

Indomethacin

COX-1 inhibitor. PPAR activator.

Non-steroidal anti-inflammatory and analgesic agent. Inhibits cyclooxygenase ($IC_{50}=0.1\mu\text{M}$) selectively over lipoxygenases ($IC_{50}=100\mu\text{M}$ for 5-, 12- and 15-LO). Preferentially inhibits PGH synthase-1 over PGH synthase-2 ($ID_{50}=4.9-8.1$ and $130-160\mu\text{M}$ respectively). Activates PPAR α and γ receptors and induces adipocyte differentiation ($EC_{50}=8\mu\text{M}$). A clinically useful NSAID.

Citations: 11

[View Online »](#)

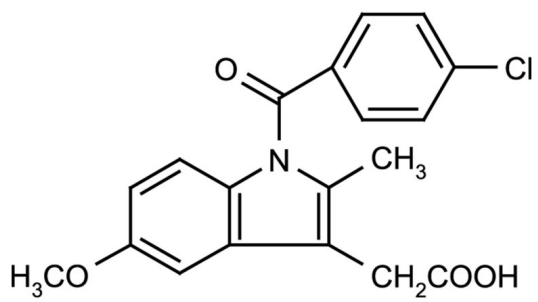
Ordering Information

[Order Online »](#)

ALX-270-086-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature – Dangerous Good

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to light yellow to light orange powder.

CAS 53-86-1

Couple Target COX, PPAR

Couple Type Activator, Inhibitor

Formula $C_{19}H_{16}ClNO_4$

MI **14:** 4968

MW 357.8

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in acetone or methanol; insoluble in 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-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)