

Compound A

Glucocorticoid receptor modulator

Cell permeable, nonsteroidal glucocorticoid receptor modulator with high binding affinity. Exerts anti-inflammatory potential by down-modulating TNF-induced pro-inflammatory gene expression, such as IL-6 and E-selectin. Unlike dexamethasone (Prod. No. BML-EI126) it does not enhance glucocorticoid response element (GRE)-driven gene expression or induce glucocorticoid receptor (GR) binding to GRE-dependent genes *in vivo*.

Citations: 6

[View Online »](#)

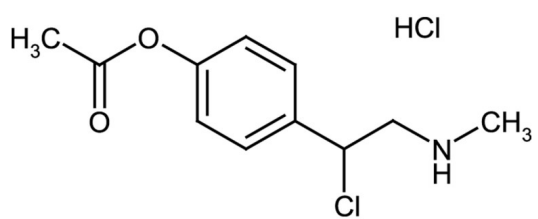
Ordering Information

[Order Online »](#)

ALX-550-516-M005	5mg
ALX-550-516-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in ethanol are stable for up to 1 month when stored at -80°C.
Handling	Protect from moisture. After reconstitution, prepare aliquots and store at -80°C.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CpdA, 2-((4-Acetophenyl)-2-chloro-N-methyl)ethylammonium chloride
Appearance	White to off-white solid.
CAS	14593-25-0
Formula	$C_{11}H_{14}ClNO_2 \cdot HCl$
MW	227.7 . 36.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol or water.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)