PRODUCT INFORMATION



ONO-8711 (dicyclohexylamine salt)

Item No. 38432

Formal Name:	6-[(2R,3S)-3-[[[(4-chloro-2-methylphenyl) sulfonyl]amino]methyl]bicyclo[2.2.2] oct-2-yl]-5Z-hexenoic acid, dicyclohexylamine salt	CI
MF:	$C_{22}H_{30}CINO_4S \bullet C_{12}H_{23}N$	
FW:	621.3	•
Purity:	≥98%	
Supplied as:	A solid	но
Storage:	-20°C	Соон
Stability:	≥4 years	

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

Laboratory Procedures

ONO-8711 (dicyclohexylamine salt) is supplied as a solid. A stock solution may be made by dissolving the ONO-8711 (dicyclohexylamine salt) in the solvent of choice, which should be purged with an inert gas. ONO-8711 (dicyclohexylamine salt) is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

ONO-8711 is an antagonist of the prostaglandin E_2 (PGE₂; Item No. 14010) receptor subtype EP_1 (K_is = 0.6 and 1.7 nM for human and mouse receptors, respectively, expressed in CHO cells).¹ It is selective for EP₁ over EP₂, EP₃, and EP₄ (K_is = >1,000, 67, and >1,000 nM, respectively, for mouse receptors expressed in CHO cells). ONO-8711 inhibits PGE_2 -induced increases in intracellular calcium concentrations in cells expressing human or mouse EP₁ (EC₅₀s = 50 and 210 nM, respectively). Dietary administration of ONO-8711 (400 and 800 ppm) reduces the incidence of squamous cell carcinoma and severe dysplasia in a rat model of tongue carcinogenesis induced by 4-nitroquinoline 1-oxide.²

References

- 1. Watanabe, K., Kawamori, T., Nakatsugi, S., et al. Role of the prostaglandin E receptor subtype EP₁ in colon carcinogenesis. Cancer Res. 59(20), 5093-5096 (1999).
- 2. Makita, H., Mutoh, M., Maruyama, T., et al. A prostaglandin E₂ receptor subtype EP₁-selective antagonist, ONO-8711, suppresses 4-nitroquinoline 1-oxide-induced rat tongue carcinogenesis. Carcinogenesis 28(3), 677-684 (2007).

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

uyer 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, 03/23/2023

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