

SC-560

COX inhibitor

Selective COX-1 inhibitor (IC_{50} =9nM). Shows 700-fold more selectivity over COX-2 (IC_{50} =6.3 μ M). Inhibits COX-1-derived platelet thromboxane B_2 , gastric PGE_2 , and dermal PGE_2 production. Orally active in rat.

Citations: 5

[View Online »](#)

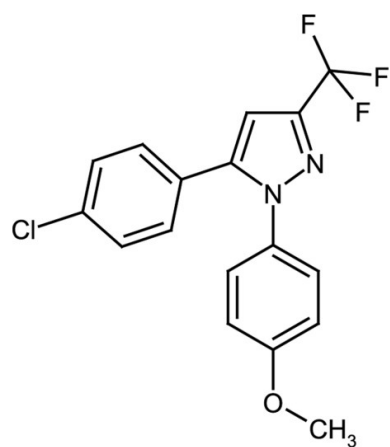
Ordering Information

[Order Online »](#)

ALX-270-427-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-(4-Chlorophenyl)-1-(4-methoxyphenyl)-3-(trifluoromethyl)-1H-pyrazole
Appearance	White to off-white crystalline solid.
CAS	188817-13-2
Couple Target	COX
Couple Type	Inhibitor
Formula	$C_{17}H_{12}N_2OCIF_3$
MW	352.7
Purity	≥98%
Solubility	Soluble in 100% ethanol, DMSO or dimethyl formamide; sparingly soluble in aqueous buffers.

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)