

Sulindac sulfone

Sulindac metabolite

Induces apoptosis but does not inhibit cyclooxygenase. Inhibits growth and induces apoptosis in human cancer cell lines. Metabolite of Sulindac (Prod. No. ALX-430-105).

Citations: 2

[View Online »](#)

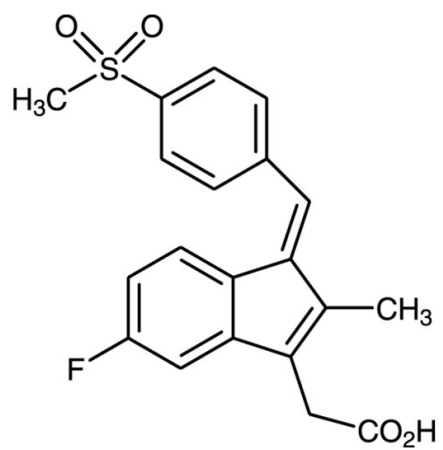
Ordering Information

[Order Online »](#)

BML-AP201-0005	5mg
BML-AP201-0025	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-(methyl-sulfonyl)benzylidene]indene-3-acetic acid
Appearance	Yellow solid.
CAS	59864-04-9
Formula	$C_{20}H_{17}FO_4S$
MW	372.4
Purity	≥94% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (3mg/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-fr@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