

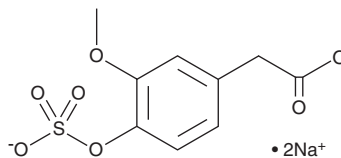
# PRODUCT INFORMATION



## Homovanillic Acid sulfate (sodium salt)

Item No. 20877

**Formal Name:** 3-methoxy-4-(sulfoxy)-benzeneacetic acid, disodium salt  
**MF:** C<sub>9</sub>H<sub>8</sub>O<sub>7</sub>S • 2Na  
**FW:** 306.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 202, 277 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

Homovanillic acid sulfate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the homovanillic acid sulfate (sodium salt) in the solvent of choice. Homovanillic acid sulfate (sodium salt) is soluble in the organic solvent methanol, which should be purged with an inert gas. Homovanillic acid sulfate (sodium salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Homovanillic acid sulfate is a phenol found in olive oil and a dopamine metabolite.<sup>1-3</sup> It is a component of low-density lipoprotein in humans that consume olive oil.<sup>1</sup> Homovanillic acid sulfate levels are increased in response to acute renal injury in humans and rats, and it has been used as a biomarker for early detection of kidney disease in pediatric nephrology.<sup>2,3</sup>

### References

1. de la Torre-Carbot, K., Chávez-Servín, J.L., Jaúregui, O., *et al.* Presence of virgin olive oil phenolic metabolites in human low density lipoprotein fraction: Determination by high-performance liquid chromatography-electrospray ionization tandem mass spectrometry. *Anal. Chim. Acta* **583(2)**, 402-410 (2007).
2. Sun, J., Bhattacharyya, S., Schnackenberg, L.K., *et al.* Discovery of early urinary biomarkers in preclinical study of gentamicin-induced kidney injury and recovery in rats. *Metabolomics* **8(6)**, 1181-1193 (2012).
3. Hanna, M.H. and Brophy, P.D. Metabolomics in pediatric nephrology: Emerging concepts. *Pediatr. Nephrol.* **30(6)**, 881-887 (2015).

#### 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, 12/08/2022

#### 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