

SLU-PP-332

pan-ERR agonist/Exercise mimetic

Pan-agonist of the estrogen receptor-related receptors with highest potency for ERR α . SLU-PP-332 increased mitochondrial function and cellular respiration in a skeletal muscle cell line. It also increased fast oxidative skeletal muscle fibers (Type IIa) and exercise endurance in mice in an ERR α -specific manner. SLU-PP-332 activates an acute aerobic exercise genetic response via inducing the expression of DDIT4.

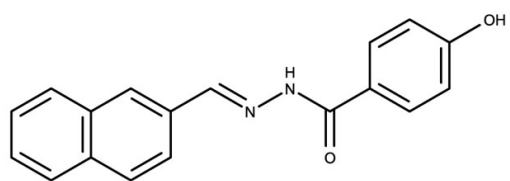
Ordering Information

[Order Online »](#)

ENZ-CHM474-0050	50mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (E)-4-Hydroxy-N'-(naphthalen-2-ylmethylene)benzohydrazide

Appearance White solid.

CAS 303760-60-3

Couple Target Estrogen-related receptor alpha

Couple Type Activator

Formula $C_{18}H_{14}N_2O_2$

Identity Determined by NMR.

MW 290.32

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (75 mg/ml).

Last modified: February 27, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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