

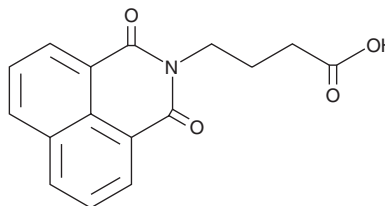
PRODUCT INFORMATION



Virstatin

Item No. 21176

CAS Registry No.: 88909-96-0
Formal Name: 1,3-dioxo-1H-benz[de]
isoquinoline-2(3H)-butanoic acid
MF: $C_{16}H_{13}NO_4$
FW: 283.3
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 214, 234 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Virstatin is supplied as a crystalline solid. A stock solution may be made by dissolving the virstatin in the solvent of choice, which should be purged with an inert gas. Virstatin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of virstatin in these solvents is approximately 33 mg/ml.

Virstatin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, virstatin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Virstatin has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Virstatin is an inhibitor of the ToxT transcriptional regulator of *V. cholerae*, which regulates transcription of virulence factors that enable intestinal colonization.¹ It inhibits ToxT dimerization and decreases expression of cholera toxin (CT) and toxin coregulated pilus (TCP) when used at a concentration of 50 μM but virstatin does not inhibit growth of *V. cholerae* at this concentration (MBCs = 600 and 1,200 μM for O395 and C6706 strains, respectively).^{1,2} Virstatin administration protects infant mice (5 to 6 days old) from intestinal colonization by ToxT-dependent *V. cholerae* but not from strains that colonize *via* non-ToxT-dependent mechanisms. Virstatin (100 μM) inhibits biofilm formation by *A. baumannii* by 38% under static conditions, which is at a lower concentration than that which inhibits growth (MIC = 1.6 mM).³ It decreases the motility of 60% of 30 mobile *A. baumannii* strains. Virstatin also binds to the accessory cholera enterotoxin (Ace) from *V. cholerae* ($K_a = 9 \times 10^4 M^{-1}$; $K_d = 11 \mu M$).⁴

References

1. Hung, D.T., Shakhnovich, E.A., Pierson, E.E., *et al.* Small-molecule inhibitor of *Vibrio cholerae* virulence and intestinal colonization. *Science* **310**(5748), 670-674 (2005).
2. Shakhnovich, E.A., Hung, D.T., Pierson, E., *et al.* Virstatin inhibits dimerization of the transcriptional activator ToxT. *Proc. Nat. Acad. Sci. USA* **104**(7), 2372-2377 (2007).
3. Chabane, Y.N., Mlouka, M.B., Alexandre, S., *et al.* Virstatin inhibits biofilm formation and motility of *Acinetobacter baumannii*. *BMC Microbiol.* **14**(62) (2014).
4. Chatterjee, T.K., Mukherjee, D., Dey, S., *et al.* Accessory cholera enterotoxin, Ace, from *Vibrio cholerae*: Structure, unfolding, and virstatin binding. *Biochemistry* **50**(14), 2962-2972 (2011).

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, 11/14/2022

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