

PRODUCT INFORMATION

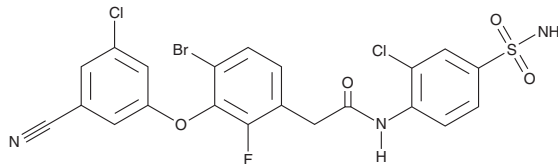


Ro 0335

Item No. 36124

CAS Registry No.: 867365-76-2
Formal Name: N-[4-(aminosulfonyl)-2-chlorophenyl]-4-bromo-3-(3-chloro-5-cyanophenoxy)-2-fluorobenzeneacetamide

Synonym: VM 1500A
MF: C₂₁H₁₃BrCl₂FN₃O₄S
FW: 573.2
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

Ro 0335 is supplied as a solid. A stock solution may be made by dissolving the Ro 0335 in the solvent of choice, which should be purged with an inert gas. Ro 0335 is soluble in DMSO.

Description

Ro 0335 is a non-nucleoside reverse transcriptase inhibitor (NNRTI) and an active metabolite of the antiviral prodrug el sulfavirine.¹ It inhibits HIV-1 reverse transcriptase in a cell-free assay and the replication of HIV-1 in MT-4 cells (IC₅₀s = 8.1 and 1.1 nM, respectively). Ro 0335 also selectively inhibits human carbonic anhydrase VII (CAVII; K_i = 9.7 nM) over human CAI and CAIII-VI (K_is = >125 nM for all) but also inhibits human CAII, CAIX, CAXII, CAXIII, and CAXIV (K_is = 35.5, 72.1, 62.3, 52.2, and 15.4 nM, respectively).²

References

- Javanbakht, H., Ptak, R., G., Chow, E., *et al.* *In vitro* resistance development for RO-0335, a novel diphenylether nonnucleoside reverse transcriptase inhibitor. *Antiviral Res.* **86(2)**, 212-219 (2010).
- Supuran, C.T., Nocentini, A., Yakubova, E., *et al.* Biochemical profiling of anti-HIV prodrug El sulfavirine (Elpida®) and its active form VM1500A against a panel of twelve human carbonic anhydrase isoforms. *J. Enzyme Inhib. Med. Chem.* **36(1)**, 1056-1060 (2021).

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, 12/19/2022

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