

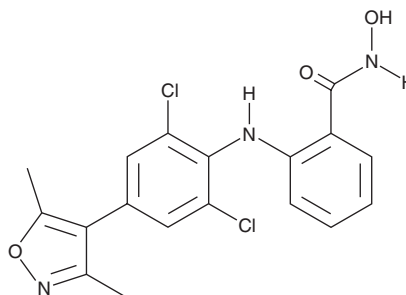
# PRODUCT INFORMATION



## FB23-2

Item No. 36606

**CAS Registry No.:** 2243736-45-8  
**Formal Name:** 2-[[2,6-dichloro-4-(3,5-dimethyl-4-isoxazolyl)phenyl]amino]-N-hydroxy-benzamide  
**MF:** C<sub>18</sub>H<sub>15</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 392.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216 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

FB23-2 is supplied as a solid. A stock solution may be made by dissolving the FB23-2 in the solvent of choice, which should be purged with an inert gas. FB23-2 is slightly soluble in methanol, DMSO, and dimethyl formamide.

### Description

FB23-2 is an inhibitor of the mRNA N<sup>6</sup>-methyladenosine (m<sup>6</sup>A) demethylase fat mass and obesity-associated (FTO) protein (IC<sub>50</sub> = 2.6 μM).<sup>1</sup> It reduces the proliferation of NB4 and Mono-Mac-6 cells (IC<sub>50</sub>s = 0.8 and 1.5 μM, respectively). FB23-2 increases m<sup>6</sup>A levels in isolated human bone marrow cells, as well as suppresses the proliferation of isolated human bone marrow cells and a panel of acute myeloid leukemia (AML) cells (IC<sub>50</sub>s = 1.9-5.2 μM). *In vivo*, FB23-2 (2 mg/kg) delays disease onset and increases survival in a Mono-Mac-6 mouse xenograft model.

### Reference

- Huang, Y., Su, R., Sheng, Y., *et al.* Small-molecule targeting of oncogenic FTO demethylase in acute myeloid leukemia. *Cancer Cell* **35**(4), 677-691.e10 (2019).

#### 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, 12/20/2022

#### 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