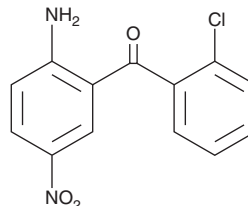


PRODUCT INFORMATION

2-Amino-2'-chloro-5-nitrobenzophenone

Item No. 37885

CAS Registry No.: 2011-66-7
Formal Name: (2-amino-5-nitrophenyl)(2-chlorophenyl)-methanone
Synonyms: 2-Amino-5-nitro-2'-chlorobenzophenone, BF-CLO, BF-Clonazepam, Clonazepam Benzophenone, CLO-Benzophenone
MF: C₁₃H₉ClN₂O₃
FW: 276.7
Purity: 276.7
UV/Vis.: λ_{max}: 228, 353 nm
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

2-Amino-2'-chloro-5-nitrobenzophenone (Item No. 37885) is an analytical reference standard categorized as a precursor in the synthesis of benzodiazepines.¹ 2-Amino-2'-chloro-5-nitrobenzophenone is also an acid hydrolysis product of clonazepam (Item No. 14263).² This product is intended for research and forensic applications.

References

1. Dowling, G., Kavanagh, P.V., Eckhardt, H.-G., *et al.* An approach to shortening the timeframe between the emergence of new compounds on the drugs market and the availability of reference standards: The microscale syntheses of nitrazolam and clonazolam for use as reference materials, utilizing polymer-supported reagents. *Drug Test Anal.* **10**(7), 1198-1208 (2018).
2. Roškar, R. and Sollner Dolenc, M. Determination of benzodiazepines in urine via benzophenone derivatives using liquid chromatography-tandem mass spectrometry. *Arh. Hig. Rada Toksikol.* **61**(4), 381-388 (2010).

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, 07/13/2023

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