

# PRODUCT INFORMATION

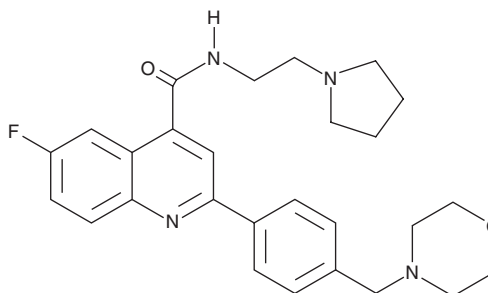


**DDD107498**

Item No. 18305

**CAS Registry No.:** 1469439-69-7  
**Formal Name:** 6-fluoro-2-[4-(4-morpholinylmethyl)phenyl]-N-[2-(1-pyrrolidinyl)ethyl]-4-quinolinecarboxamide

**Synonym:** DDD 498  
**MF:** C<sub>27</sub>H<sub>31</sub>FN<sub>4</sub>O<sub>2</sub>  
**FW:** 462.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 268, 336 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

## Laboratory Procedures

DDD107498 is supplied as a solid. A stock solution may be made by dissolving the DDD107498 in the solvent of choice, which should be purged with an inert gas. DDD107498 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DDD107498 in ethanol and DMF is approximately 30 mg/ml and approximately 10 mg/ml in DMSO.

## Description

DDD107498 is an antimalarial.<sup>1</sup> It is an inhibitor of *P. falciparum* translation elongation factor 2 (PfeEF2) that inhibits protein synthesis. DDD107498 is active against *P. falciparum* strain 3D7 (EC<sub>50</sub> = 1 nM) and clinical isolates of *P. falciparum* and *P. vivax* isolated from patients in areas where multidrug-resistant malaria is endemic (median EC<sub>50</sub>s = 0.81 and 0.51 nM, respectively). It reduces parasitemia in a mouse model of *P. berghei* infection (ED<sub>90</sub> = 0.57 mg/kg), indicating activity against the blood-stage form of the parasite. DDD107498 is also active against liver schizont forms of *P. berghei* and *P. yoelii* (EC<sub>50</sub> = ~1 nM) and eradicates *P. berghei* infection in mice when administered prior to infection at a dose of 3 mg/kg.

## Reference

1. Baragaña, B., Hallyburton, I., Lee, M.C.S., *et al.* A novel multiple-stage antimalarial agent that inhibits protein synthesis. *Nature* 522(7556), 315-331 (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, 10/04/2021

## 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