982).

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, 12/12/2022

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