

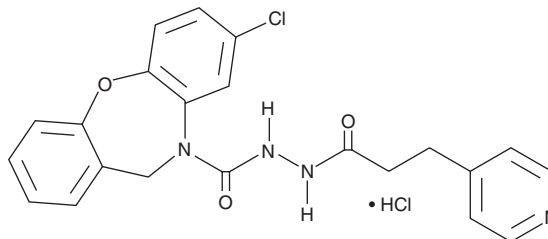
Product Information



SC-51089

Catalog No. 10011561

CAS Registry No.: 146033-02-5
Formal Name: 8-chloro-2-[1-oxo-3-(4-pyridinyl)propyl]hydrazide-dibenz[b,f][1,4]oxazepine-10(11H)-carboxylic acid, monohydrochloride
Synonym: CID132748
MF: C₂₂H₁₉ClN₄O₃ • HCl
FW: 459.3
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that SC-51089 be stored as supplied at -20°C. It should be stable for at least two years.

SC-51089 is supplied as a crystalline solid. A stock solution may be made by dissolving the SC-51089 in the solvent of choice. SC-51089 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SC-51089 in these solvents is approximately 20 mg/ml.

SC-51089 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SC-51089 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SC-51089 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The prostaglandin E₂ (PGE₂) receptor EP₁ is involved in triggering PGE₂-mediated pain as well as neuronal survival and growth. SC-51089 is a selective EP₁ antagonist, first shown to have *in vivo* analgesic activity in mice (ED₅₀ = 6.3 mg/kg when given subcutaneously) and in the rat.^{1,2} SC-51089 also inhibits the growth of glioma cell lines *in vitro* (IC₅₀ = -1 μM) and slows tumor growth *in vivo*.³ In addition, SC-51089 is neuroprotective, attenuating neuronal cell death in response to oxidative stress, an effect that is also found in EP₁^{-/-} mice and mediated by PGE₂.^{4,5}

References

- Hallinan, E.A., Hagen, T.J., Husa, R.K., *et al.* N-substituted dibenzoxazepines as analgesic PGE₂ antagonists. *J. Med. Chem.* **36**, 3293-3299 (1993).
- Malmberg, A.B., Rafferty, M.F., and Yaksh, T.L. Antinociceptive effect of spinally delivered prostaglandin E receptor antagonists in the formalin test on the rat. *Neurosci. Lett.* **173**, 193-196 (1994).
- Matsuo, M., Yoshida, N., Zaitso, M., *et al.* Inhibition of human glioma cell growth by a PHS-2 inhibitor, NS398, and a prostaglandin E receptor subtype EP₁-selective antagonist, SC51089. *Journal of Neuro-Oncology* **66**, 285-292 (2004).
- Carrasco, E., Casper, D., and Werner, P. PGE₂ receptor EP₁ renders dopaminergic neurons selectively vulnerable to low-level oxidative stress and direct PGE₂ neurotoxicity. *J. Neurosci. Res.* **85**, 3109-3117 (2007).
- Saleem, S., Li, R., Wei, G., *et al.* Effects of EP₁ receptor on cerebral blood flow in the middle cerebral artery occlusion model of stroke in mice. *J Neurosci Res.* **85(11)**, 2433-2440 (2007).

Related Products

AH 6809 - Cat. No. 14050 • SC-19220 - Cat. No. 14060 • GW 848687X - Cat. No. 10010410

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

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 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, 02/09/2010