

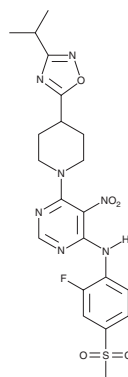
PRODUCT INFORMATION



AR231453

Item No. 34898

CAS Registry No.: 733750-99-7
Formal Name: N-[2-fluoro-4-(methylsulfonyl)phenyl]-6-[4-[3-(1-methylethyl)-1,2,4-oxadiazol-5-yl]-1-piperidinyl]-5-nitro-4-pyrimidinamine
MF: C₂₁H₂₄FN₇O₅S
FW: 505.5
Purity: ≥95%
UV/Vis.: λ_{max}: 285, 358 nm
Supplied as: A 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

AR231453 is supplied as a solid. A stock solution may be made by dissolving the AR231453 in the solvent of choice, which should be purged with an inert gas. AR231453 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AR231453 in these solvents is approximately 3 and 25 mg/ml, respectively.

Description

AR231453 is an agonist of GPR119.¹ It increases cAMP accumulation in HIT-T15 β cells that endogenously express high levels of GPR119 (EC₅₀ = 4.7 nM). AR231453 increases insulin release from HIT-T15 cells (EC₅₀ = 3.5 nM) and increases glucose-induced insulin release in isolated rat and mouse islets. It reduces plasma glucose levels in wild-type, KK/Ay, and db/db diabetic mice when administered at a dose of 20 mg/kg.

Reference

1. Chu, Z.L., Jones, R.M., He, H., *et al.* A role for β-cell-expressed GPR119 in glycemic control by enhancing glucose-dependent insulin release. *Endocrinology* **148**(6), 2601-2609 (2007).

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, 02/17/2026

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