

PRODUCT INFORMATION



JW 55

Item No. 16260

CAS Registry No.: 664993-53-7
Formal Name: N-[4-[[[tetrahydro-4-(4-methoxyphenyl)-2H-pyran-4-yl]methyl]amino]carbonyl]phenyl]-2-furancarboxamide

Synonym: Tankyrase 1/2 Inhibitor IV

MF: C₂₅H₂₆N₂O₅

FW: 434.5

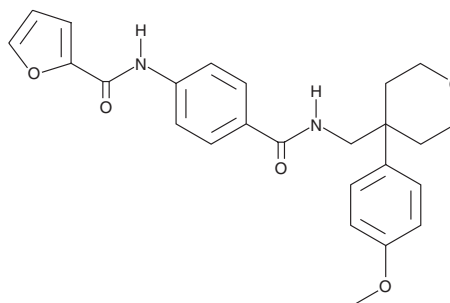
Purity: ≥98%

UV/Vis.: λ_{max}: 291 nm

Supplied as: A crystalline 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

JW 55 is supplied as a crystalline solid. A stock solution may be made by dissolving the JW 55 in the solvent of choice, which should be purged with an inert gas. JW 55 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JW 55 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

JW 55 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JW 55 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JW 55 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

JW 55 is an inhibitor of the PARP domain of tankyrases (TNKS) 1 and 2 (IC₅₀s = 1.9 and 0.83 μM, respectively).¹ By inhibiting TNKS1/2, JW 55 stabilizes Axin2 and increases degradation of β-catenin, inhibiting the β-catenin signaling pathway (IC₅₀ = 470 nM).¹ It has been shown to decrease canonical Wnt signaling in SW480 and HCT15 colon carcinoma cell lines and to reduce proliferation of SW480 cells *in vitro*.¹

Reference

1. Waaler, J., Machon, O., Tumova, L., *et al.* A novel tankyrase inhibitor decreases canonical Wnt signaling in colon carcinoma cells and reduces tumor growth in conditional APC mutant mice. *Cancer Res.* **72(11)**, 2822-2832 (2012).

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, 11/14/2022

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