

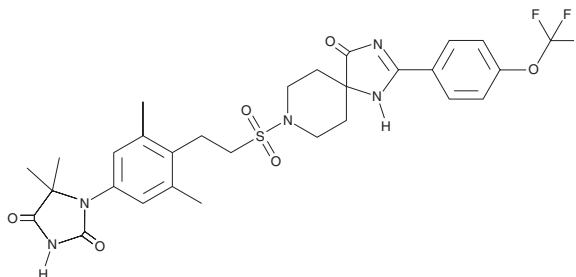
# PRODUCT INFORMATION



## PCO371

Item No. 39182

**CAS Registry No.:** 1613373-33-3  
**Formal Name:** 1-[3,5-dimethyl-4-[2-[[4-oxo-2-[4-(trifluoromethoxy)phenyl]-1,3,8-triazaspiro[4.5]dec-1-en-8-yl]sulfonyl]ethyl]phenyl]-5,5-dimethyl-2,4-imidazolidinedione  
**MF:** C<sub>29</sub>H<sub>32</sub>F<sub>3</sub>N<sub>5</sub>O<sub>6</sub>S  
**FW:** 635.7  
**Purity:** ≥95%  
**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

PCO371 is supplied as a solid. A stock solution may be made by dissolving the PCO371 in the solvent of choice, which should be purged with an inert gas. PCO371 is soluble in methanol.

### Description

PCO371 is a parathyroid hormone receptor type 1 (PTH1R) agonist.<sup>1</sup> It selectively induces cAMP production in COS-7 cells expressing human PTH1R (EC<sub>50</sub> = 2.4 μM) over COS-7 cells expressing human PTH2R (EC<sub>50</sub> = > 100 μM). PCO371 (1 and 3 μM) induces calcium release from isolated fetal rat long bones. It increases lumbar spine and proximal femur bone mineral density (BMD), lumbar spine bone strength, serum levels of osteocalcin, and bone formation in the lumbar spine without affecting serum calcium levels in an ovariectomized (OVX) rat model of osteopenia when administered at a dose of 10 mg/kg per day. PCO371 (9 mg/kg) increases serum, but not urinary, calcium levels in a hypocalcemic thyroparathyroidectomized (TPTX) rat model of hypothyroidism.

### References

1. Tamura, T., Noda, H., Joyashiki, E., *et al.* Identification of an orally active small-molecule PTHR1 agonist for the treatment of hypoparathyroidism. *Nat. Commun.* **7**, 13384 (2016).

#### 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, 02/22/2024

#### 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