

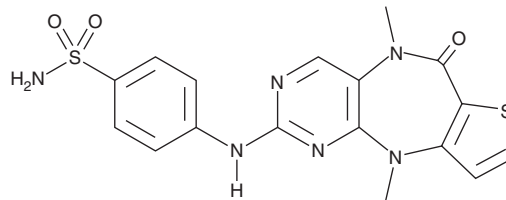
# PRODUCT INFORMATION



## XMU-MP-1

Item No. 22083

**CAS Registry No.:** 2061980-01-4  
**Formal Name:** 4-[(6,10-dihydro-5,10-dimethyl-6-oxo-5H-pyrimido[5,4-b]thieno[3,2-e][1,4]diazepin-2-yl)amino]-benzenesulfonamide  
**MF:** C<sub>17</sub>H<sub>16</sub>N<sub>6</sub>O<sub>3</sub>S<sub>2</sub>  
**FW:** 416.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 299 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



**Special Conditions:** Requires sonication

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

XMU-MP-1 is supplied as a solid. A stock solution may be made by dissolving the XMU-MP-1 in the solvent of choice, which should be purged with an inert gas. XMU-MP-1 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of XMU-MP-1 in these solvents is approximately 0.5 and 2 mg/ml, respectively.

### Description

XMU-MP-1 is an inhibitor of the pro-apoptotic, sterile 20-like kinases MST1 and 2 (IC<sub>50</sub>s = 71.1 and 38.1 nM, respectively).<sup>1</sup> It is selective for MST1 and 2 over a panel of 468 kinases at a concentration of 1 μM. XMU-MP-1 inhibits phosphorylation of MOB1, LATS1 and 2, and YAP in HepG2 cells as well as MST1 and 2 autophosphorylation in RAW 264.7, U2OS, SW480, RPE1, SNU423, and HepG2 cells in a concentration-dependent manner. It also prevents cell death induced by overexpression of MST2. XMU-MP-1 (1 mg/kg) enhances liver regeneration in mice post partial hepatectomy and reduces liver fibrosis in a mouse model of chronic liver injury.

### Reference

1. Fan, F., He, Z., Kong, L.-L., *et al.* Pharmacological targeting of kinases MST1 and MST2 augments tissue repair and regeneration. *Sci. Transl. Med.* **8**(352), 352ra108 (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, 12/12/2022

#### 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