

# NS-398

## COX inhibitor

Cyclooxygenase inhibitor. Selectively inhibits PGH synthase-2 ( $IC_{50}=1\text{ }\mu\text{M}$ ) with no inhibition of PGH synthase-1 at concentrations up to  $100\text{ }\mu\text{M}$ . Protects neuronal cultures from LPS-induced neurotoxicity.

Citations: 14

[View Online »](#)

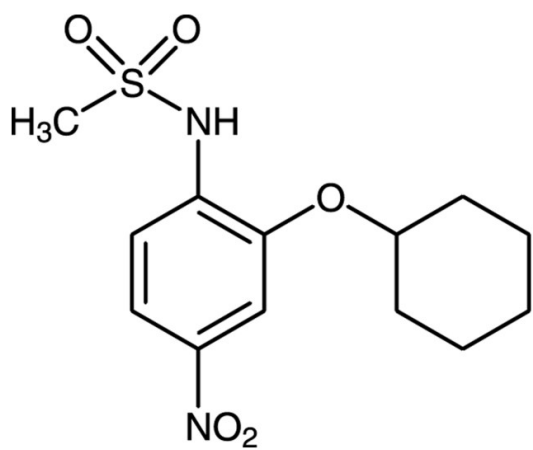
## Ordering Information

[Order Online »](#)

|                |      |
|----------------|------|
| BML-EI261-0005 | 5mg  |
| BML-EI261-0025 | 25mg |

## 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

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-[2-(Cyclohexyloxy)-4-nitrophenyl]-methanesulfonamide

**CAS** 123653-11-2

**Couple Target** COX

**Couple Type** Inhibitor

**Formula**  $C_{13}H_{18}N_2O_5S$

**MW** 314.4

**Purity**  $\geq 98\%$  (HPLC)

**Solubility** Soluble in 100% ethanol (warm; sonication), DMSO (25mg/ml) or dimethyl formamide. 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)