

Mycophenolate mofetil

Immunosuppressant

Mycophenolate mofetil (CellCept) is an immunosuppressant prodrug of mycophenolic acid, a fermentation product of *Penicillium brevicompactum*. Immunosuppression occurs via inhibition of type II inosine-5'-monophosphate dehydrogenase, depleting guanosine preferentially in T and B lymphocytes. In addition, mycophenolic acid displays anti-inflammatory activity by preventing NO production specifically from iNOS and by blocking glycosylation and expression of adhesion molecules and therefore recruitment of lymphocytes and monocytes to sites of inflammation.

Citations: 5

[View Online »](#)

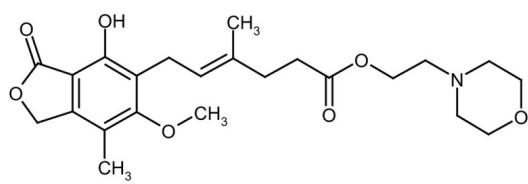
Ordering Information

[Order Online »](#)

BML-IM101-0020	20mg
BML-IM101-0100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Solutions can be stored for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	128794-94-5
Couple Target	Inosine-5'-monophosphate dehydrogenase, NOS
Couple Type	Inhibitor
Formula	$C_{23}H_{31}NO_7$
MW	433.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml, warm).
Source	Isolated from <i>Penicillium brevicompactum</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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