

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 γ -secretase. NF- κ B inhibitor. Inhibits Ras signaling.

Citations: 6

[View Online »](#)

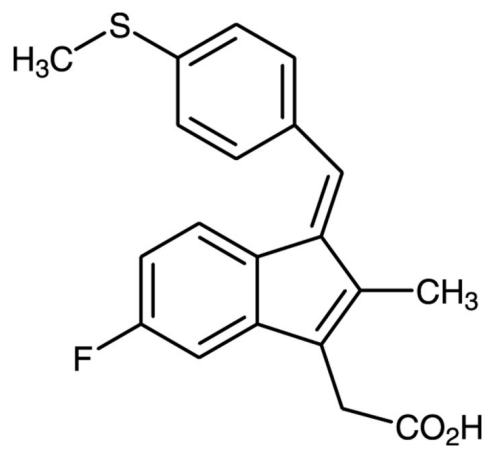
Ordering Information

[Order Online »](#)

BML-AP200-0005	5mg
BML-AP200-0025	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 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Z-5-Fluoro-2-methyl-1-[p-(methylthio)benzylidene]indene-3-acetic acid
Appearance	Yellow-orange solid.
CAS	32004-67-4
Couple Target	COX
Couple Type	Inhibitor
Formula	$C_{20}H_{17}FO_2S$
MW	340.4
Purity	≥99% (TLC)
Solubility	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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com