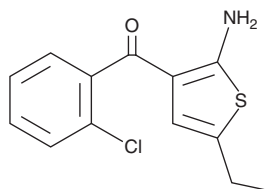


PRODUCT INFORMATION

2-Amino-3-(2-chlorobenzoyl)-5-ethylthiophene

Item No. 36070

CAS Registry No.: 50508-60-6
Formal Name: (2-amino-5-ethyl-3-thienyl)(2-chlorophenyl)-methanone
Synonyms: 2-Amino-5-ethyl-3-(2-chlorobenzoyl)thiophene, 2-Amino-3-(o-chlorobenzoyl)-5-ethylthiophene, Etizolam Related Compound
MF: C₁₃H₁₂ClNOS
FW: 265.8
Purity: ≥98%
UV/Vis.: λ_{max}: 243 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-3-(2-chlorobenzoyl)-5-ethylthiophene (Item No. 36070) is an analytical reference standard categorized as a benzophenone.¹ 2-Amino-3-(2-chlorobenzoyl)-5-ethylthiophene is a precursor in the synthesis of etizolam (Item Nos. 20991 | 12063). It has been found in seized etizolam samples. This product is intended for research and forensic applications.

Reference

- Downey, C., O'Donnell, A., McLaughlin, G., et al. An unusual detection of 2-amino-3-(2-chlorobenzoyl)-5-ethylthiophene and 2-methylamino-5-chlorobenzophenone in illicit yellow etizolam tablets marked "5617" seized in the Republic of Ireland. *Drug Test. Anal.* (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, 04/16/2024

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