

# PRODUCT INFORMATION

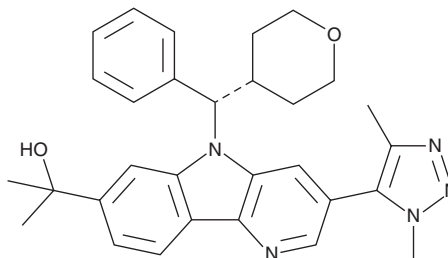


## BMS 986158

Item No. 27789

**CAS Registry No.:** 1800340-40-2  
**Formal Name:** 3-(1,4-dimethyl-1H-1,2,3-triazol-5-yl)- $\alpha,\alpha$ -dimethyl-5-[(S)-phenyl(tetrahydro-2H-pyran-4-yl)methyl]-5H-pyrido[3,2-b]indole-7-methanol

**MF:** C<sub>30</sub>H<sub>33</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 495.6  
**Purity:** ≥98%  
**UV/Vis.:**  $\lambda_{\text{max}}$ : 214, 270, 322 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

BMS 986158 is supplied as a solid. A stock solution may be made by dissolving the BMS 986158 in the solvent of choice, which should be purged with an inert gas. BMS 986158 is soluble in the organic solvent DMSO at a concentration of approximately 50 mM.

### Description

BMS 986158 is a BET bromodomain inhibitor with anticancer activity.<sup>1</sup> It decreases c-Myc expression and induces cancer cell death in c-Myc-driven cancer cell lines *in vitro*. BMS 986158 inhibits tumor growth in triple-negative breast, colorectal, and lung cancer patient-derived mouse xenograft models.

### Reference

1. Gavai, A.V., Norris, D., Tortolani, D., *et al.* Abstract 5789: Discovery of clinical candidate BMS-986158, an oral BET inhibitor, for the treatment of cancer. *Cancer Res.* **78(Suppl. 13)**, 5789-5789 (2018).

#### 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, 11/22/2022

#### 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