

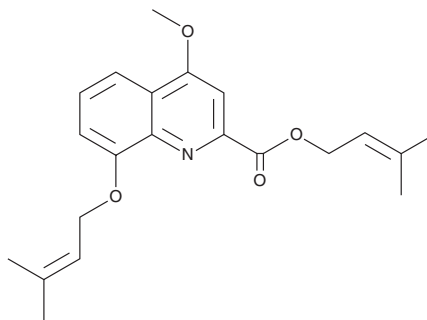
PRODUCT INFORMATION



Ppc-1

Item No. 42352

CAS Registry No.: 1245818-17-0
Formal Name: 4-methoxy-8-[(3-methyl-2-buten-1-yl)oxy]-2-quinolinecarboxylic acid, 3-methyl-2-buten-1-yl ester
MF: $C_{21}H_{25}NO_4$
FW: 355.4
Purity: $\geq 98\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/*P. pseudo-candidum*



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

Laboratory Procedures

Ppc-1 is supplied as a solid. A stock solution may be made by dissolving the Ppc-1 in the solvent of choice, which should be purged with an inert gas. Ppc-1 is soluble in DMSO.

Description

Ppc-1 is a prenylated quinolinecarboxylic acid originally isolated from *P. pseudo-candidum* that has diverse biological activities.¹⁻³ It reduces the proliferation of K562 chronic myelogenous leukemia (CML) cells and 3T3-L1 mouse fibroblasts when used at a concentration of 15 μM .¹ Ppc-1 (1 and 10 μM) inhibits the production of IL-2 induced by concanavalin A (Item No. 14951) in Jurkat T cells.² *In vivo*, Ppc-1 (0.8 mg/kg) reduces body weight gain and increases serum fatty acid levels in mice fed a commercial diet.³

References

1. Kikuchi, H., Nakamura, K., Kubohara, Y., *et al.* Novel prenylated and geranylated aromatic compounds isolated from *Polysphondylium* cellular slime molds. *Tetrahedron* **66**(32), 6000-6007 (2010).
2. Ogura, M., Kikuchi, H., Suzuki, T., *et al.* Prenylated quinolinecarboxylic acid derivative suppresses immune response through inhibition of PAK2. *Biochem Pharmacol.* **105**, 55-65 (2016).
3. Suzuki, T., Kikuchi, H., Ogura, M., *et al.* Weight loss by Ppc-1, a novel small molecule mitochondrial uncoupler derived from slime mold. *PLoS One* **10**(2), e0117088 (2015).

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, 07/01/2025

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