

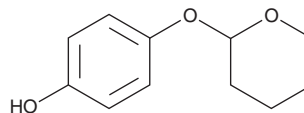
PRODUCT INFORMATION



(S)-(-)-Deoxyarbutin

Item No. 9001104

Formal Name: 4-[(tetrahydro-2H-pyran-2-yl)oxy]-phenol
MF: C₁₁H₁₄O₃
FW: 194.2
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

(S)-(-)-Deoxyarbutin is supplied as a solid. A stock solution may be made by dissolving the (S)-(-)-deoxyarbutin in the solvent of choice, which should be purged with an inert gas. (S)-(-)-Deoxyarbutin is soluble in methanol.

Description

(S)-(-)-Deoxyarbutin is a tyrosinase inhibitor and an isomer of deoxyarbutin (Item No. 21077).¹ It inhibits tyrosinase activity when used at a concentration of 74 μM.

Reference

1. Delong, M.A., Kvalnes, K.L., and Boissy, R. Chiral compounds, compositions, products and methods employing same. **WO2013103873A1** (2013).

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, 10/11/2023

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