

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 »](#)

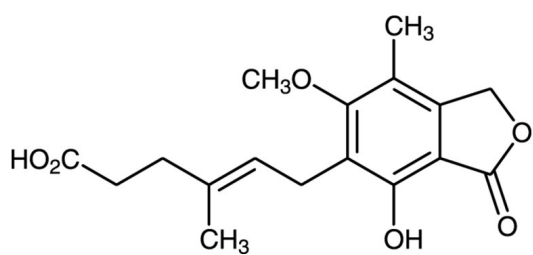
Ordering Information

[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> |



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-be@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