

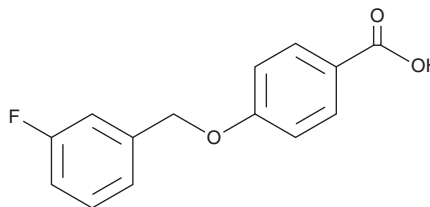
PRODUCT INFORMATION



Safinamide Metabolite NW 1689

Item No. 35395

CAS Registry No.: 405-85-6
Formal Name: 4-[(3-fluorophenyl)methoxy]-benzoic acid
MF: $C_{14}H_{11}FO_3$
FW: 246.2
Purity: $\geq 95\%$
UV/Vis.: $\lambda_{\max}$: 246 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

Safinamide metabolite NW 1689 is supplied as a solid. A stock solution may be made by dissolving the safinamide metabolite NW 1689 in the solvent of choice, which should be purged with an inert gas. Safinamide metabolite NW 1689 is slightly soluble in ethanol.

Description

Safinamide metabolite NW 1689 is a potential impurity found in commercial preparations of safinamide.¹ It is also a degradation product of safinamide (Item No. 21546) formed under conditions of thermal and oxidative stress.

Reference

1. Zou, L., Sun, L., Zhang, H., *et al.* Identification, characterization, and quantification of impurities of safinamide mesilate: Process-related impurities and degradation products. *J. AOAC Int.* **100**(4), 1029-1037 (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, 06/06/2023

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