

DuP-697

COX inhibitor

DuP-697 is a highly selective and irreversible inhibitor of cyclooxygenase-2 (COX-2). (COX-2 IC_{50} =10nM, COX-1 IC_{50} =800nM). It is an orally effective anti-inflammatory, analgesic and antipyretic agent which is not ulcerogenic. Inhibition of COX-2 by DuP-697 and other inhibitors exacerbates ischemia-induced myocardial dysfunction in the rabbit. It inhibits angiogenesis and induces apoptosis in human endothelial cells.

Citations: 4

[View Online »](#)

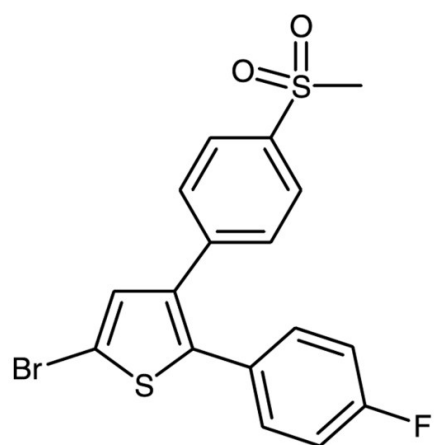
Ordering Information

[Order Online »](#)

BML-EI358-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (5-Bromo-2[4-fluorophenyl]-3-[4-methylsulfonyl-phenyl]thiophene)

Appearance White to off-white solid.

CAS 88149-94-4

Couple Target COX

Couple Type Inhibitor

Formula $C_{17}H_{12}BrFO_2S_2$

Identity Determined by NMR.

MW 411.3

Purity ≥98% (TLC)

Solubility Soluble in DMSO (10mg/ml) or 100% ethanol (5mg/ml).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)