

# 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 »](#)

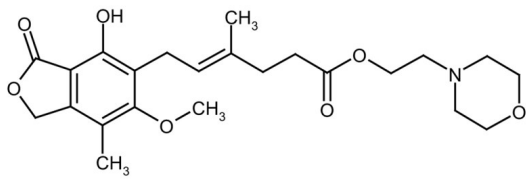
Ordering Information

[Order Online »](#)

BML-IM101-0020	20mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Appearance</b>	White to off-white solid.
<b>CAS</b>	128794-94-5
<b>Couple Target</b>	Inosine-5'-monophosphate dehydrogenase, NOS
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{23}H_{31}NO_7$
<b>MW</b>	433.5
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml, warm).
<b>Source</b>	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-  
be@enzolifesciences.com](mailto:info-be@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)