

Dexamethasone

Glucocorticoid receptor ligand

Anti-inflammatory glucocorticoid with anti-inflammatory activity. Induces the production of phospholipase A2 inhibitory protein (lipocortin). Inhibits induction of nitric oxide synthase (NOS) ($IC_{50}=5nM$). Induces apoptosis in eosinophils, but inhibits it in neutrophils. Both effects are blocked by RU-486. Modulates osteogenesis of mesenchymal stem cells. Inhibits constitutive OPG mRNA steady-state levels in a dose- and time-dependent fashion, promoting osteoclast formation. Oral dexamethasone administration induces an increase in insulin resistance in humans. Furthermore, it inhibits differentiation of immature to mature dendritic cells.

Citations: 36

[View Online »](#)

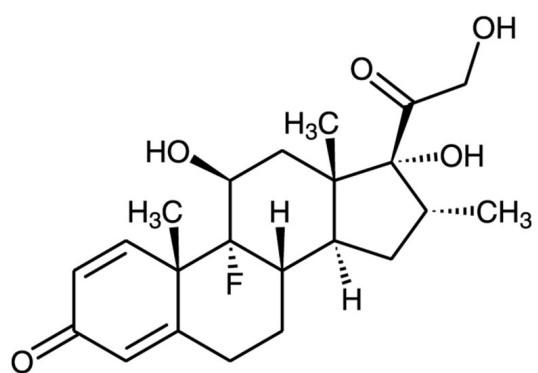
Ordering Information

[Order Online »](#)

BML-EI126-0001	1g
BML-EI126-0005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 5 years after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	9α-Fluoro-16α-methylprednisolone
Appearance	White to off-white solid.
CAS	50-02-2
Couple Target	Glucocorticoid receptor, IDO
Couple Type	Activator, Ligand
Formula	$C_{22}H_{29}FO_5$
MI	14: 2943
MW	392.5
Purity	≥98% (HPLC)
RTECS	TU3980000
Solubility	Soluble in DMSO (25mg/ml), 100% ethanol (25mg/1.6ml) or methanol; insoluble in water.



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)