

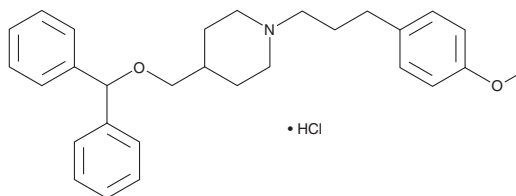
PRODUCT INFORMATION



UK 78282

Item No. 45311

CAS Registry No.: 136647-02-4
Formal Name: 4-[[diphenylmethoxy)methyl]-1-[3-(4-methoxyphenyl)propyl]-piperidine, monohydrochloride
MF: C₂₉H₃₅NO₂ • HCl
FW: 466.1
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

UK 78282 is supplied as a solid. A stock solution may be made by dissolving the UK 78282 in the solvent of choice, which should be purged with an inert gas. UK 78282 is soluble (≥10 mg/ml) in DMSO.

Description

UK 78282 is an inhibitor of voltage-gated potassium channel 1.3 (K_v1.3) and K_v1.4 (IC₅₀s = 0.28 and 0.17 μM, respectively).¹ It is selective for K_v1.3 and K_v1.4 over K_v1.1-1.2 and K_v1.5-1.6 (IC₅₀s = 2.9-70 μM), K_v3.1, -3.2, and -3.4 (IC₅₀s = 11-127 μM), K_v4.2 (IC₅₀ = 10 μM), and intermediate conductance calcium-activated potassium channel (K_{Ca}3.1/IK; IC₅₀ = >30 μM). UK 78282 inhibits phytohemagglutinin-induced proliferation and IL-2 production in T cells (IC₅₀s = 2 and 2.9 μM, respectively).

Reference

1. Hanson, D.C., Nguyen, A., Mather, R.J., *et al.* UK-78,282, a novel piperidine compound that potently blocks the Kv1.3 voltage-gated potassium channel and inhibits human T cell activation. *Br. J. Pharmacol.* **126**(8), 1707-1716 (1999).

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, 03/27/2026

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