

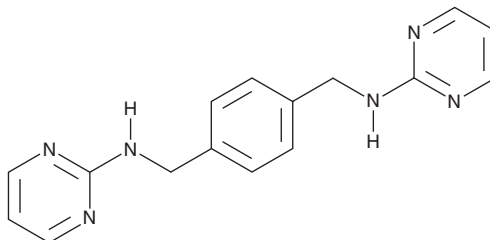
PRODUCT INFORMATION



MSX-122

Item No. 23426

CAS Registry No.: 897657-95-3
Formal Name: N¹,N⁴-di-2-pyrimidinyl-1,4-benzenedimethanamine
MF: C₁₆H₁₆N₆
FW: 292.3
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 302 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

MSX-122 is supplied as a crystalline solid. A stock solution may be made by dissolving the MSX-122 in the solvent of choice, which should be purged with an inert gas. MSX-122 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of MSX-122 in these solvents is approximately 0.2 mg/ml.

Description

MSX-122 is a small molecule partial antagonist of C-X-C chemokine 4 (CXCR4; IC₅₀ = ~10 nM).¹ It reduces inflammation and infiltration of CXCR4⁺ cells in a dextran sulfate (Item No. 23250) model of experimental colitis in mice. Specifically, it prevents the production of TNF-α, IL-1β, and IL-6. MSX-122 also reduces TNF-α secretion by macrophages infected with patient-derived invasive *E. coli*. It decreases metastases in mouse models utilizing MDA-MB-231 breast cancer or 686LN-Ms squamous cell carcinoma cells injected intravenously or OMM2.3 melanoma cells inoculated into the eye.

Reference

1. Liang, Z., Zhan, W., Zhu, A., *et al.* Development of a unique small molecule modulator of CXCR4. *PLoS One* 7(4), e34038 (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, 10/10/2022

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