

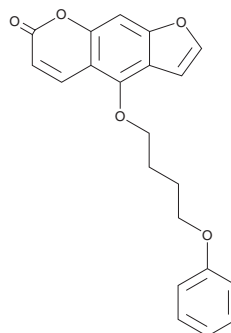
PRODUCT INFORMATION



PAP-1

Item No. 30105

CAS Registry No.: 870653-45-5
Formal Name: 4-(4-phenoxybutoxy)-7H-furo[3,2-g][1]benzopyran-7-one
Synonym: 5-(4-Phenoxybutoxy)psoralen
MF: C₂₁H₁₈O₅
FW: 350.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 249, 268, 309 nm
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

PAP-1 is supplied as a solid. A stock solution may be made by dissolving the PAP-1 in the solvent of choice, which should be purged with an inert gas. PAP-1 is soluble in the organic solvent DMSO at a concentration of approximately 9 mg/ml.

Description

PAP-1 is an inhibitor of K_v1.3 voltage-gated potassium channels (EC₅₀ = 2 nM at -80 mV in cells overexpressing the human channel).¹ It is selective for K_v1.3 over a panel of 22 ion channels, including potassium, sodium, calcium, and chloride channels (EC₅₀s = 45-15,000 nM for all). PAP-1 inhibits anti-CD3 antibody-induced proliferation of isolated human CCR7⁺ effector memory T cells. *In vivo*, PAP-1 (3 mg/kg) inhibits ovalbumin-induced delayed-type hypersensitivity in rats.

Reference

1. Schmitz, A., Sankaranarayanan, A., Azam, P., *et al.* Design of PAP-1, a selective small molecule Kv1.3 blocker, for the suppression of effector memory T cells in autoimmune diseases. *Mol. Pharmacol.* **68**(5), 1254-1270 (2005).

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, 10/27/2022

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