

A-769662

AMPK activator

A-769662 is a potent activator of AMP-activated protein kinase (AMPK), with an EC50 of 0.8 μ M in vitro. It specifically targets AMPK heterotrimers containing the β 1 subunit.

Key features and applications include:

- **High Selectivity:** Selectively activates AMPK β 1 subunit-containing heterotrimers.
- **Potency:** Demonstrates strong activation with an EC50 of 0.8 μ M.
- **Stimulate Fatty Acid Oxidation:** Promotes fatty acid oxidation in various cell types.
- **Inhibit Adipocyte Differentiation:** Reduces differentiation of preadipocytes into adipocytes.
- **Enhance Glucose Uptake:** Increases glucose uptake in skeletal muscle cells.
- **Promote Endothelial Cell Survival:** Supports endothelial cell survival during metabolic stress.

Relevant disease states include:

- **Type 2 Diabetes:** Improves glucose homeostasis and reduces plasma glucose levels.
- **Obesity:** Inhibits adipogenesis, making it useful in obesity research.
- **Metabolic Syndrome:** Addresses key components of metabolic syndrome by activating AMPK.
- **Cardiovascular Diseases:** Enhances endothelial cell survival, which is beneficial in cardiovascular research.

- GMP format available

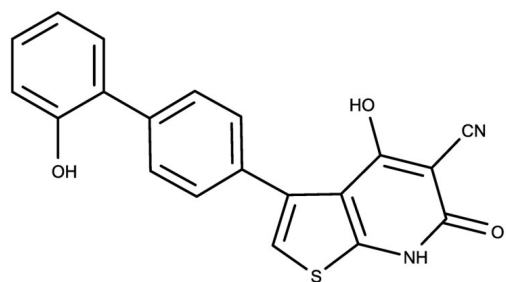
Ordering Information

[Order Online »](#)

ENZ-CHM360-0025	25mg
-----------------	------

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO or ethanol may be stored at -20°C for up to 3 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-Hydroxy-3-(2'-hydroxy-[1,1'-biphenyl]-4-yl)-6-oxo-6,7-dihydrothieno[2,3-b]pyridine-5-carbonitrile
Appearance	Beige solid.
CAS	844499-71-4
Couple Target	AMPK
Couple Type	Activator
Formula	$C_{20}H_{12}N_2O_3S$
Identity	Determined by NMR.
MW	360.39
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 30 mg/ml) or in ethanol (up to 5 mg/ml).



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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@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)