

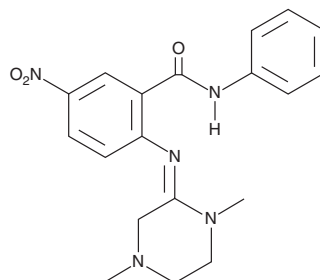
# PRODUCT INFORMATION



## ML-336

Item No. 9001920

**CAS Registry No.:** 1613465-33-0  
**Formal Name:** 2-[(E)-(1,4-dimethyl-2-piperazinylidene)amino]-5-nitro-N-phenyl-benzamide  
**MF:** C<sub>19</sub>H<sub>21</sub>N<sub>5</sub>O<sub>3</sub>  
**FW:** 367.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 274, 345 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

ML-336 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-336 in the solvent of choice. ML-336 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of ML-336 in these solvents is approximately 0.2, 10, and 20 mg/ml, respectively.

ML-336 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-336 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ML-336 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Venezuelan Equine Encephalitis Virus (VEEV) is a mosquito-borne alphavirus that can cause encephalomyelitis in all equine species and also in humans. ML-336 is a quinazolinone-based compound that has been shown to inhibit a VEEV-induced cytopathic effect in three strains of the virus (IC<sub>50</sub>s = 32, 20, and 41 nM for TC-83, V3526, and Trinidad donkey strains, respectively).<sup>1</sup> It is reported to target the VEEV non-structural protein 2, which is necessary for transcription and replication of viral RNA.<sup>1</sup>

### Reference

1. Chung, D., Schroeder, C.E., Sotsky, J., *et al.* ML336: Development of quinazolinone-based inhibitors against Venezuelan Equine Encephalitis Virus (VEEV), in *Probe Reports from the NIH Molecular Libraries Program*, 1 (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, 11/09/2022

#### 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