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

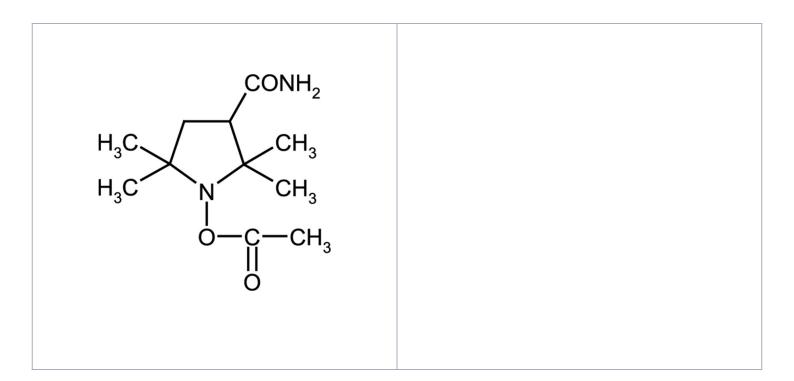
Order Online »

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

