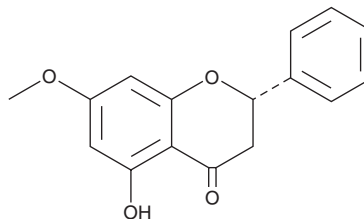


# PRODUCT INFORMATION

## Pinostrobin

Item No. 22436

**CAS Registry No.:** 480-37-5  
**Formal Name:** 2,3-dihydro-5-hydroxy-7-methoxy-2S-phenyl-4H-1-benzopyran-4-one  
**MF:** C<sub>16</sub>H<sub>14</sub>O<sub>4</sub>  
**FW:** 270.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 213, 288 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

Pinostrobin is supplied as a crystalline solid. A stock solution may be made by dissolving the pinostrobin in the solvent of choice. Pinostrobin is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of pinostrobin in these solvents is approximately 12 and 20 mg/ml, respectively.

Pinostrobin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pinostrobin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pinostrobin has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pinostrobin is a flavonoid with diverse biological activities, including antioxidant, anti-inflammatory, and anticancer properties.<sup>1</sup> It induces quinone reductase (QR) in murine hepatoma cells with a QR doubling concentration of 500 nM.<sup>2</sup> Pinostrobin inhibits TNF-α and IL-1β production in RAW 264.7 macrophages (IC<sub>50</sub>s = 17.28 and 23.5 μM, respectively) and in LPS-stimulated rats (48.6 and 55% reduction, respectively).<sup>3</sup> Pinostrobin also shows selective cytotoxicity for CCRF-CEM leukemia cells (IC<sub>50</sub> = 10.2 μM) in a panel of eight cancer cell lines (IC<sub>50</sub>s = >30 μM).<sup>4</sup>

### References

1. Patel, N.K., Jaiswal, G., and Bhutani, K.K. A review on biological sources, chemistry and pharmacological activities of pinostrobin. *Nat. Prod. Res.* **30(18)**, 2017-2027 (2016).
2. Fahey, J.W. and Stephenson, K.K. Pinostrobin from honey and thai ginger (*Boesenbergia pandurata*): A potent flavonoid inducer of mammalian phase 2 chemoprotective and antioxidant enzymes. *J. Agric. Food Chem.* **50(25)**, 7472-7476 (2002).
3. Patel, N.K. and Bhutani, K.K. Pinostrobin and Cajanus lactone isolated from *Cajanus cajan* Millsp. leaves inhibits TNF-α and IL-1 production: *In vitro* and *in vivo* experimentation. *Phytomedicine* **21(7)**, 946-953 (2014).
4. Ashidi, J.S., Houghton, P.J., Hylands, P.J., et al. Ethnobotanical survey and cytotoxicity testing of plants of South-western Nigeria used to treat cancer, with isolation of cytotoxic constituents from *Cajanus cajan* Millsp. leaves. *J. Ethnopharmacol.* **128(2)**, 501-512 (2010).

#### 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, 12/22/2022

#### 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