

# Celecoxib

## COX-2 inhibitor

Celecoxib is a non-steroidal anti-inflammatory drug (NSAID) that selectively inhibits cyclooxygenase-2 (COX-2) with an  $IC_{50}$  of 40 nM. It is commonly used to treat pain and inflammation associated with conditions like osteoarthritis and rheumatoid arthritis. Celecoxib is known for its lower risk of gastrointestinal side effects compared to non-selective NSAIDs.

Key features and applications include:

- **Selective COX-2 Inhibition:** Targets the COX-2 enzyme specifically, reducing pain and inflammation with fewer gastrointestinal side effects compared to non-selective NSAIDs.

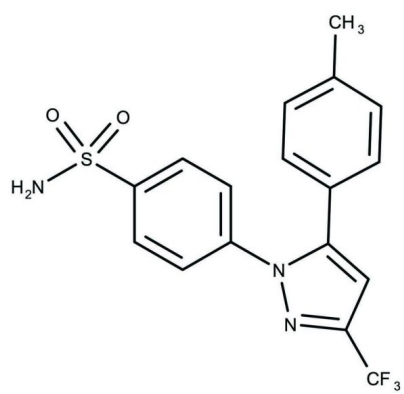
## Ordering Information

[Order Online »](#)

ENZ-CHM332-0250	250mg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	Ambient
Long Term Storage	Ambient
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	4-[5-(4-Methylphenyl)-3-(trifluoromethyl)-1H-pyrazol-1-yl]benzenesulfonamide, CJ-016377, CP-598107, PF-00345549, PHA-00846533, SC-58635 YM-177
Appearance	White solid.
CAS	169590-42-5
Couple Target	COX
Couple Type	Inhibitor
Formula	$C_{17}H_{14}F_3N_3O_2S$
Identity	Determined by NMR.
MW	381.38
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 100 mg/mL) or in ethanol (up to 100 mg/mL).

Last modified: July 28, 2025



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)