

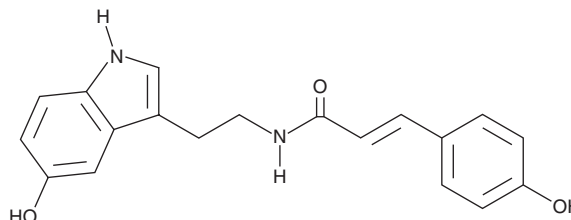
PRODUCT INFORMATION



N-(p-Coumaroyl) Serotonin

Item No. 17797

CAS Registry No.: 68573-24-0
Formal Name: N-[2-(5-hydroxy-1H-indol-3-yl)ethyl]-3-(4-hydroxyphenyl)-2-propenamide
Synonym: NSC 369503
MF: C₁₉H₁₈N₂O₃
FW: 322.4
Purity: ≥95%
UV/Vis.: λ_{max}: 212, 225, 294 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

N-(p-Coumaroyl) serotonin is supplied as a crystalline solid. A stock solution may be made by dissolving the N-(p-coumaroyl) serotonin in the solvent of choice, which should be purged with an inert gas. N-(p-Coumaroyl) serotonin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-(p-coumaroyl) serotonin in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

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

Description

N-(p-Coumaroyl) serotonin is an antioxidative phenolic naturally found in plants, including safflower seed and millet grain.^{1,2} It ameliorates atherosclerosis *in vivo*, reducing lesion area in mice and rabbits.^{1,3} N-(p-Coumaroyl) serotonin also suppresses inflammatory cytokine generation by human monocytes, augments fibroblast proliferation, and promotes relaxation of vascular smooth muscle cells *in vitro*.⁴⁻⁶

References

1. Katsuda, S., Suzuki, K., Koyama, N., et al. *Hypertens. Res.* **32**(11), 944-949 (2009).
2. Watanabe, M. *J. Agric. Food Chem.* **47**(11), 4500-4505 (1999).
3. Koyama, N., Kuribayashi, K., Seki, T., et al. *J. Agric. Food Chem.* **54**(14), 4970-4976 (2015).
4. Kawashima, S., Hayashi, M., Takii, T., et al. *J. Interferon Cytokine Res.* **18**(6), 423-428 (1998).
5. Takii, T., Hayashi, M., Hiroma, H., et al. *J. Biochem.* **125**(5), 910-915 (1999).
6. Takimoto, T., Suzuki, K., Arisaka, H., et al. *Mol. Nutr. Food Res.* **55**(10), 1561-1571 (2011).

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, 11/16/2022

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