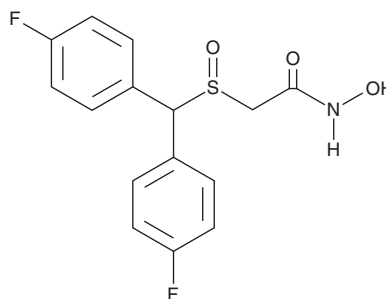


PRODUCT INFORMATION



CRL-40,941
Item No. 19138

CAS Registry No.: 90212-80-9
Formal Name: 2-[[bis(4-fluorophenyl)methyl]sulfinyl]-N-hydroxy-acetamide
Synonyms: bis(4-fluorophenyl)-Adrafinil, Fladrafinil, Fluorafinil
MF: C₁₅H₁₃F₂NO₃S
FW: 325.3
Purity: ≥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

CRL-40,941 (Item No. 19138) is an analytical reference standard categorized as a nootropic.¹ This product is intended for research and forensic applications.

Reference

1. Dowling, G., Kavanagh, P.V., Talbot, B., *et al.* Outsmarted by nootropics? An investigation into the thermal degradation of modafinil, modafinic acid, adrafinil, CRL-40,940 and CRL-40,941 in the GC injector: Formation of 1,1,2,2-tetraphenylethane and its tetra fluoro analog. *Drug Test Anal.* **9(3)**, 518-528 (2017).

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, 04/17/2024

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