

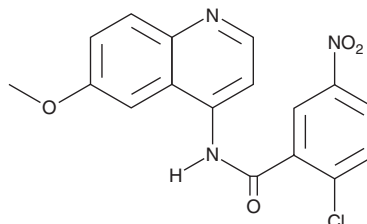
PRODUCT INFORMATION



SR 16832

Item No. 27632

CAS Registry No.: 2088135-12-8
Formal Name: 2-chloro-N-(6-methoxy-4-quinoliny)-5-nitro-benzamide
MF: C₁₇H₁₂ClN₃O₄
FW: 357.7
Purity: ≥98%
UV/Vis.: λ_{max}: 233 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

SR 16832 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 16832 in the solvent of choice, which should be purged with an inert gas. SR 16832 is soluble in the organic solvent DMSO.

Description

SR 16832 is a dual-site covalent antagonist of peroxisome proliferator-activated receptor γ (PPAR γ).¹ It inhibits MRL20-induced allosteric activation of PPAR γ in a reporter assay using HEK293T cells when used at a concentration of 5 μ M. SR 16832 also reduces basal activity of PPAR γ and inhibits binding of docosahexaenoic acid (DHA; Item No. 90310) to PPAR γ in a time-resolved FRET (TR-FRET) assay.

Reference

1. Brust, R., Lin, H., Fuhrmann, J., *et al.* Modification of the orthosteric PPAR γ covalent antagonist scaffold yields an improved dual-site allosteric inhibitor. *ACS Chem. Biol.* **12**(4), 969-978 (2017).

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