

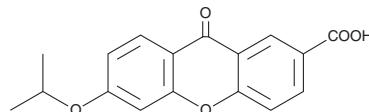
# PRODUCT INFORMATION



**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%  
**UV/Vis.:** λ<sub>max</sub>: 244, 305 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

AH 6809 is supplied as a crystalline solid. A stock solution may be made by dissolving the AH 6809 in the solvent of choice, which should be purged with an inert gas. 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 solid in aqueous buffers. The solubility of AH 6809 in PBS (pH 7.2) is approximately 0.34 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

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**(2), 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**(10), 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**(35), 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**(2), 217-224 (1997).
6. Keery, R.J. and Lumley, P. AH6809, a prostaglandin DP-receptor blocking drug on human platelets. *Br. J. Pharmacol.* **94**(3), 745-754 (1988).

### 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, 11/09/2023

## 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