

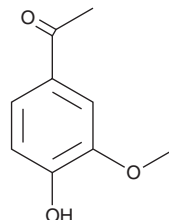
# PRODUCT INFORMATION



## Apocynin

Item No. 11976

**CAS Registry No.:** 498-02-2  
**Formal Name:** 1-(4-hydroxy-3-methoxyphenyl)-ethanone  
**Synonyms:** Acetoguaiacone, Acetovanillone, NSC 2146, NSC 209524  
**MF:** C<sub>9</sub>H<sub>10</sub>O<sub>3</sub>  
**FW:** 166.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 276, 304 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

Apocynin is supplied as a crystalline solid. A stock solution may be made by dissolving the apocynin in the solvent of choice, which should be purged with an inert gas. Apocynin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of apocynin in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

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

### Description

Apocynin is an acetophenone originally isolated from *P. kurroa* that has antioxidant and anti-inflammatory activities.<sup>1,2</sup> It inhibits glucose production in H-4-II-E-C3 cells (IC<sub>50</sub> = 25 μM).<sup>3</sup> Apocynin inhibits the production of superoxide anions induced by PMA (Item No. 10008014) in isolated human leukocytes in a concentration-dependent manner.<sup>1</sup> It increases nitric oxide (NO) bioavailability in rat carotid arterial rings isolated from spontaneously hypertensive stroke-prone (SHRSP) rats when used at concentrations of 0.3 or 3 mM.<sup>4</sup> Apocynin (5 mg/kg) reduces lung injury and pulmonary TNF-α and IL-1β levels in a mouse model of pleurisy induced by carrageenan.<sup>2</sup> It inhibits IκBα degradation and reduces NF-κB p65 and inducible nitric oxide synthase (iNOS) protein levels in the same model. Formulations containing apocynin have been used in the treatment of rheumatoid arthritis, osteoarthritis, bronchial asthma, and chronic obstructive pulmonary disease (COPD).

### References

1. Heumüller, S., Wind, S., Barbosa-Sicard, E., *et al.* Apocynin is not an inhibitor of vascular NADPH oxidases but an antioxidant. *Hypertension* **51(2)**, 211-217 (2008).
2. Hashimoto, J., Motohashi, K., Sakamoto, K., *et al.* Screening and evaluation of new inhibitors of hepatic glucose production. *J. Antibiot. (Tokyo)* **62(11)**, 625-629 (2009).
3. Hamilton, C.A., Brosnan, M.J., Al-Benna, S., *et al.* NAD(P)H oxidase inhibition improves endothelial function in rat and human blood vessels. *Hypertension* **40(5)**, 755-762 (2002).

#### 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, 10/12/2022

#### 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