Pyocyanin

ROS inducer

Cytotoxic pigment secreted by *Pseudomonas aeruginosa*, which frequently infects the lungs of immuno-suppressed patients, as well as those with cystic fibrosis. Pyocyanin toxicity results from its ability to undergo reduction by NAD(P)H and subsequent generation of superoxide and $\rm H_2O_2$ directly in the lungs.

Citations: 6

View Online »

Ordering Information

Order Online »

ENZ-53001-C001

1µmol

Manuals, SDS & CofA

View Online »

Figure: Pyocyanin is reduced by NAD(P)H and generates superoxide and $\rm H_2O_2$.

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -80°C.

Handling Avoid freeze/thaw cycles.

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Sanasin, 1-hydroxy-N-methylphenazine

CAS 85-66-5

Couple Target ROS

Couple Type Dye

Formula $C_{13}H_{10}N_2O$

Formulation Lyophilized.

MW 210.2

Purity ≥98% (HPLC)

Quantity 1 μmol

Reconstitution Reconstitute the contents of the vial in 100µl solvent to

yield a 10mM stock solution.

Solubility Soluble in DMSO or dimethyl formamide.

Technical Info / Product Notes

Pyocyanin is a member of the CELLESTIAL[®] product line, reagents and assay kits comprising fluorescent molecular probes that have been extensively benchmarked for live cell analysis applications. $\mathsf{CELLESTIAL}^{\texttt{R}}$ reagents and kits are optimal for use in demanding imaging applications, such as confocal microscopy, flow cytometry and HCS, where consistency and reproducibility are required.

Last modified: May 29, 2024



eu@enzolifesciences.com

Phone: 800.942.0430

info-