

PRODUCT INFORMATION

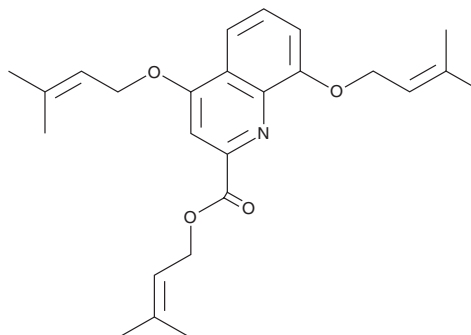


PQA-18

Item No. 42351

CAS Registry No.: 1604678-82-1
Formal Name: 4,8-bis[(3-methyl-2-buten-1-yl)oxy]-2-quinolinecarboxylic acid, 3-methyl-2-buten-1-yl ester
Synonym: Prenylated Quinolinecarboxylic Acid Compound-18

MF: C₂₅H₃₁NO₄
FW: 409.5
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

PQA-18 is supplied as a solid. A stock solution may be made by dissolving the PQA-18 in the solvent of choice, which should be purged with an inert gas. PQA-18 is soluble in DMSO.

Description

PQA-18 is an inhibitor of p21-activated kinase 2 (PAK2; IC₅₀ = 10 nM), an immunosuppressant, and a derivative of Ppc-1 (Item No. 42352).¹ It inhibits the production of IL-2 induced by concanavalin A (Item No. 14951) in Jurkat T cells (IC₅₀ = 400 nM). Topical application of PQA-18 (0.1%) decreases disease severity, serum IgE levels, and skin thickness in the NC/Nga mouse model of atopic dermatitis. It also reduces cutaneous nerve fiber density in NC/Nga mice when applied topically at 0.05%.² PQA-18 (4 mg/kg) increases graft survival rate in a rat model of small intestine transplant.³

References

1. 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).
2. Ogura, M., Endo, K., Suzuki, T., *et al.* Prenylated quinolinecarboxylic acid compound-18 prevents sensory nerve fiber outgrowth through inhibition of the interleukin-31 pathway. *PLoS One* **16(2)**, e0246630 (2021).
3. Kodama, T., Maeda, A., Lo, P.-C., *et al.* The effect of a novel immunosuppressive drug, a PAK-2 inhibitor, on macrophage differentiation/polarization in a rat small intestinal transplantation model. *Transpl. Immunol.* **57**, 101246 (2019).

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