

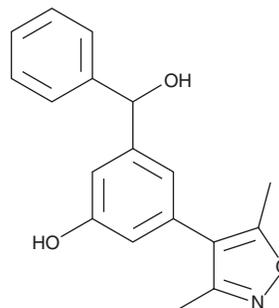
# PRODUCT INFORMATION



## OXF BD 02

Item No. 14811

**CAS Registry No.:** 1429129-68-9  
**Formal Name:** 3-(3,5-dimethyl-4-isoxazolyl)-5-hydroxy- $\alpha$ -phenyl-benzenemethanol  
**MF:** C<sub>18</sub>H<sub>17</sub>NO<sub>3</sub>  
**FW:** 295.3  
**Purity:** ≥98%  
**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

OXF BD 02 is supplied as a solid. A stock solution may be made by dissolving the OXF BD 02 in the solvent of choice, which should be purged with an inert gas. OXF BD 02 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

OXF BD 02 is an inhibitor of bromodomain 1 (BD1) in bromodomain-containing protein 4 (BRD4; IC<sub>50</sub> = 382 nM).<sup>1</sup> It is selective for BRD4 BD1 over a panel of 60 bromodomain-containing proteins but does inhibit CREB-binding protein (CBP) at 25 μM. OXF BD 02 is cytotoxic to MV4-11 acute myeloid leukemia (AML) cells (IC<sub>50</sub> = 0.794 nM) but not A549 or H1975 lung cancer cells (IC<sub>50</sub>s = >10 μM for both) or U2OS osteosarcoma or HeLa cervical cancer cells (IC<sub>50</sub>s = >100 μM for both). OXF BD 02 also binds to recombinant *S. mansoni* BRD3 (SmBRD3) that contains BD1 and BD2 (K<sub>d</sub> = 683 nM) but not recombinant BD2 of SmBRD3.<sup>2</sup>

### References

- Hewings, D.S., Fedorov, O., Filippakopoulos, P., *et al.* Optimization of 3,5-dimethylisoxazole derivatives as potent bromodomain ligands. *J. Med. Chem.* **56**(8), 3217-3227 (2013).
- Schiedel, M., McArdle, D.J.B., Padalino, G., *et al.* Small molecule ligands of the BET-like bromodomain, SmBRD3, affect *Schistosoma mansoni* survival, oviposition, and development. *J. Med. Chem.* **66**(23), 15801-15822 (2023).

#### 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, 08/12/2025

#### 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