

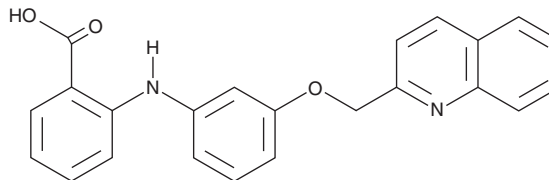
PRODUCT INFORMATION



SR 2640

Item No. 42267

CAS Registry No.: 105350-26-3
Formal Name: 2-[[3-(2-quinolinylmethoxy)phenyl]amino]-benzoic acid
MF: C₂₃H₁₈N₂O₃
FW: 370.4
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

SR 2640 is supplied as a solid. A stock solution may be made by dissolving the SR 2640 in the solvent of choice, which should be purged with an inert gas. SR 2640 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

SR 2640 is an antagonist of leukotriene D₄ (LTD₄; Item No. 20310) and LTE₄ (Item No. 20410).¹ It inhibits LTD₄- or LTE₄-induced contraction of isolated guinea pig ileum (IC₅₀s = ~0.003 μM for both), but not LTC₄-induced contractions. SR 2640 (1 mg/kg) inhibits ovalbumin-induced bronchoconstriction in anesthetized guinea pigs.

Reference

1. Ahnfelt-Rønne, I., Kirstein, D., and Kaergaard-Nielsen, C. A novel leukotriene D₄/E₄ antagonist, SR2640 (2-[[3-(2-quinolinylmethoxy)phenyl]amino]benzoic acid). *Eur. J. Pharmacol.* **155(1-2)**, 117-128 (1988).

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, 04/24/2025

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