

# Sulindac sulfide

COX inhibitor

Active metabolite of sulindac (Prod. No. ALX-430-105). Inhibits cyclooxygenase. More selective for PGHS-1 ( $ID_{50}=0.5 \mu\text{M}$ ) over PGHS-2 ( $ID_{50}=14 \mu\text{M}$ ). Induces apoptosis by a cyclooxygenase-independent mechanism. Inhibits colorectal cancer growth and  $\gamma$ -secretase. NF- $\kappa$ B inhibitor. Inhibits Ras signaling.

Citations: 6

[View Online »](#)

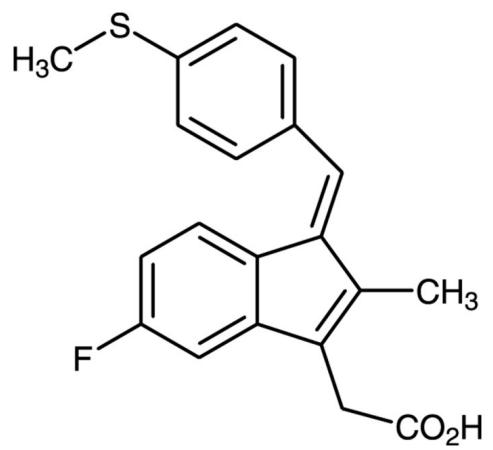
Ordering Information

[Order Online »](#)

BML-AP200-0005	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 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Z-5-Fluoro-2-methyl-1-[p-(methylthio)benzylidene]indene-3-acetic acid
<b>Appearance</b>	Yellow-orange solid.
<b>CAS</b>	32004-67-4
<b>Couple Target</b>	COX
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{20}H_{17}FO_2S$
<b>MW</b>	340.4
<b>Purity</b>	≥99% (TLC)
<b>Solubility</b>	Soluble in DMSO (>25mg/ml) or 100% ethanol (5mg/ml, warm).

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)