

# SC-236

## COX inhibitor

Highly selective and potent inhibitor of cyclooxygenase-2 ( $IC_{50}$ =10nM for COX-2 versus  $IC_{50}$ =17.8 $\mu$ 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 »](#)

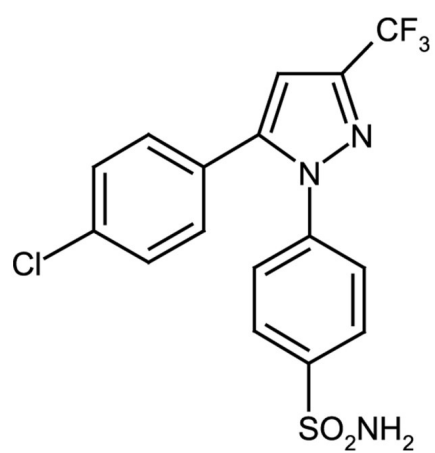
## Ordering Information

[Order Online »](#)

ALX-270-377-M005	5mg
ALX-270-377-M025	25mg

## 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. Stock solutions are stable for 3 months when stored at -20°C.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

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



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)