

ACP

Probe for oxidative stress detection

Cell permeable probe for the detection of intracellular oxidative stress. ACP can easily be deprotected with esterases to cyclic hydroxylamines, which are oxidized by ROS leading to ESR-detectable nitroxide radicals.

Citations: 4

[View Online »](#)

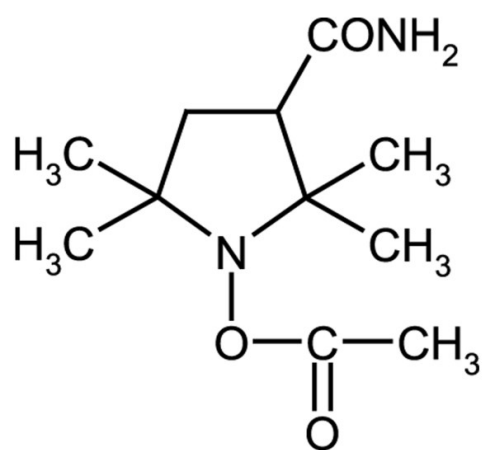
Ordering Information

[Order Online »](#)

ALX-430-118-M050	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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-Acetoxy-3-carbamoyl-2,2,5,5-tetramethylpyrrolidine

Appearance White to off-white crystalline powder.

Formula $C_{11}H_{20}N_2O_3$

MW 228.3

Purity ≤0.05% paramagnetic impurities (by EPR)

Solubility Soluble in water, methanol, 100% ethanol or DMSO.

Last modified: May 29, 2024



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)