

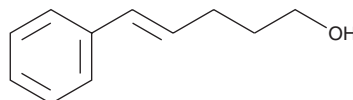
PRODUCT INFORMATION



PPA

Item No. 75751

CAS Registry No.: 13159-16-5
Formal Name: 5-phenyl-4E-pentenol
MF: C₁₁H₁₄O
FW: 162.2
Purity: ≥98%
UV/Vis.: λ_{max}: 252 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

PPA is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of PPA in ethanol and DMF is approximately 10 mg/ml and approximately 7 mg/ml in DMSO.

PPA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of PPA should be diluted with the aqueous buffer of choice. The solubility of PPA in PBS (pH 7.2) is approximately 0.54 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PPA is the product of peroxidase-catalyzed reduction of PPHP. It can be used as a reference standard for HPLC analysis of peroxidase assays.^{1,2}

References

1. Weller, P.E., Markey, C.M., and Marnett, L.J. Enzymatic reduction of 5-phenyl-4-pentenyl-hydroperoxide: Detection of peroxidases and identification of peroxidase reducing substrates. *Arch. Biochem. Biophys.* **243**(2), 633-643 (1985).
2. Shimokawa, T. and Smith, W.L. Prostaglandin endoperoxide synthase. The aspirin acetylation region. *J. Biol. Chem.* **267**(17), 12387-12392 (1992).

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, 02/01/2024

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