

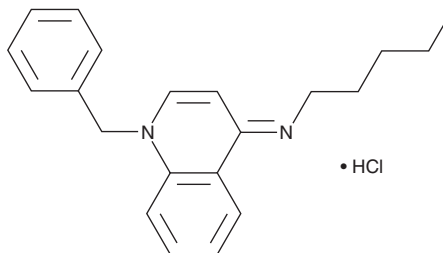
PRODUCT INFORMATION



CP 339,818 (hydrochloride)

Item No. 40766

CAS Registry No.: 478341-55-8
Formal Name: N-[1-(phenylmethyl)-4(1H)-quinolinyldiene]-1-pentanamine, monohydrochloride
MF: $C_{21}H_{24}N_2 \cdot HCl$
FW: 340.9
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

CP 339,818 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the CP 339,818 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. CP 339,818 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

CP 339,818 is an inhibitor of voltage-gated potassium channel $K_v1.3$ and $K_v1.4$ ($IC_{50}s = 0.23$ and $0.3 \mu M$, respectively).¹ It is selective for these channels over $K_v1.1$, $K_v1.2$, $K_v1.5$, and $K_v1.6$ channels ($IC_{50}s = 62$, 14 , 19 , and $20 \mu M$, respectively), as well as $K_v3.1$, $K_v3.2$, $K_v3.4$, and $K_v4.2$ channels ($IC_{50}s = 17$, 10 , 36 , and $>10 \mu M$, respectively). CP 339,818 inhibits thymidine incorporation, a marker of T cell activation, induced by an anti-CD3 antibody or a combination of ionomycin (Item Nos. 10004974 | 11932) and phorbol 12-myristate 13-acetate (PMA; Item No. 10008014) in isolated human peripheral blood T cells ($IC_{50}s = 4.7$ and $5.2 \mu M$, respectively).

Reference

1. Nguyen, A., Kath, J.C., Hanson, D.C., *et al.* Novel nonpeptide agents potently block the C-type inactivated conformation of $K_v1.3$ and suppress T cell activation. *Mol. Pharmacol.* **50**(6), 1672-1679 (1996).

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, 04/25/2024

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