

PRODUCT INFORMATION

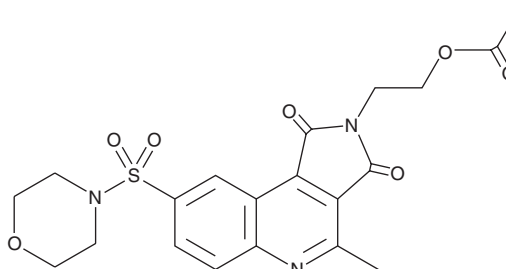


Caspase-3 Inhibitor VII

Item No. 13306

CAS Registry No.: 745046-84-8
Formal Name: 2-[2-(acetyloxy)ethyl]-4-methyl-8-(4-morpholinylsulfonyl)-1H-pyrrolo[3,4-c]quinoline-1,3(2H)-dione

Synonym: Ivachtin
MF: C₂₀H₂₁N₃O₇S
FW: 447.5
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 258, 349 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

Caspase-3 inhibitor VII is supplied as a crystalline solid. A stock solution may be made by dissolving the caspase-3 inhibitor VII in the solvent of choice, which should be purged with an inert gas. Caspase-3 inhibitor VII is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 2.5 mg/ml. Caspase-3 inhibitor VII is slightly soluble in DMSO.

Caspase-3 inhibitor VII is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, caspase-3 inhibitor VII should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Caspase-3 inhibitor VII has a solubility of approximately 0.03 mg/ml in a 1:30 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Caspase-3 inhibitor VII is a 4-methylated pyrroloquinoline that reversibly inhibits caspase-3 (IC₅₀ = 23 nM).¹ It displays anti-apoptotic activity in Jurkat T cells treated with staurosporine (Item No. 81590).¹

Reference

1. Kravchenko, D.V., Kuzovkova, Y.A., Kysil, V.M., *et al.* Synthesis and structure-activity relationship of 4-substituted 2-(2-acetyloxyethyl)-8-(morpholine-4-sulfonyl)pyrrolo[3,4-c]quinoline-1,3-diones as potent caspase-3 inhibitors. *J. Med. Chem.* **48**(11), 3680-3683 (2005).

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, 10/25/2022

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