

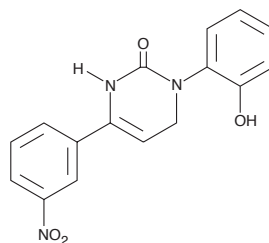
Product Information



Icilin

Catalog No. 10137

CAS Registry No.: 36945-98-9
Formal Name: 3,6-dihydro-1-(2-hydroxyphenyl)-4-(3-nitrophenyl)-2(1H)-pyrimidinone
Synonym: AG 3-5
MF: C₁₆H₁₃N₃O₄
FW: 311.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

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

Icilin is supplied as a crystalline solid. A stock solution may be made by dissolving the icilin in an organic solvent purged with an inert gas. Icilin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of icilin in these solvents is approximately 25 mg/ml.

Icilin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, icilin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Icilin has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. Therefore, further dilutions of the organic solvent 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. We do not recommend storing the aqueous solution for more than one day.

Nociceptive sensory peripheral neurons often express both heat sensitive (VR₁) and cold-sensitive (CMR1 (rat); TRPM8 (human)) receptors that are part of the transient receptor potential (TRP) superfamily. Icilin (AG 3-5) is a synthetic CMR1/TRPM8 super agonist that is 2.5-fold more efficacious and nearly 200-fold more potent than the reference cold thermosensory agonist *l*-menthol.¹ Icilin induces sensations of intense cold when applied orally in humans, and induces 'wet dog shakes', a behavioral marker of cold sensation, when given to rats. Icilin should serve as the reference cold nociceptive agonist for TRP-type ion channels in the future.²

References

1. McKemy, D.D., Neuhauser, W.M., and Julius, D. Identification of a cold receptor reveals a general role for TRP channels in thermosensation. *Nature* **416**, 52-58 (2002).
2. Wei, E.T. and Seid, D.A. AG-3-5: a chemical producing sensations of cold. *J. Pharm. Pharmacol.* **35**, 110-112 (1983).

Related Products

Capsaicin - Cat. No. 92350 • Dihydrocapsaicin - Cat. No. 92355

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

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL 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 Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty of any kind, expressed or implied, including, but not limited to, the warranties of fitness for a particular purpose and merchantability, which extends beyond the description of the chemicals on the face hereof, except that the material will meet our specifications at the time of delivery. Buyer's exclusive remedy and Cayman Chemical Company's sole liability hereunder shall be limited to refund of the purchase price of, or at Cayman Chemical Company's option the replacement of, all material that does not meet our specifications. Cayman Chemical Company shall not be liable otherwise or for incidental or consequential damages, including, but not limited to, the costs of handling. Said refund or replacement is conditioned on Buyer giving written notice to Cayman Chemical Company within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within said thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material. Copyright Cayman Chemical Company, 03/20/2006