

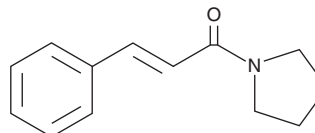
PRODUCT INFORMATION



1-Cinnamoylpyrrolidine

Item No. 36483

CAS Registry No.: 52438-21-8
Formal Name: (2E)-3-phenyl-1-(1-pyrrolidinyl)-2-propen-1-one
MF: $C_{13}H_{15}NO$
FW: 201.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years
Item Origin: Plant/Alkaloids



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

Laboratory Procedures

1-Cinnamoylpyrrolidine is supplied as a solid. A stock solution may be made by dissolving the 1-cinnamoylpyrrolidine in the solvent of choice, which should be purged with an inert gas. 1-Cinnamoylpyrrolidine is soluble in organic solvents such as acetone, chloroform, ethyl acetate, DMSO, and dichloromethane.

Description

1-Cinnamoylpyrrolidine is an amide that has been found in *P. taiwanense* and has antiplatelet and DNA-cleaving activities.^{1,2} It inhibits arachidonic acid- or collagen-induced aggregation of isolated washed rabbit platelets (IC_{50} s = 30.5 and 37.3 μM , respectively).¹ 1-Cinnamoylpyrrolidine (250 μM) cleaves DNA in a cell-free DNA strand scission assay in the presence, but not absence, of copper ions.²

References

1. Chen, I.-S., Chen, Y.-C., and Liao, C.-H. Amides with anti-platelet aggregation activity from *Piper taiwanense*. *Fitoterapia* **78**(6), 414-419 (2007).
2. Ma, J., Jones, S.H., and Hecht, S.M. Phenolic acid amides: A new type of DNA strand scission agent from *Piper caninum*. *Bioorg. Med. Chem.* **12**(14), 3885-3889 (2004).

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, 10/21/2022

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