

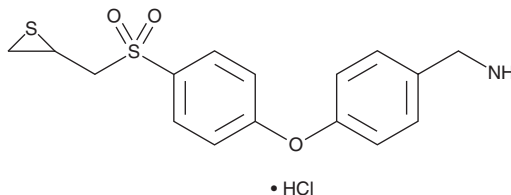
PRODUCT INFORMATION



ND-336

Item No. 19207

CAS Registry No.: 1807453-83-3
Formal Name: 4-[4-[(2-thiiranylmethyl)sulfonyl]phenoxy]-benzenemethanamine, monohydrochloride
MF: C₁₆H₁₇NO₃S₂ • HCl
FW: 371.9
Purity: ≥95%
UV/Vis.: λ_{max}: 245, 333 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

ND-336 is supplied as a crystalline solid. A stock solution may be made by dissolving the ND-336 in the solvent of choice, which should be purged with an inert gas. ND-336 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ND-336 in these solvents is approximately 30 and 5 mg/ml, respectively.

ND-336 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ND-336 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ND-336 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ND-336 is a selective inhibitor of the gelatinases matrix metalloproteinase-2 (MMP-2) and MMP-9, as well as MMP-14 (K_is = 85, 150, and 120 nM, respectively).¹ It has minimal effect on other MMPs. ND-336 accelerates diabetic wound healing in mice by lowering inflammation and by enhancing angiogenesis and re-epithelialization of the wound.¹

Reference

1. Gao, M., Nguyen, T.T., Suckow, M.A., *et al.* Acceleration of diabetic wound healing using a novel protease-anti-protease combination therapy. *Proc. Natl. Acad. Sci. USA* **112**(49) (2015).

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, 02/06/2025

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