

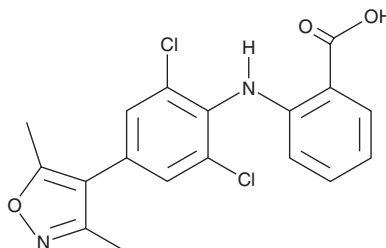
# PRODUCT INFORMATION



## FB23

Item No. 37383

**CAS Registry No.:** 2243736-35-6  
**Formal Name:** 2-[[2,6-dichloro-4-(3,5-dimethyl-4-isoxazolyl)phenyl]amino]-benzoic acid  
**MF:** C<sub>18</sub>H<sub>14</sub>Cl<sub>2</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 377.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221 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 is supplied as a solid. A stock solution may be made by dissolving the FB23 in the solvent of choice, which should be purged with an inert gas. FB23 is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

### Description

FB23 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> = 0.06 μM) and an active metabolite of FB23-2 (Item No. 36606).<sup>1</sup> It reduces the proliferation of NB4 and Mono-Mac-6 acute myeloid leukemia (AML) cells (IC<sub>50</sub>s = 44.8 and 23.6 μM, respectively).

### Reference

1. 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, 09/25/2023

#### 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