

PRODUCT INFORMATION



N-Boc Norfentanyl

Item No. 36073

CAS Registry No.: 2665695-47-4
Formal Name: 4-[(1-oxopropyl)phenylamino]-

1-piperidinecarboxylic acid,
1,1-dimethylethyl ester
Synonym: t-BOC Norfentanyl

MF: $C_{19}H_{28}N_2O_3$

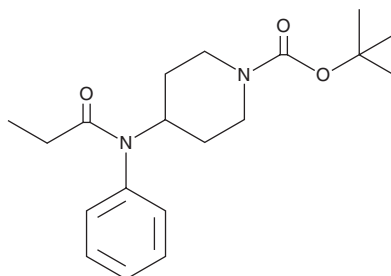
FW: 332.4

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 5 years



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

Description

N-Boc Norfentanyl (Item No. 36073) is an analytical reference standard that is structurally similar to known opioids. N-Boc Norfentanyl is an intermediate in the synthesis of fentanyl-based haptens for use in fentanyl vaccines.¹ This product is intended for research and forensic applications.

Reference

1. Torres, O., Matyas, G.R., Rice, K.C., *et al.* Fentanyl haptens for the preparation of a fentanyl vaccine. *The Henry M. Jackson Foundation for the Advancement of Military Medicine, Inc.* **WO 2021/146334 A1** (2021).

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, 12/07/2023

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