

AdipoRon

Novel and orally active adiponectin receptor agonist for AdipoR1 and AdipoR2.

AdipoR agonist that binds to both AdipoR1 and AdipoR2.

Produced by adipocytes, adiponectin has circulating levels that are inversely related to the degree of adiposity. Adiponectin concentrations correlate negatively with glucose, insulin, triglyceride concentrations, liver fat content and body mass index, while correlating positively with high-density lipoprotein-cholesterol levels, hepatic insulin sensitivity and insulin-stimulated glucose disposal. Adiponectin controls many metabolic processes and is known to contribute to metabolic pathologies including diabetes, atherosclerosis, and obesity. It has been shown to increase insulin sensitivity and decrease plasma glucose by increasing tissue fat oxidation. Adiponectin receptor activation is a promising potential therapy for diabetes and other obesity-associated diseases.

AdipoRon is a small-molecule adiponectin mimetic. At low micromolar concentrations, it binds to the adiponectin receptors AdipoR1 and AdipoR2. In addition, AdipoRon is orally administered and readily absorbed and delivered to target tissues, making it potentially clinically relevant. AdipoRon has been shown to induce insulin sensitivity, inhibit lipid synthesis and increase glucose metabolism.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-CHM101-0005	5mg
ENZ-CHM101-0010	10mg

Manuals, SDS & CofA

[View Online »](#)

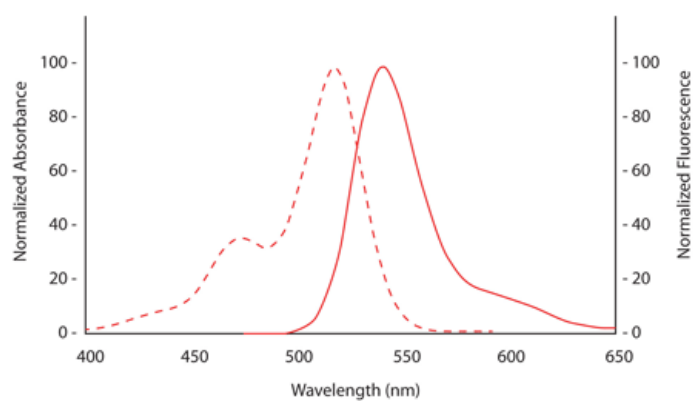
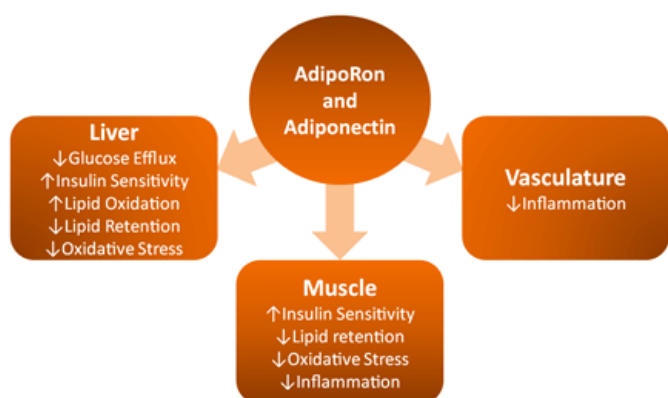
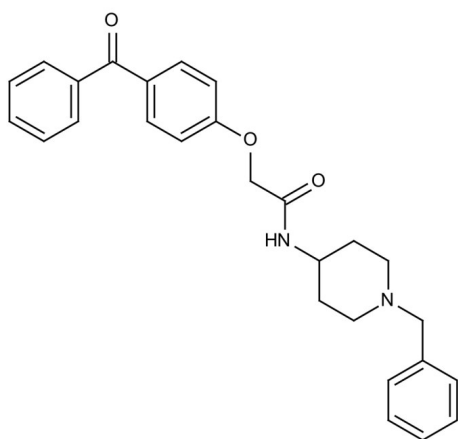


Fig 1. AdipoRon and Adiponectin have similar downstream effects.



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, up to 3 years at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(4-benzoylphenoxy)-N-(1-benzylpiperidin-4-yl)acetamide
CAS	924416-43-3
Couple Target	Adiponectin receptor
Couple Type	Activator, Ligand
Formula	$C_{27}H_{28}N_2O_3$
MW	428.52
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (86mg/ml) or ethanol (50mg/ml). Slightly soluble in water (<1mg/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-be@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