

## PRODUCT INFORMATION

## MELK-8a (hydrochloride)

Item No. 37025

**CAS Registry No.:** 2096992-20-8

**Formal Name:** 1-methyl-4-[4-[4-[3-(4-piperidinylmethoxy)-4-pyridinyl]-1H-pyrazol-1-yl]phenyl]-piperazine, monohydrochloride

**Synonym:** NVS-MELK8a

MF:  $C_{25}H_{22}N_4O \bullet HCl$

FW: 469.0

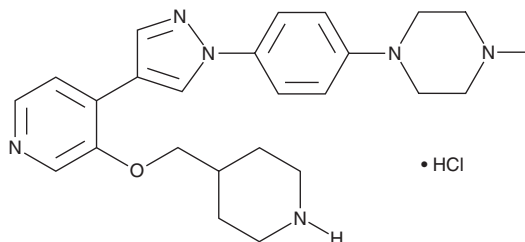
Purity:  $\geq 95\%$ 

UV/Vis.:  $\lambda_{\text{max}}$ : 307 nm

Supplied as: <sup>max</sup>  
A solid

**Storage:** -20°C

**Stability:**  $\geq 4$  years



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

## Laboratory Procedures

MELK-8a (hydrochloride) is supplied as a solid. Aqueous solutions of MELK-8a (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of MELK-8a (hydrochloride) in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

MELK-8a is an inhibitor of maternal embryonic leucine zipper kinase (MELK;  $IC_{50} = 0.002 \mu M$ ).<sup>1</sup> It is selective for MELK over FLT3, haspin, KIT, and PDGFR $\alpha$ , ( $IC_{50}$ s = 0.18, 0.19, >10, and  $0.42 \mu M$ , respectively), as well as a panel of 456 kinases at  $1 \mu M$ . MELK-8a inhibits the proliferation of MDA-MB-468 and MCF-7 breast cancer cells ( $IC_{50}$ s = 0.11 and  $3.68 \mu M$ ).

## Reference

1. Touré, B.B., Giraldez, J., Smith, T., *et al.* Toward the validation of maternal embryonic leucine zipper kinase: Discovery, optimization of highly potent and selective inhibitors, and preliminary biology insight. *J. Med. Chem.* **59**(10), 4711-4723 (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, 12/21/2022

**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