

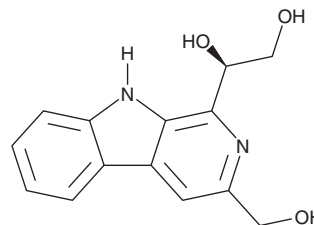
PRODUCT INFORMATION



Pyridindolol

Item No. 32914

CAS Registry No.: 55812-46-9
Formal Name: α¹R-(hydroxymethyl)-9H-pyrindo[3,4-b]indole-1,3-dimethanol
Synonym: NSC 266530
MF: C₁₄H₁₄N₂O₃
FW: 258.3
Purity: ≥90%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces alboverticillatus* MD401-C5



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

Laboratory Procedures

Pyridindolol is supplied as a crystalline solid. A stock solution may be made by dissolving the pyridindolol in the solvent of choice. Pyridindolol is soluble in organic solvents such as ethanol, methanol, 1-butanol, DMSO, dimethyl formamide, and 0.5M HCl, which should be purged with an inert gas.

Pyridindolol is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Pyridindolol is a bacterial metabolite that has been found in *S. alboverticillatus*.¹ It inhibits neutral β-galactosidase by 50% under acidic, but not neutral, conditions when used at a concentration of 2 μg/ml. It is selective for β-galactosidase isolated from bovine liver over β-galactosidases isolated from human, bovine, pig, and rat tissues and sialidases isolated from *C. perfringens*, *Streptomyces*, and the H3N2 strain of influenza virus (IC₅₀s = >250 μg/ml for all).^{1,2}

References

1. Aoyagi, T., Kumagai, M., Hazato, T., *et al.* Pyridindolol, a new β-galactosidase inhibitor produced by actinomycetes. *J. Antibiot. (Tokyo)* **28(7)**, 555-557 (1975).
2. Kumagai, M., Aoyagi, T., and Umezawa, H. Inhibitory activity of pyridindolol on β-galactosidase. *J. Antibiot. (Tokyo)* **29(7)**, 696-703 (1976).

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