

AA147

ATF6 activator

AA147 is an endoplasmic reticulum (ER) proteostasis regulator. It activates the ATF6-dependent ER stress response and the NRF2 oxidative stress response, providing protection against oxidative damage in neuronal cells.

Key features and applications include:

- **ER Proteostasis Regulation:** Activates ATF6 and NRF2 pathways.
- **Oxidative Stress Protection:** Reduces reactive oxygen species (ROS) and protects against glutamate-induced cytotoxicity.
- **Broad-Spectrum Antiviral Activity:** Inhibits dengue and Zika virus infections in an ATF6-independent manner.

Relevant disease states include:

- **Neurodegenerative Diseases:** Potential therapeutic for conditions like ischemic stroke and Alzheimer's disease.
- **Multiple Sclerosis:** Alleviates symptoms in experimental autoimmune encephalitis (EAE) models by reducing ER stress and oligodendrocyte loss.
- **Viral Infections:** Effective against dengue and Zika viruses.

Ordering Information

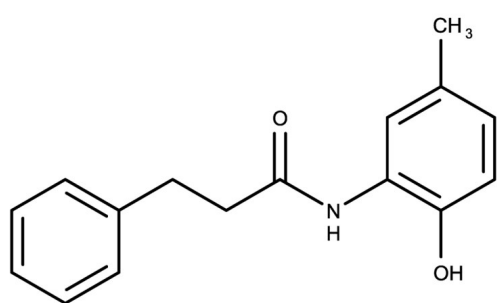
[Order Online »](#)

ENZ-CHM350-0050	50mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- GMP format available



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	N-(2-Hydroxy-5-methylphenyl)-3-phenylpropanamide, Compound 147
Appearance	Beige to pale orange solid.
CAS	393121-74-9
Couple Target	ATF
Couple Type	Activator
Formula	$C_{16}H_{17}NO_2$
Identity	Determined by NMR.
MW	255.32
Purity	≥98% HPLC
Solubility	Soluble in DMSO (>25 mg/ml) and ethanol (>25 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-
be@enzolifesciences.com](mailto:info-be@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)