

CERTIFICATE of ANALYSIS



4-Anilino-1-Boc-piperidine

4-(phenylamino)-1-piperidinecarboxylic acid, 1,1-dimethylethyl ester

Item No. 29119 • Batch No. 0572608

Purity Specification: $\geq 98\%$

Molecular Formula : C₁₆H₂₄N₂O₂

CAS Number: 125541-22-2

Formula Weight : 276.37404

Expiry date: 20JAN2023

Tests	Results
HPLC	Purity: 99.0 %
IR	Conforms
Mass spec	MH ⁺ : 276.8
TLC	Purity: 100 %
UV	λ max: 249 nm
NMR	Conforms

Reviewed and approved by: Jennifer LaBrecque

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

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