

# Meclofenamate sodium

COX inhibitor

Dual 5-lipoxygenase ( $IC_{50}=46.8\text{ }\mu\text{M}$ ) and cyclooxygenase I and II inhibitor ( $IC_{50}=0.6\text{ }\mu\text{M}$ ). Clinically useful NSAID.

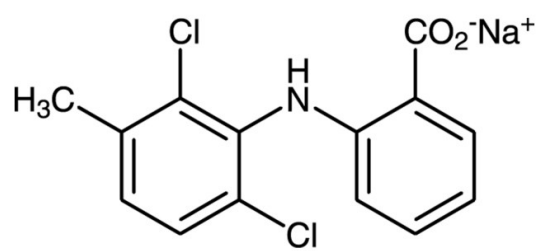
## Ordering Information

[Order Online »](#)

|                |    |
|----------------|----|
| BML-EI163-0005 | 5g |
|----------------|----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                                                                                                                         |
|-------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months. |
| Long Term Storage | Ambient                                                                                                                                                                 |
| Shipping          | Ambient Temperature                                                                                                                                                     |

## Regulatory Status

RUO - Research Use Only

## Product Details

|               |                                             |
|---------------|---------------------------------------------|
| Appearance    | White solid.                                |
| CAS           | 6385-02-0                                   |
| Couple Target | COX, Lipxygenase                            |
| Couple Type   | Inhibitor                                   |
| Formula       | $C_{14}H_{10}Cl_2NO_2 \cdot Na$             |
| MW            | 318.1                                       |
| Purity        | ≥98% (HPLC)                                 |
| Solubility    | Soluble in water (15mg/ml slightly cloudy). |

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)