

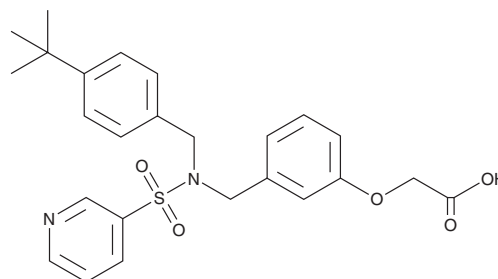
PRODUCT INFORMATION



Evatanepag

Item No. 42727

CAS Registry No.: 223488-57-1
Formal Name: 2-[3-[[[4-(1,1-dimethylethyl)phenyl]methyl](3-pyridinylsulfonyl)amino]methyl]phenoxy]-acetic acid
Synonym: CP 533,536
MF: C₂₅H₂₈N₂O₅S
FW: 468.6
Purity: ≥98%
Supplied as: A 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

Evatanepag is supplied as a solid. A stock solution may be made by dissolving the evatanepag in the solvent of choice, which should be purged with an inert gas. Evatanepag is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Evatanepag is an agonist of prostaglandin E₂ (PGE₂) receptor subtype EP₂.¹ It induces cAMP accumulation in cells expressing rat EP₂ (EC₅₀ = 0.3 nM). Evatanepag (3 mg/kg) increases total bone area, mineral content, and mineral density in rat tibia.

Reference

1. Cameron, K.O., Lefker, B.A., Ke, H.Z., *et al.* Discovery of CP-533536: An EP₂ receptor selective prostaglandin E₂ (PGE₂) agonist that induces local bone formation. *Bioorg. Med. Chem. Lett.* **19**(7), 2075-2078 (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, 02/04/2025

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