

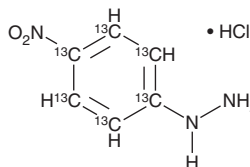
PRODUCT INFORMATION



¹³C₆-4-Nitrophenylhydrazine (hydrochloride)

Item No. 40479

Formal Name:	(4-nitrophenyl-1,2,3,4,5,6- ¹³ C ₆ -4-)hydrazine, monohydrochloride
Synonyms:	¹³ C ₆ -4-NPH, ¹³ C ₆ - <i>p</i> -Nitrophenylhydrazine, ¹³ C ₆ - <i>para</i> -Nitrophenylhydrazine
MF:	[¹³ C] ₆ H ₇ N ₃ O ₂ • HCl
FW:	195.6
Purity:	≥95%
Supplied as:	A solid
Storage:	-20°C
Stability:	≥3 years



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

Laboratory Procedures

¹³C₆-4-Nitrophenylhydrazine is intended for use as an internal standard for the quantification of 4-nitrophenylhydrazine by GC- or LC-MS.

¹³C₆-4-Nitrophenylhydrazine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ¹³C₆-4-nitrophenylhydrazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ¹³C₆-4-Nitrophenylhydrazine (hydrochloride) is soluble in methanol.

Description

¹³C₆-4-Nitrophenylhydrazine is a byproduct in the synthesis of ¹³C₆-3-nitrophenylhydrazine (Item No. 20744). 4-Nitrophenylhydrazine has been used as an intermediate in the synthesis of benzo-indazole derivatives with anti-uterotrophic activity *in vivo* and pyrazolobenzothiazine derivatives with antiviral activity *in vitro*.^{1,2}

References

1. Ragab, M.A., Elagawany, M., Daabees, H., *et al.* Structure-based design and synthesis of conformationally constrained derivatives of methyl-piperidinopyrazole (MPP) with estrogen receptor (ER) antagonist activity. *Bioorg. Chem.* **119**, 105554 (2022).
2. Barreca, M.L., Manfroni, G., Leyssen, P., *et al.* Structure-based discovery of pyrazolobenzothiazine derivatives as inhibitors of hepatitis C virus replication. *J. Med. Chem.* **56**(6), 2270-2282 (2013).

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, 03/20/2024

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