

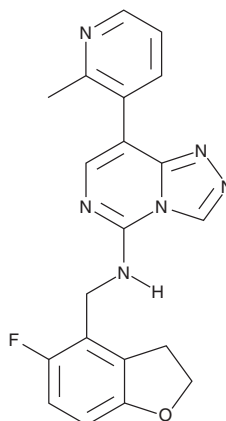
# PRODUCT INFORMATION



## MAK-683

Item No. 29316

**CAS Registry No.:** 1951408-58-4  
**Formal Name:** N-[(5-fluoro-2,3-dihydro-4-benzofuranyl)methyl]-8-(2-methyl-3-pyridinyl)-1,2,4-triazolo[4,3-c]pyrimidin-5-amine  
**MF:** C<sub>20</sub>H<sub>17</sub>FN<sub>6</sub>O  
**FW:** 376.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 298 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

MAK-683 is supplied as a solid. A stock solution may be made by dissolving the MAK-683 in the solvent of choice, which should be purged with an inert gas. MAK-683 is soluble in the organic solvent DMSO at a concentration of approximately 15 mg/ml.

### Description

MAK-683 is an inhibitor of trimethylation of histone 3 lysine 27 (H3K27) by polycomb repressive complex 2 (PRC2), a complex that includes embryonic ectoderm development protein (EED).<sup>1</sup> MAK-683 has an IC<sub>50</sub> value of 5.9 nM in an EED-H3K27me3 peptide competition binding assay and inhibits H3K27 trimethylation in G401 cells (IC<sub>50</sub> = 2.6 nM). It inhibits proliferation of KARPAS422 B cell lymphoma cells with an IC<sub>50</sub> value of 3 nM.

### Reference

1. Chan, H.M., Gu, X.-J., Huang, Y., *et al.* Triazolopyrimidine compounds and uses thereof. *Novartis AG. US20160176882A1* (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, 12/08/2022

#### 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