

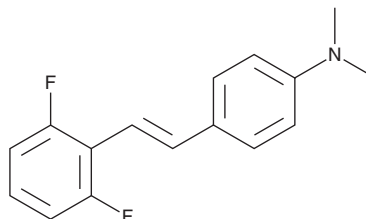
# PRODUCT INFORMATION



## FIDAS-3

Item No. 37716

CAS Registry No.: 1266684-01-8  
Formal Name: 4-[(1E)-2-(2,6-difluorophenyl)ethenyl]-  
N,N-dimethyl-benzenamine  
MF:  $C_{16}H_{15}F_2N$   
FW: 259.3  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 242, 356 nm  
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

FIDAS-3 is supplied as a solid. A stock solution may be made by dissolving the FIDAS-3 in the solvent of choice, which should be purged with an inert gas. FIDAS-3 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

FIDAS-3 is an inhibitor of S-adenosylmethionine synthase isoform type-2 (AdoMet synthase 2;  $IC_{50} = 4.9 \mu M$ ).<sup>1</sup> It also is an inhibitor of human-ether-a-go-go (hERG;  $IC_{50} = 32.1 \mu M$  in a voltage-clamp assay using HEK293 cells expressing hERG).<sup>2</sup> FIDAS-3 reduces the proliferation of LS 174T colorectal cancer cells ( $IC_{50} = 26.9 \mu M$ ) and the levels of Axin2 and c-Myc in the same cells when used at a concentration of 10  $\mu M$ .<sup>1,3</sup> Intraperitoneal administration of FIDAS-3 (20 mg/kg per day) decreases tumor growth in an LS 174T mouse xenograft model.<sup>3</sup>

### References

1. Sviripa, V.M., Zhang, W., Balia, A.G., *et al.* 2',6'-Dihalostyrylanilines, pyridines, and pyrimidines for the inhibition of the catalytic subunit of methionine S-adenosyltransferase-2. *J. Med. Chem.* **57**(14), 6083-6091 (2014).
2. Zhang, W., Sviripa, V., Chen, X., *et al.* Fluorinated N,N-dialkylaminostilbenes repress colon cancer by targeting methionine S-adenosyltransferase 2A. *ACS Chem. Biol.* **8**(4), 796-803 (2013).
3. Zhang, W., Sviripa, V., Kril, L.M., *et al.* Fluorinated N,N-dialkylaminostilbenes for Wnt pathway inhibition and colon cancer repression. *J. Med. Chem.* **54**(5), 1288-1297 (2011).

#### 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, 03/17/2023

#### 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