

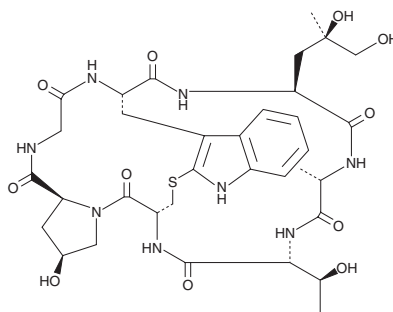
PRODUCT INFORMATION



Phalloidin

Item No. 18039

CAS Registry No.: 17466-45-4
Formal Name: phalloidin
Synonym: NSC 523214
MF: C₃₅H₄₈N₈O₁₁S
FW: 788.9
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 291 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

Phalloidin is supplied as a crystalline solid. A stock solution may be made by dissolving the phalloidin in the solvent of choice. Phalloidin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of phalloidin in these solvents is approximately 10 mg/ml. It is also soluble in warm water. The solubility of phalloidin in warm water is approximately 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Phalloidin is a natural mycotoxin first isolated from the death cap mushroom, *A. phalloides*. It binds filamentous actin (F-actin) and stabilizes actin fibers.^{1,2} Fluorescently labeled phalloidin is commonly used to stain F-actin in cells.^{3,4}

References

1. Lengsfeld, A.M., Löw, I., Wieland, T., *et al.* Interaction of phalloidin with actin. *Proc. Natl. Acad. Sci. USA* **71(7)**, 2803-2807 (1974).
2. Löw, I., Dancker, P., and Wieland, T. Stabilization of F-actin by phalloidin. Reversal of the destabilizing effect of cytochalasin B. *FEBS Lett.* **54(2)**, 263-265 (1975).
3. Wulf, E., Deboben, A., Bautz, F.A., *et al.* Fluorescent phalloidin, a tool for the visualization of cellular actin. *Proc. Natl. Acad. Sci. USA* **76(9)**, 4498-4502 (1979).
4. Jumat, M.R., Nguyen Hong, T., Wong, P., *et al.* Imaging analysis of human metapneumovirus-infected cells provides evidence for the involvement of F-actin and the raft-lipid microdomains in virus morphogenesis. *Virology* **11:198** (2014).

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, 11/17/2022

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