

# PRODUCT INFORMATION



**MF498**

Item No. 15973

**CAS Registry No.:** 915191-42-3

**Formal Name:** N-[[[4-(5,9-diethoxy-6-oxo-7H-pyrrolo[3,4-g]quinolin-7-yl)-3-methylphenyl]methyl]sulfonyl]-2-methoxy-benzeneacetamide

**MF:** C<sub>32</sub>H<sub>33</sub>N<sub>3</sub>O<sub>7</sub>S

**FW:** 603.7

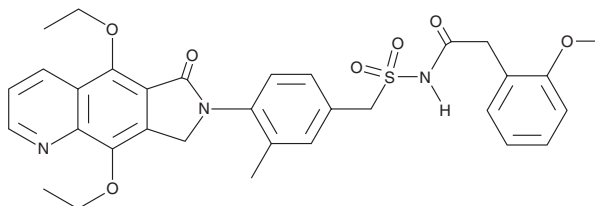
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 204, 255 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

MF498 is supplied as a crystalline solid. A stock solution may be made by dissolving the MF498 in the solvent of choice, which should be purged with an inert gas. MF498 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MF498 in ethanol is approximately 20 mg/ml and approximately 10 mg/ml in DMSO and DMF.

MF498 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MF498 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. MF498 has a solubility of approximately 0.09 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) activates four E prostanoid (EP) receptors, EP<sub>1-4</sub>. EP<sub>4</sub> is a Gs protein-coupled receptor that, by elevating the second messenger cAMP, plays important roles in bone formation and resorption, inflammation, cancer, and atherosclerosis.<sup>1-3</sup> MF498 is a selective EP<sub>4</sub> receptor antagonist (K<sub>i</sub> = 0.7 nM versus a K<sub>i</sub> > 1 μM for other EP receptors).<sup>4</sup> In HEK293 cells expressing the human EP<sub>4</sub> receptor, MF498 inhibits EP ligand induced activity with an IC<sub>50</sub> value of 1.7 nM.<sup>4</sup> In various animal models for arthritis, MF498 has been shown to inhibit inflammation without gastrointestinal toxicity (ED<sub>50</sub> values as low as 0.02 mg/kg/day).<sup>4</sup>

## References

1. Li, M., Thompson, D.D., and Paralkar, V.M. Prostaglandin E<sub>2</sub> receptors in bone formation. *Int. Orthop.* **31(6)**, 767-772 (2007).
2. Hawcroft, G., Ko, C.W.S., and Hull, M.A. Prostaglandin E<sub>2</sub>-EP4 receptor signalling promotes tumorigenic behaviour of HT-29 human colorectal cancer cells. *Oncogene* **26(21)**, 3006-3019 (2007).
3. Babaev, V.R., Chew, J.D., Ding, L., et al. Macrophage EP4 deficiency increases apoptosis and suppresses early atherosclerosis. *Cell Metab.* **8(6)**, 492-501 (2008).
4. Clark, P., Rowland, S.E., Denis, D., et al. MF498 [N-[[[4-(5,9-diethoxy-6-oxo-6,8-dihydro-7H-pyrrolo[3,4-g]quinolin-7-yl)-3-methylbenzyl]sulfonyl]-2-(2-methoxyphenyl)acetamide], a selective E prostanoid receptor 4 antagonist, relieves joint inflammation and pain in rodent models of rheumatoid and osteoarthritis. *J. Pharmacol. Exp. Ther.* **325(2)**, 425-434 (2008).

### 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, 11/30/2022

## 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