

AR231453

GPR119 activator

Synthetic agonist of the orphan GPR119. Although endogenous lipid amides such as oleoyl ethanolamide have been reported to be agonists at GPR119, AR231453 is the first reported high potency agonist ($EC_{50} = 0.68\text{nM}$). The compound displayed *in vivo* activity in rodents and was active in an oral glucose tolerance test in mice following oral administration. AR231453 stimulated the release of GLP-1 via GPR119 receptors in mouse models. It was shown to be a highly selective ligand with no off-target activity across a 76 receptor and enzyme profiling panel (CEREP) and no activity against 140 known and orphan GPCRs in melanophore dispersion assays.

Citations: 5

[View Online »](#)

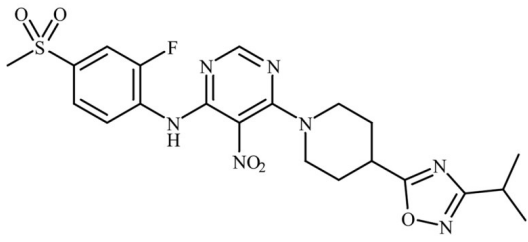
Ordering Information

[Order Online »](#)

BML-GP100-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2-Fluoro-4-methanesulfonyl-phenyl)-{6-[4-(3-isopropyl-[1,2,4]oxadiazol-5-yl)-piperidin-1-yl]-5-nitro-pyrimidin-4-yl]-amine
Appearance	Yellow solid.
CAS	733750-99-7
Couple Target	GPR119
Couple Type	Activator, Ligand
Formula	$C_{21}H_{24}FN_7O_5S$
Identity	Determined by NMR.
MW	505.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com