

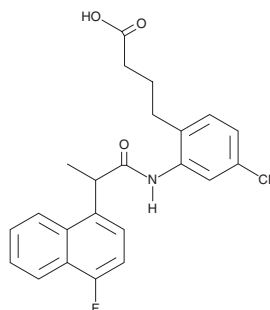
# PRODUCT INFORMATION



## ONO-AE3-208

Item No. 14522

**CAS Registry No.:** 402473-54-5  
**Formal Name:** 4-cyano-2-[[2-(4-fluoro-1-naphthalenyl)-1-oxopropyl]amino]-benzenebutanoic acid  
**MF:** C<sub>24</sub>H<sub>21</sub>FN<sub>2</sub>O<sub>3</sub>  
**FW:** 404.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 286 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

ONO-AE3-208 is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-AE3-208 in the solvent of choice, which should be purged with an inert gas. ONO-AE3-208 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ONO-AE3-208 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

ONO-AE3-208 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ONO-AE3-208 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ONO-AE3-208 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

ONO-AE3-208 is an antagonist of the EP<sub>4</sub> receptor (K<sub>i</sub> = 1.3 nM) which less potently affects EP<sub>3</sub>, FP, and TP receptors (K<sub>i</sub>s = 30, 790, and 2,400 nM, respectively) and is without effect on other prostanoid receptors.<sup>1,2</sup> In wild type mice, it mimics deletion of EP<sub>4</sub> by producing severe colitis, with epithelial loss, crypt damage, and inflammation, after treatment with 3% dextran sodium sulfate.<sup>1</sup> ONO-AE3-208 has also been used to implicate EP<sub>4</sub> signaling in immune and autoimmune responses, inflammation, and cancer.<sup>3-6</sup>

### References

1. Kabashima, K., Saji, T., Murata, T., *et al.* The prostaglandin receptor EP<sub>4</sub> suppresses colitis, mucosal damage and CD4 cell activation in the gut. *J. Clin. Invest.* **109**(7), 883-893 (2002).
2. Jones, R.L., Giembycz, M.A., and Woodward, D.F. Prostanoid receptor antagonists: Development strategies and therapeutic applications. *Br. J. Pharmacol.* **158**(1), 104-145 (2009).
3. Ma, X., Kundu, N., Rifat, S., *et al.* Prostaglandin E receptor EP<sub>4</sub> antagonism inhibits breast cancer metastasis. *Cancer Res.* **66**(6), 2923-2927 (2006).
4. Kabashima, K., Sakata, D., Nagamachi, M., *et al.* Prostaglandin E<sub>2</sub>-EP<sub>4</sub> signaling initiates skin immune responses by promoting migration and maturation of Langerhans cells. *Nat. Med.* **9**(6), 744-749 (2003).
5. Esaki, Y., Li, Y., Sakata, D., *et al.* Dual roles of PGE<sub>2</sub>-EP<sub>4</sub> signaling in mouse experimental autoimmune encephalomyelitis. *Proc. Natl. Acad. Sci. USA* **107**(27), 12233-12238 (2010).
6. Komatsu, H., Enjouji, S., Ito, A., *et al.* Prostaglandin E<sub>2</sub> inhibits proteinase-activated receptor 2-signal transduction through regulation of receptor internalization. *J. Vet. Med. Sci.* **75**(3), 255-261 (2013).

#### 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, 04/02/2024

#### 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