

Mycophenolic acid

Inosine-5'-monophosphate dehydrogenase inhibitor

Antiviral and immunosuppressive activity. Clinically useful antipsoriatic agent. Mycophenolic acid prodrugs are effective in rheumatoid arthritis and for the prevention of organ transplant rejection. Mechanism of action is inhibition of inosine-5'-monophosphate dehydrogenase.

Citations: 4

[View Online »](#)

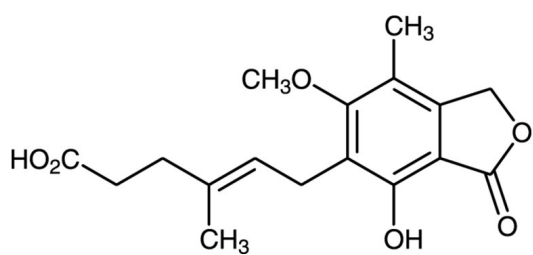
Ordering Information

[Order Online »](#)

BML-A249-0500	500mg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-(4-Hydroxy-6-methoxy-7-methyl-3-oxo-5-phthalanyl)-4-methyl-4-hexanoic acid
Appearance	White solid.
CAS	24280-93-1
Couple Target	Inosine-5'-monophosphate dehydrogenase
Couple Type	Inhibitor
Formula	$C_{17}H_{20}O_6$
MW	320.3
Purity	≥98% (HPLC)
RTECS	MP8050000
Solubility	Soluble in DMSO (clear colorless solution at 50 mg/ml) or 100% ethanol.
Source	Isolated from <i>Penicillium brevi-compactum</i>

Last modified: May 29, 2024



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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com