

# SC-236

## COX inhibitor

Highly selective and potent inhibitor of cyclooxygenase-2 ( $IC_{50}=10\text{nM}$  for COX-2 versus  $IC_{50}=17.8\mu\text{M}$  for COX-1) with antitumor properties. Exhibits longer half-life time and reduced gastric toxicity in fasting rat model. Has antitumor properties and has been shown to induce apoptosis and bFGF and VEGF-driven angiogenesis. Potent antimetastatic activity against both spontaneous metastases arising following primary tumour excision and experimental metastases.

Citations: 14

[View Online »](#)

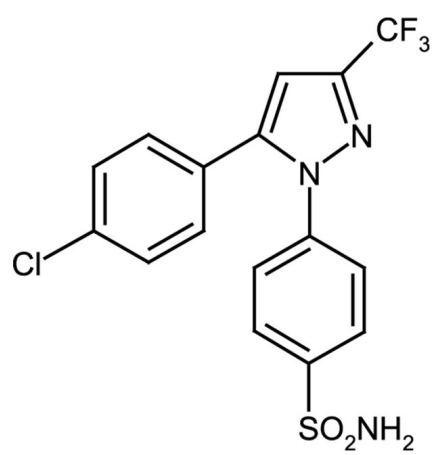
## Ordering Information

[Order Online »](#)

ALX-270-377-M005	5mg
------------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C. Stock solutions are stable for 3 months when stored at -20°C.
<b>Handling</b>	Protect from light. Packaged under inert gas.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	4-[5-(4-Chlorophenyl)-3-(trifluoromethyl)-1H-pyrazol-1-yl]benzenesulfonamide
<b>Appearance</b>	White to off-white crystalline powder.
<b>CAS</b>	170569-86-5
<b>Couple Target</b>	COX
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{16}H_{11}ClF_3N_3O_2S$
<b>MW</b>	401.8
<b>Purity</b>	≥95% (HPLC)
<b>Solubility</b>	Soluble in DMSO.

**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-fr@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)