

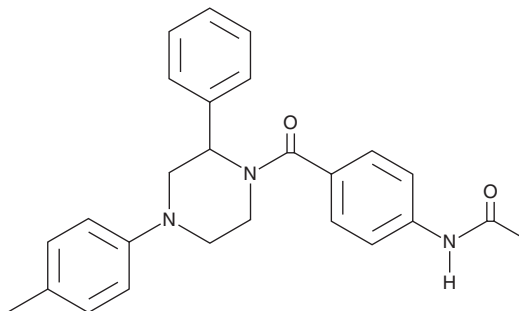
PRODUCT INFORMATION



SBC-110736

Item No. 23103

CAS Registry No.: 1629166-02-4
Formal Name: N-[4-[[4-(4-methylphenyl)-2-phenyl-1-piperazinyl]carbonyl]phenyl]-acetamide
MF: C₂₆H₂₇N₃O₂
FW: 413.5
Purity: ≥98%
UV/Vis.: λ_{max}: 253 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

SBC-110736 is supplied as a crystalline solid. A stock solution may be made by dissolving the SBC-110736 in the solvent of choice, which should be purged with an inert gas. SBC-110736 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SBC-110736 in these solvents is approximately 10 and 15 mg/ml, respectively.

SBC-110736 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SBC-110736 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SBC-110736 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SBC-110736 is an inhibitor of proprotein convertase subtilisin kexin type 9 (PCSK9).¹ It increases LDL receptor (LDLR) levels in HepG2 cells expressing PCSK9 when used at a concentration of 400 nM. SBC-110736 (8 mg/kg per day) reduces plasma total cholesterol levels by a mean of 38% in mice fed a high-fat diet. It also reduces plasma LDL cholesterol levels in mice fed a high-fat diet and exhibits an additive effect when administered in combination with atorvastatin (Item No. 10493).

Reference

1. Abdel-Meguid, S.S., Abou-Gharbia, M., Blass, B., *et al.* Anti-pcsk9 compounds and methods for the treatment and/or prevention of cardiovascular diseases. *Shifa Biomedical Corp.* **US20150376139A1** (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, 10/04/2022

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