

PRODUCT INFORMATION



ONO-AE3-208 (sodium salt)

Item No. 16108

CAS Registry No.: 2309931-05-1
Formal Name: 4-cyano-2-[[2-(4-fluoro-1-naphthalenyl)-1-oxopropyl]amino]-benzenebutanoic acid, monosodium salt

MF: C₂₄H₂₀FN₂O₃ • Na

FW: 426.4

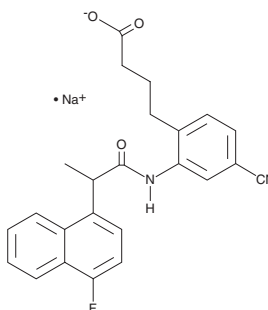
Purity: ≥95%

UV/Vis.: λ_{max}: 223, 285 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 (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-AE3-208 (sodium salt) in the solvent of choice, which should be purged with an inert gas. ONO-AE3-208 (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ONO-AE3-208 (sodium salt) in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ONO-AE3-208 (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ONO-AE3-208 (sodium salt) in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ONO-AE3-208 is an antagonist of the EP₄ receptor (K_i = 1.3 nM) that less potently affects EP₃, FP, and TP receptors (K_is = 30, 790, and 2,400 nM, respectively) and is without effect on other prostanoid receptors.^{1,2} In wild type mice, it mimics deletion of EP₄ by producing severe colitis, with epithelial loss, crypt damage, and inflammation, after treatment with 3% dextran sodium sulfate.¹ ONO-AE3-208 has also been used to implicate EP₄ signaling in immune and autoimmune responses, inflammation, and cancer.³⁻⁶

References

1. Kabashima, K., Saji, T., Murata, T., *et al.* The prostaglandin receptor EP4 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 EP4 antagonism inhibits breast cancer metastasis. *Cancer Res.* **66**(6), 2923-2927 (2006).
4. Kabashima, K., Sakata, D., Nagamachi, M., *et al.* Prostaglandin E₂-EP4 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 PGE₂-EP4 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 E₂ 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, 02/06/2024

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