

# PRODUCT INFORMATION

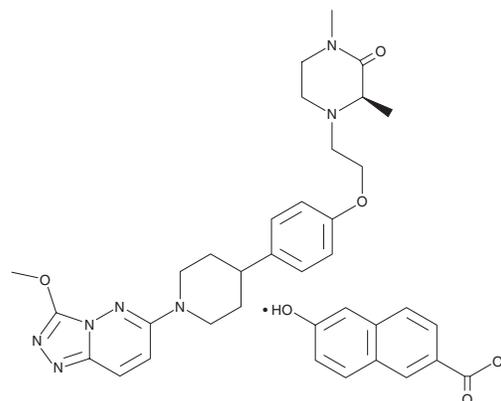


## AZD 5153 (6-hydroxy-2-naphthoic acid)

Item No. 42341

**CAS Registry No.:** 1869912-40-2  
**Formal Name:** 6-hydroxy-2-naphthalenecarboxylic acid, compd. with (3R)-4-[2-[4-[1-(3-methoxy-1,2,4-triazolo[4,3-b]pyridazin-6-yl)-4-piperidinyl]phenoxy]ethyl]-1,3-dimethyl-2-piperazinone (1:1)

**Synonym:** AZD 5153 HNT salt  
**MF:** C<sub>25</sub>H<sub>33</sub>N<sub>7</sub>O<sub>3</sub> • C<sub>11</sub>H<sub>8</sub>O<sub>3</sub>  
**FW:** 667.8  
**Purity:** ≥95%  
**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

AZD 5153 (6-hydroxy-2-naphthoic acid) is supplied as a solid. A stock solution may be made by dissolving the AZD 5153 (6-hydroxy-2-naphthoic acid) in the solvent of choice, which should be purged with an inert gas. AZD 5153 (6-hydroxy-2-naphthoic acid) is soluble (≥10 mg/ml) in DMSO.

### Description

AZD 5153 is an orally available, bivalent inhibitor of the bromodomain and extraterminal (BET) protein BRD4 (IC<sub>50</sub> = 5 nM).<sup>1</sup> It can simultaneously bind two bromodomains in BRD4, which has been shown to increase antitumor activity in multiple xenograft models of acute myeloid leukemia, multiple myeloma, and diffuse large B cell lymphoma.

### Reference

1. Rhyasen, G.W., Hattersley, M.M., Yao, Y., *et al.* AZD5153: A novel bivalent BET bromodomain inhibitor highly active against hematologic malignancies. *Mol. Cancer Ther.* **15(11)**, 2563-2574 (2016).

**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, 11/15/2024

### 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