

# PRODUCT INFORMATION



## iBET-BD2

Item No. 31766

**CAS Registry No.:** 2474876-09-8  
**Formal Name:** 4-(acetylamino)-3-fluoro-N-(trans-4-hydroxycyclohexyl)-5-[(1S)-1-phenylethoxy]-benzamide

**Synonym:** GSK046  
**MF:** C<sub>23</sub>H<sub>27</sub>FN<sub>2</sub>O<sub>4</sub>

**FW:** 414.5

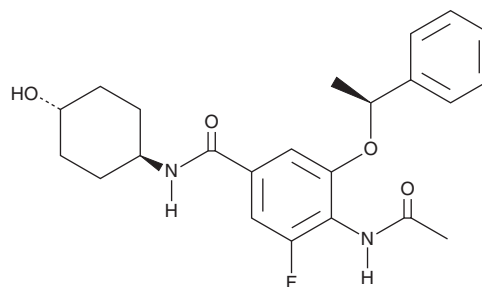
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 253 nm

**Supplied as:** A crystalline 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

iBET-BD2 is supplied as a crystalline solid. A stock solution may be made by dissolving the iBET-BD2 in the solvent of choice, which should be purged with an inert gas. iBET-BD2 is soluble in the organic solvent DMSO.

### Description

iBET-BD2 is an inhibitor of the bromodomain and extra-terminal domain (BET) family protein bromodomain 2 (BD2).<sup>1</sup> It inhibits BD2 in bromodomain-containing protein 2 (BRD2), BRD3, BRD4, and BRDT (IC<sub>50</sub>s = 264, 98, 49, and 214 nM, respectively) and is selective for BD2 over BD1 (IC<sub>50</sub>s = 10.96, 36.31, 70.55, and >50.11 μM for BRD2, BRD3, BRD4, and BDT, respectively) in a time-resolved FRET (TR-FRET) assay. iBET-BD2 (1 μM) inhibits IFN-γ-induced protein expression of MHC class I in K562 cells. It reduces the production of anti-KLH IgM antibodies in mice immunized with keyhole limpet hemocyanin (KLH) when administered at a dose of 40 mg/kg.

### Reference

1. Gilan, O., Rioja, I., Knezevic, K., et al. Selective targeting of BD1 and BD2 of the BET proteins in cancer and immuno-inflammation. *Science* **368**(6489), 387-394 (2020).

#### 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, 01/03/2023

#### 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