

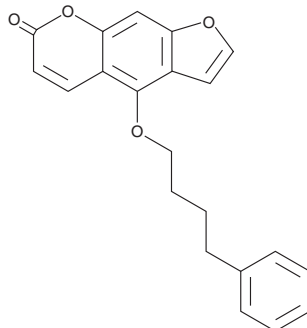
PRODUCT INFORMATION



Psora 4

Item No. 28389

CAS Registry No.: 724709-68-6
Formal Name: 4-(4-phenylbutoxy)-7H-furo[3,2-g]
[1]benzopyran-7-one
MF: C₂₁H₁₈O₄
FW: 334.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 251, 269, 310 nm
Supplied as: A crystalline 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

Psora 4 is supplied as a crystalline solid. A stock solution may be made by dissolving the psora 4 in the solvent of choice, which should be purged with an inert gas. Psora 4 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

Psora 4 is an inhibitor of voltage-gated potassium channels (K_v1.3; EC₅₀ = 3 nM).¹ It is selective for K_v1.3 over K_v1.1, K_v1.2, K_v1.4, K_v1.7, and K_v3.1 channels (EC₅₀s = 62, 49, 202, 100, and 1,500 nM, respectively) but does inhibit the K_v1.5 channel (EC₅₀ = 7.7 nM). Psora 4 selectively binds to the C-type inactivated state of the K_v1.3 channel and inhibits potassium current in a use-dependent manner. It has immunosuppressive activity, inhibiting proliferation of human and rat myelin-specific effector memory T cells *in vitro* (TEM cells; EC₅₀s = 25 and 60 nM, respectively).

Reference

1. Vennekamp, J., Wulff, H., Beeton, C., *et al.* K_v1.3-blocking 5-phenylalkoxy-psoralens: A new class of immunomodulators. *Mol. Pharmacol.* **65**(6), 1364-1374 (2004).

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, 11/10/2022

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