

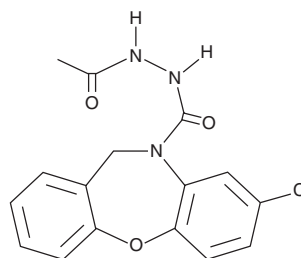
# PRODUCT INFORMATION



## SC-19220

Item No. 14060

**CAS Registry No.:** 19395-87-0  
**Formal Name:** 8-chloro-dibenz[b,f][1,4]oxazepine-10(11H)-carboxy-(2-acetyl)hydrazide  
**MF:** C<sub>16</sub>H<sub>14</sub>ClN<sub>3</sub>O<sub>3</sub>  
**FW:** 331.8  
**Purity:** ≥96%  
**UV/Vis.:** λ<sub>max</sub>: 203 nm  
**Supplied as:** A crystalline 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

SC-19220 is supplied as a crystalline solid. A stock solution may be made by dissolving the SC-19220 in an organic solvent, which should be purged with an inert gas. SC-19220 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SC-19220 in these solvents is approximately 14 mg/ml.

SC-19220 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SC-19220 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SC-19220 has a solubility of approximately 300 µg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

The prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) receptor EP<sub>1</sub> is involved in triggering PGE<sub>2</sub>-mediated pain, neuronal survival and growth as well as the suppression of tumor metastasis. SC-19220 is a dibenzoxazepine that acts as a selective antagonist of PGE<sub>2</sub> at the EP<sub>1</sub> receptor.<sup>1,2</sup> At doses between 0.3-300 µM, SC-19220 acts as a competitive antagonist of PGE<sub>2</sub>-induced smooth muscle contractions of guinea pig ileum and stomach.<sup>3,4</sup> SC-19220 also acts as a PGE<sub>2</sub> antagonist in the EP<sub>1</sub> receptor-mediated contraction of guinea pig trachea.<sup>5</sup> SC-19220 displaces radiolabeled PGE<sub>2</sub> from the cloned human EP<sub>1</sub> receptor with an IC<sub>50</sub> of 6.7 µM and exhibits no binding at the human EP<sub>2</sub> receptor.<sup>6,7</sup> It binds with apparent low affinity to the cloned mouse EP<sub>1</sub> receptor expressed in CHO cells.<sup>8</sup> In a model of bone resorption, 3-150 µM SC-19220 inhibited both 1,25 dihydroxy vitamin D<sub>3</sub>- and PGE<sub>2</sub>-stimulated osteoclastogenesis in mouse osteoclast precursors.<sup>9</sup> Antagonism of EP<sub>1</sub> with 1 µM SC-19220 has been shown to promote metastasis in a mouse model of metastatic breast cancer.<sup>10</sup> In a rodent model of Parkinson's disease, 1.5 µM SC-19220 prevented PGE<sub>2</sub>-mediated loss of dopaminergic neurons from rat substantia nigra.<sup>11</sup>

### References

1. Coleman, R.A., Eglen, R.M., Jones, R.L., et al. *IUPHAR Compendium* 1-12 (1997).
2. Zeng, L., An, S., and Goetzl, E.J. *Biochemistry* **35**, 7159-7164 (1996).
3. Rakovska, A. and Milenov, K. *Arch. Int. Pharmacodyn.* **268**, 59-69 (1984).
4. Sanner, J. *Ann. N. Y. Acad. Sci.* **180**, 396-409 (1971).
5. Coleman, R.A. and Kennedy, I. *Prostaglandins* **29**, 363-375 (1985).
6. Funk, C.D., Furci, L., Fitzgerald, G.A., et al. *J. Biol. Chem.* **268**, 26767-26772 (1993).
7. Bastien, L., Sawyer, N., Grygorczyk, R., et al. *J. Biol. Chem.* **269**, 11873-11877 (1994).
8. Kiriya, M., Ushikubi, F., Kobayashi, T., et al. *Br. J. Pharmacol.* **122**, 217-224 (1997).
9. Inoue, H., Tsujisawa, T., Fukuizumi, T., et al. *J. Endocrinol.* **161**(2), 231-236 (1999).
10. Ma, X., Kundu, N., Ioffe, O.B., et al. *Mol. Cancer Res.* **8**(10), 1310-1318 (2010).
11. Carrasco, E., Casper, D., and Werner, P. J. *Neurosci. Res.* **85**, 3109-3117 (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, 11/21/2022

#### 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