

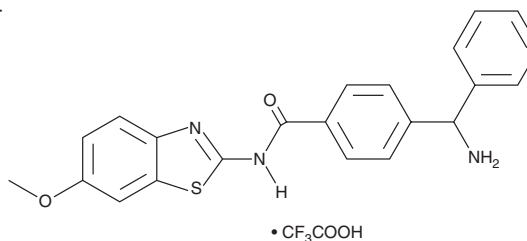
PRODUCT INFORMATION



SW203668 (trifluoroacetate salt)

Item No. 19379

CAS Registry No.: 2117405-48-6
Formal Name: 4-(aminophenylmethyl)-N-(6-methoxy-2-benzothiazolyl)-benzamide, trifluoroacetate salt
MF: C₂₂H₁₉N₃O₂S • CF₃COOH
FW: 503.5
Purity: ≥98%
UV/Vis.: λ_{max}: 316 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

SW203668 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the SW203668 (trifluoroacetate salt) in the solvent of choice. SW203668 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of SW203668 (trifluoroacetate salt) in these solvents is approximately 20 mg/ml.

SW203668 (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SW203668 (trifluoroacetate salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SW203668 (trifluoroacetate salt) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SW203668 is an irreversible inhibitor of stearoyl-CoA desaturase (IC₅₀ = 54 nM).¹ It is selectively cytotoxic to H2122, H460, HCC44, and HCC95 cell lines that express cytochrome P450 (CYP) isoform CYP4F11 over eight other cancer cell lines that lack CYP4F11 *in vitro* (IC₅₀s = 22-116 and >10,000 nM, respectively) and ectopic expression of CYP4F11 in SW203668-insensitive H1155 cells results in sensitization to SW203668. *In vivo*, SW203668 reduces tumor growth rate without reducing sebocyte production in the H2122 wild-type and nonobese diabetic severe combined immunodeficiency (NOD-SCID) mouse xenograft models when administered at doses of 20 and 6 mg/kg, respectively.

Reference

1. Theodoropoulos, P.C., Gonzales, S.S., Winterton, S.E., *et al.* Discovery of tumor-specific irreversible inhibitors of stearoyl CoA desaturase. *Nat. Chem. Biol.* **12**(4), 218-225 (2016).

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, 07/08/2019

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