

PRODUCT INFORMATION



CJ-023423

Item No. 10010355

CAS Registry No.: 415903-37-6
Formal Name: N-[[2,4-(2-ethyl-4,6-dimethyl-1H-imidazo[4,5-c]pyridin-1-yl)phenylethylamino]carbonyl]-4-methyl-benzenesulfonamide

Synonyms: Grapiprant, RQ-00000007

MF: C₂₆H₂₉N₅O₃S

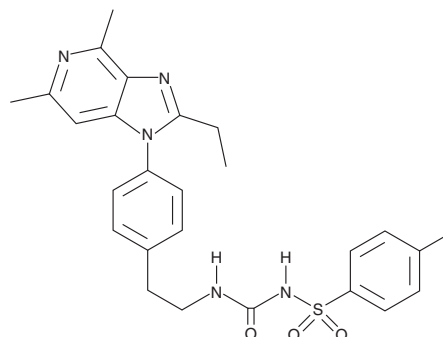
FW: 491.6

Purity: ≥98%

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

CJ-023423 is supplied as a crystalline solid. A stock solution may be made by dissolving the CJ-023423 in the solvent of choice, which should be purged with an inert gas. CJ-023423 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CJ-023423 in ethanol and DMSO is approximately 10 mg/ml and approximately 20 mg/ml in DMF.

CJ-023423 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CJ-023423 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CJ-023423 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Prostaglandin E₂ (PGE₂) activates four E prostanoid (EP) receptors, EP₁₋₄. EP₄ is a G_s protein-coupled receptor that, by elevating the second messenger cAMP, plays important roles in bone formation and resorption, cancer, and atherosclerosis.¹⁻³ CJ-023423 is a potent and selective antagonist of the EP₄ receptor (K_i = 13 and 20 nM for human and rat EP₄, respectively).^{4,5} It inhibits PGE₂-evoked elevation in intracellular cAMP in cells and, *in vivo*, reduces thermal hyperalgesia induced by intraplantar injection of PGE₂.⁵ CJ-023423 reduces acute and chronic inflammatory pain in different mouse models.⁵

References

1. 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).
2. Li, M., Thompson, D.D., and Paralkar, V.M. Prostaglandin E₂ receptors in bone formation. *Int. Orthop.* **31(6)**, 767-772 (2007).
3. Hawcroft, G., Ko, C.W.S., and Hull, M.A. Prostaglandin E₂-EP4 receptor signalling promotes tumorigenic behaviour of HT-29 human colorectal cancer cells. *Oncogene* **26(21)**, 3006-3019 (2007).
4. Nakao, K., Murase, A., Ohshiro, H., *et al.* CJ-023,423, a novel, potent and selective prostaglandin EP4 receptor antagonist with antihyperalgesic properties. *J. Pharmacol. Exp. Ther.* **322(2)**, 686-694 (2007).
5. 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).

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, 11/04/2022

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