

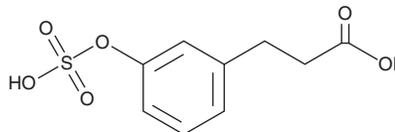
PRODUCT INFORMATION



3-(3-Hydroxyphenyl)propionic Acid sulfate

Item No. 9004103

CAS Registry No.: 86321-32-6
Formal Name: 3-(sulfooxy)-benzenepropanoic acid
Synonym: 3-HPPA sulfate
MF: C₉H₁₀O₆S
FW: 246.2
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

3-(3-Hydroxyphenyl)propionic acid (3-HPPA) sulfate is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-HPPA sulfate in the solvent of choice, which should be purged with an inert gas. 3-HPPA sulfate is soluble in methanol.

Description

3-HPPA sulfate is a metabolite of a variety of phenols, including ferulic acid (Item No. 19871), dihydroferulic acid, and dihydrocaffeic acid (Item No. 28390), and the glycosides echinacoside (Item No. 29927) and acteoside.^{1,2} Urinary excretion of 3-HPPA sulfate is reduced in protein-deficient mice infected with *C. parvum* and increased in mice fed wheat bran or functionalized wheat bran.³

References

1. Ma, Y., Lu, Y., Petrofsky, K., *et al.* Double-edged metabolic effects from short-term feeding of functionalized wheat bran to mouse revealed by metabolomic profiling. *J. Agric. Food Chem* **69**(23), 6543-6555 (2021).
2. Cui, Q., Pan, Y., Bai, X., *et al.* Systematic characterization of the metabolites of echinacoside and acteoside from *Cistanche tubulosa* in rat plasma, bile, urine and feces based on UPLC-ESI-Q-TOF-MS. *Biomed. Chromatogr.* **30**(9), 1406-1415 (2015).
3. Bolick, D.T., Mayneris-Perxachs, J., Medlock, G.L., *et al.* Increased urinary trimethylamine n-oxide following cryptosporidium infection and protein malnutrition independent of microbiome effects. *J. Infect. Dis.* **216**(1), 64-71 (2017).

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, 11/17/2025

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