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) (IC50=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 »

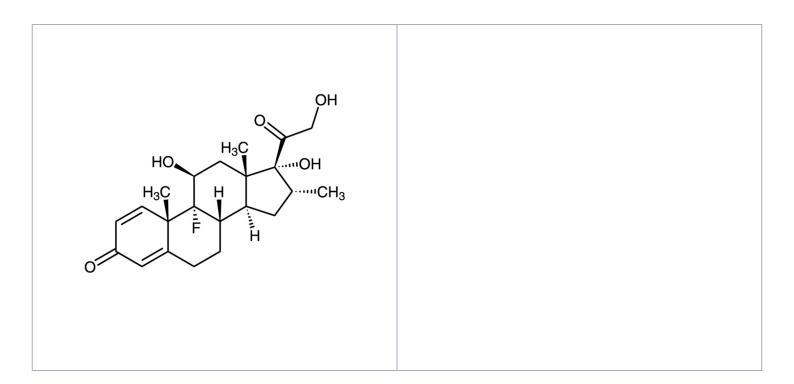
Ordering Information

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-16 α -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

Soluble in DMSO (25mg/ml), 100% ethanol (25mg/1.6ml)

or methanol; insoluble in water.