

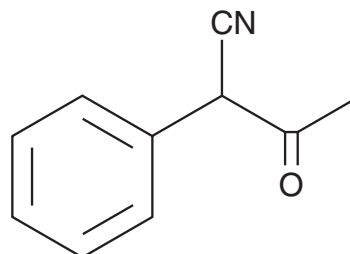
PRODUCT INFORMATION



α -Phenylacetonitrile

Item No. 21105

CAS Registry No.: 4468-48-8
Formal Name: α -acetyl-benzeneacetonitrile
Synonyms: APAAN, 2-Phenylacetoacetonitrile, NSC 11777, NSC 25183, NSC 55206
MF: C₁₀H₉NO
FW: 159.2
Purity: \geq 98%
UV/Vis.: λ_{max} : 264 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

α -Phenylacetonitrile (Item No. 21105) is an analytical reference standard categorized as a precursor in the synthesis of amphetamines.^{1,2} This product is intended for research and forensic applications.

References

1. Power, J.D., Barry, M.G., Scott, K.R., *et al.* An unusual presentation of a customs importation seizure containing amphetamine, possibly synthesized by the APAAN-P2P-Leuckart route. *Forensic Sci. Int.* **234**, 10-13 (2014).
2. Power, J.D., Kavanagh, P., McLaughlin, G., *et al.* Forensic analysis of P2P derived amphetamine synthesis impurities: Identification and characterization of indene by-products. *Drug Test Anal.* Epub ahead of print (2016).

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, 06/22/2023

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