

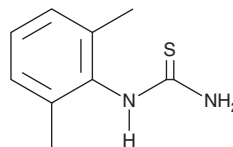
PRODUCT INFORMATION



1-(2,6-Xylyl)-2-thiourea

Item No. 39373

CAS Registry No.: 6396-76-5
Formal Name: N-(2,6-dimethylphenyl)-thiourea
Synonyms: 2,6-Dimethylphenylthiourea,
NSC 202179
MF: C₉H₁₂N₂S
FW: 180.3
Purity: ≥98%
UV/Vis.: λ_{max}: 249 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

1-(2,6-Xylyl)-2-thiourea (Item No. 39373) is an analytical reference standard categorized as an anesthetic metabolite.¹ 1-(2,6-Xylyl)-2-thiourea is a metabolite of xylazine (Item Nos. 37854 | 22641). 1-(2,6-Xylyl)-2-thiourea has been used as a marker for xylazine doping in horses.² This product is intended for research and forensic applications.

References

- Mutlib, A.E., Chui, Y.C., Young, L.M., et al. Characterization of metabolites of xylazine produced in vivo and in vitro by LC/MS/MS and by GC/MS. *Drug Metab. Dispos.* **20(6)**, 840-848 (1992).
- Spyrodaki, M.-H., Lyris, E., Georgoulakis, I., et al. Determination of xylazine and its metabolites by GC-MS in equine urine for doping analysis. *J. Pharm. Biomed. Anal.* **35(1)**, 107-116 (2004).

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, 10/02/2023

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