

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 H_2O_2 directly in the lungs.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-53001-C001	1µmol
----------------	-------

Manuals, SDS & CofA

[View Online »](#)

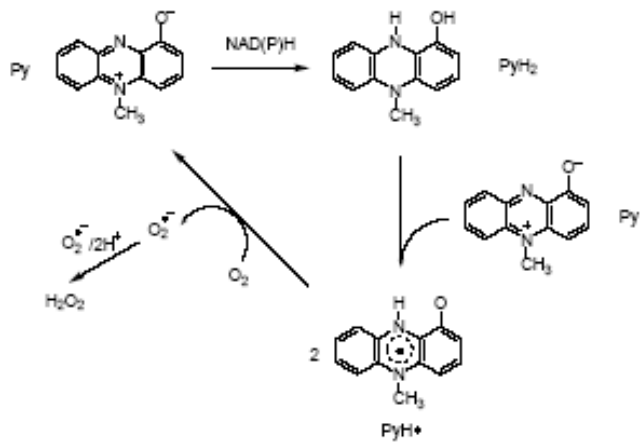
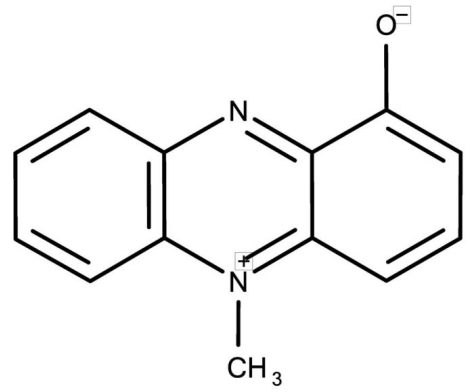


Figure: Pyocyanin is reduced by NAD(P)H and generates superoxide and H_2O_2 .



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 ₂ O
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. CELLESTIAL[®] 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.

ENZO LIFE SCIENCES, INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office (ELS) AG
info-eu@enzolifesciences.com

Belgium, The Netherlands
Phone: +32 3 466 0420
be@enzolifesciences.com

France
Phone: +33 4 67 40 60 60
info-fr@enzolifesciences.com

Germany
Phone: +49 622 760 630
info-de@enzolifesciences.com

UK & Ireland
Phone: 0845 601 1488
info-uk@enzolifesciences.com