

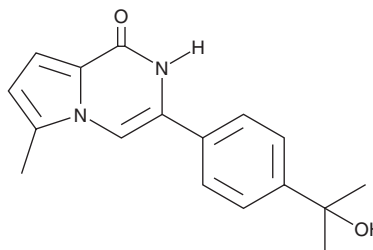
PRODUCT INFORMATION



M2912

Item No. 45586

CAS Registry No.: 1460286-21-8
Formal Name: 3-[4-(1-hydroxy-1-methylethyl)phenyl]-6-methyl-pyrrolo[1,2-a]pyrazin-1(2H)-one
Synonym: MSC2504877
MF: C₁₇H₁₈N₂O₂
FW: 282.3
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

M2912 is supplied as a solid. A stock solution may be made by dissolving the M2912 in the solvent of choice, which should be purged with an inert gas. M2912 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

M2912 is an inhibitor of tankyrase 1 (TNKS1) and TNKS2 (IC₅₀s = 0.6 and 0.9 nM, respectively).¹ It is selective for TNKS1 and TNKS2 over poly(ADP-ribose) polymerase 1 (PARP1; IC₅₀ = 928 nM). M2912 increases the levels of Axin2 (EC₅₀ = 17 nM) and reduces Wnt reporter activity in DLD-1 cells. It reduces the growth of a variety of colorectal cancer cells *in vitro* at 0.01-10 μM and reduces tumor volume in a COLO 320 mouse xenograft model at 20 mg/kg.

Reference

1. Buchstaller, H.P., Anlauf, U., Dorsch, D., *et al.* Optimization of a screening hit toward M2912, an oral tankyrase inhibitor with antitumor activity in colorectal cancer models. *J. Med. Chem.* **64**(14), 10371-10392 (2021).

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, 05/29/2026

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