

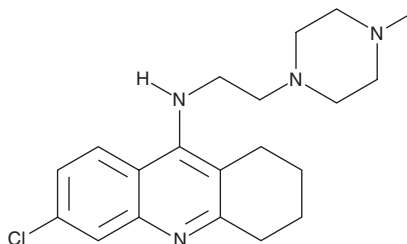
# PRODUCT INFORMATION



## Quinacrine analog 34

Item No. 17733

**CAS Registry No.:** 1411646-44-0  
**Formal Name:** 6-chloro-1,2,3,4-tetrahydro-N-[2-(4-methyl-1-piperazinyl)ethyl]-9-acridinamine  
**MF:** C<sub>20</sub>H<sub>27</sub>ClN<sub>4</sub>  
**FW:** 358.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 226, 252, 330 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

Quinacrine analog 34 is supplied as a crystalline solid. A stock solution may be made by dissolving the quinacrine analog 34 in the solvent of choice, which should be purged with an inert gas. Quinacrine analog 34 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Quinacrine analog 34 in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Quinacrine analog 34 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, quinacrine analog 34 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Quinacrine analog 34 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

Quinacrine analog 34 is an inhibitor of autophagy with a minimum effective concentration (MEC) of 0.5 μM in U2OS cells.<sup>1</sup> It induces lysosomal deacidification and accumulation of LC3-II in the same cells when used at concentrations of 0.5 or 3 μM.

### Reference

1. Wang, T., Goodall, M.L., Gonzales, P., *et al.* Synthesis of improved lysomotropic autophagy inhibitors. *J. Med. Chem.* **58**(7), 3025-3035 (2015).

#### 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, 02/15/2024

#### 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