

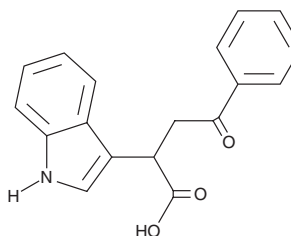
PRODUCT INFORMATION



PEO-IAA

Item No. 40665

CAS Registry No.: 6266-66-6
Formal Name: α -(2-oxo-2-phenylethyl)-1H-indole-3-acetic acid
Synonym: NSC 36840
MF: $C_{18}H_{15}NO_3$
FW: 293.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

PEO-IAA is supplied as a solid. A stock solution may be made by dissolving the PEO-IAA in the solvent of choice, which should be purged with an inert gas. PEO-IAA is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

PEO-IAA is an auxin antagonist.¹ It slows shoot growth and suppresses auxin-induced apical dominance in *C. sativa* seedlings when used at a concentration of 10 $\mu g/L$ in growth media.

Reference

1. Smýkalová, I., Vrbová, M., Cvečková, M., *et al.* The effects of novel synthetic cytokinin derivatives and endogenous cytokinins on the in vitro growth responses of hemp (*Cannabis sativa* L.) explants. Plant Cell Tissue Organ Cult. 139, 381-394 (2019).

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, 04/26/2024

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