

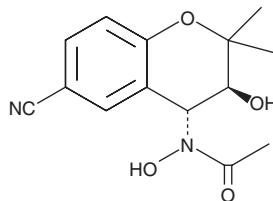
PRODUCT INFORMATION



Y-26763

Item No. 21459

CAS Registry No.: 127408-31-5
Formal Name: N-[(3S,4R)-6-cyano-3,4-dihydro-3-hydroxy-2,2-dimethyl-2H-1-benzopyran-4-yl]-N-hydroxy-acetamide
MF: C₁₄H₁₆N₂O₄
FW: 276.3
Purity: ≥98%
UV/Vis.: λ_{max}: 253, 290 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

Y-26763 is supplied as a crystalline solid. A stock solution may be made by dissolving the Y-26763 in the solvent of choice, which should be purged with an inert gas. Y-26763 is soluble in organic solvents such as ethanol and DMSO. The solubility of Y-26763 in these solvents is approximately 100 mM.

Description

Y-26763 is an activator of ATP-sensitive potassium (K_{ATP}) channels.¹ It increases the frequency of K_{ATP} channel openings in the presence of ATP and activates K_{ATP} channel currents in isolated human β-cells when used at a concentration of 200 μM. It also inhibits increases in intracellular calcium induced by histamine or norepinephrine (Item No. 16673) in isolated rabbit mesenteric artery smooth muscle in a concentration-dependent manner.² Y-26763 induces relaxation in isolated precontracted rat aortic rings (IC₅₀ = 0.027 μM).³ It decreases mean blood pressure in spontaneously hypertensive rats when administered intravenously at a dose of 0.03 mg/kg, an effect that can be reversed by glibenclamide (Item No. 15009).

References

1. Cosgrove, K.E., Straub, S.G., Barnes, P.D., *et al.* Y-26763: ATP-sensitive K⁺ channel activation and the inhibition of insulin release from human pancreatic beta-cells. *Eur. J. Pharmacol.* **486**(2), 133-139 (2004).
2. Watanabe, Y., Suzuki, A., Suzuki, H., *et al.* Effect of membrane hyperpolarization induced by a K⁺ channel opener on histamine-induced Ca²⁺ mobilization in rabbit arterial smooth muscle. *Br. J. Pharmacol.* **117**(6), 1302-1308 (1996).
3. Nakajima, T., Shinohara, T., Yaoka, O., *et al.* Y-27152, a long-acting K⁺ channel opener with less tachycardia: Antihypertensive effects in hypertensive rats and dogs in conscious state. *J. Pharmacol. Exp. Ther.* **261**(2), 730-736 (1992).

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, 11/07/2022

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