

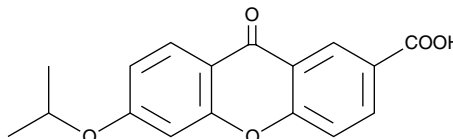
# Product Information



## AH 6809

Item No. 14050

**CAS Registry No.:** 33458-93-4  
**Formal Name:** 6-isopropoxy-9-oxoxanthene-2-carboxylic acid  
**MF:** C<sub>17</sub>H<sub>14</sub>O<sub>5</sub>  
**FW:** 298.3  
**Purity:** ≥98%  
**Stability:** ≥2 years at 4°C  
**Supplied as:** A crystalline solid  
**UV/Vis.:** λ<sub>max</sub>: 244, 305 nm



### Laboratory Procedures

For long term storage, we suggest that AH 6809 be stored as supplied at 4°C. It should be stable for at least two years.

AH 6809 is supplied as a crystalline solid. A stock solution may be made by dissolving the AH 6809 in an organic solvent. AH 6809 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AH 6809 in these solvents is approximately 0.5, 1, and 8 mg/ml, respectively.

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 AH 6809 can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of AH 6809 in PBS (pH 7.2) is approximately 0.34 mg/ml. For greater aqueous solubility, AH 6809 can be directly dissolved in 0.1 M Na<sub>2</sub>CO<sub>3</sub> (4 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

AH 6809 is an EP and DP receptor antagonist with nearly equal affinity for the cloned human EP<sub>1</sub>, EP<sub>2</sub>, EP<sub>3-III</sub>, and DP<sub>1</sub> receptors.<sup>1</sup> AH 6809 blocks the PGE<sub>2</sub>-induced accumulation of cAMP in COS cells transfected with the human EP<sub>2</sub> receptor.<sup>2</sup> It also blocks the accumulation of Ca<sup>2+</sup> in *Xenopus* oocytes expressing the human EP<sub>1</sub> receptor.<sup>3</sup> In the human and guinea pig, the activity profile of AH 6809 is similar to that of SC-19220, but somewhat more potent.<sup>4</sup> In the mouse, AH 6809 has the highest affinity for the EP<sub>2</sub> receptor,<sup>5</sup> but also acts as a weak antagonist at the murine EP<sub>1</sub> and DP<sub>1</sub> receptors. In whole human platelets, AH 6809 is an effective antagonist of the antiaggregatory actions of PGD<sub>2</sub>, but not PGI<sub>2</sub>, with an EC<sub>50</sub> value of about 5 x 10<sup>-5</sup> M.<sup>6</sup>

### References

1. Abramovitz, M., Adam, M., Boie, Y., *et al.* The utilization of recombinant prostanoid receptors to determine the affinities and selectivities of prostaglandins and related analogs. *Biochim. Biophys. Acta* **1483**, 285-293 (2000).
2. Woodward, D.F., Pepperl, D.J., Burkey, T.H., *et al.* 6-Isopropoxy-9-oxoxanthene-2-carboxylic acid (AH 6809), a human EP<sub>2</sub> receptor antagonist. *Biochem. Pharmacol.* **50**, 1731-1733 (1995).
3. Funk, C.D., Furci, L., Fitzgerald, G.A., *et al.* Cloning and expression of a cDNA for the human prostaglandin E receptor EP<sub>1</sub> subtype. *J. Biol. Chem.* **268**, 26767-26772 (1993).
4. Coleman, R.A., Kennedy, I., and Sheldrick, R.L.G. AH6809, a prostanoid EP<sub>1</sub> receptor blocking drug. *Br. J. Pharmacol.* **85**, 273P (1985).
5. Kiriya, M., Ushikubi, F., Kobayashi, T., *et al.* Ligand binding specificities of the eight types and subtypes of the mouse prostanoid receptors expressed in Chinese hamster ovary cells. *Br. J. Pharmacol.* **122**, 217-224 (1997).
6. Keery, R.J. and Lumley, P. AH6809, a prostaglandin DP-receptor blocking drug on human platelets. *Br. J. Pharmacol.* **94**, 745-754 (1988).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/14050](http://www.caymanchem.com/catalog/14050)

**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, 06/20/2012

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)