

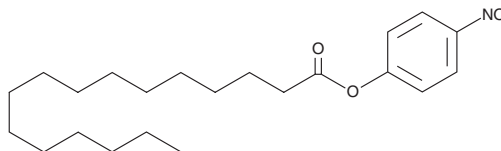
PRODUCT INFORMATION



4-Nitrophenyl Palmitate

Item No. 30870

CAS Registry No.: 1492-30-4
Formal Name: hexadecanoic acid, 4-nitrophenyl ester
Synonyms: *p*-Nitrophenyl Palmitate,
para-Nitrophenyl Palmitate, *p*NPP
MF: C₂₂H₃₅NO₄
FW: 377.5
Purity: ≥95%
UV/Vis.: λ_{max}: 212, 270 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

4-Nitrophenyl palmitate is supplied as a solid. A stock solution may be made by dissolving the 4-nitrophenyl palmitate in the solvent of choice, which should be purged with an inert gas. 4-Nitrophenyl palmitate is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

4-Nitrophenyl palmitate is a colorimetric lipase and esterase substrate.¹ Upon enzymatic hydrolysis of 4-nitrophenyl palmitate, 4-nitrophenol is released which can be quantified by colorimetric detection at 410 nm as a measure of enzyme activity. 4-Nitrophenyl palmitate has been used to characterize the activity of various bacterial and mammalian enzymes, including *Burkholderia* and porcine pancreatic lipases.

Reference

1. Gupta, N., Rathi, P., and Gupta, R. Simplified *para*-nitrophenyl palmitate assay for lipases and esterases. *Anal. Biochem.* **311**(1), 98-99 (2002).

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, 11/10/2020

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