

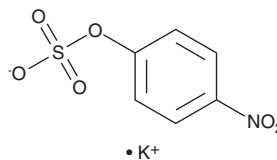
PRODUCT INFORMATION



p-Nitrophenyl Sulfate (potassium salt)

Item No. 35691

CAS Registry No.: 6217-68-1
Formal Name: sulfuric acid, mono(4-nitrophenyl) ester, monopotassium salt
Synonym: 4-Nitrophenyl Sulfate
MF: C₆H₄NO₆S • K
FW: 257.3
Purity: ≥98%
Supplied as: A 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

p-Nitrophenyl sulfate (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the *p*-nitrophenyl sulfate (potassium salt) in the solvent of choice, which should be purged with an inert gas. *p*-Nitrophenyl sulfate (potassium salt) is soluble in acetonitrile.

p-Nitrophenyl sulfate (potassium salt) is soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

p-Nitrophenyl sulfate is a chromogenic substrate for arylsulfatases.¹ Upon enzymatic cleavage by arylsulfatase, *p*-nitrophenyl is released, which can be quantified by colorimetric detection at 400 nm as a measure of arylsulfatase activity.

Reference

- 1 Leon, Y.A., Bulbrook, R.D., and Corner, E.D. Steroid sulphatase, arylsulphatase and beta-glucuronidase in the Mollusca. *Biochem. J.* **75(3)**, 612-617

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, 03/04/2024

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