

Phenylmethylsulfonyl fluoride

Inhibitor of serine proteases and FAAH

Irreversible inhibitor of serine proteases. Widely used fatty acid amid hydrolase (FAAH) inhibitor for pretreatment of brain membrane preparations when testing CB₁ receptor activity. Spleen membranes used for CB₂ studies do not require this pretreatment, as FAAH activity is absent in these cells. Inhibits also cysteine proteases like papain (reversible by DTT treatment, Prod. No. ALX-280-001), as well as internucleosomal DNA fragmentation in immature thymocytes. For a related, more stable inhibitor see AEBSF (Prod. No. ALX-270-022).

Citations: 13

[View Online »](#)

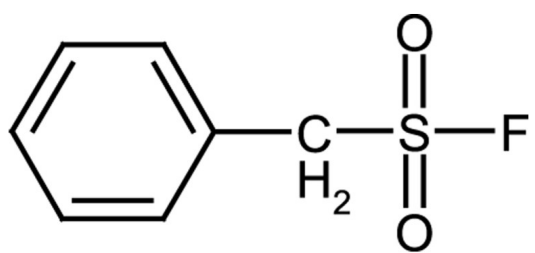
Ordering Information

[Order Online »](#)

ALX-270-184-G005	5g
ALX-270-184-G025	25g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PMSF, Benzylsulfonyl fluoride

Appearance White to off-white needle crystals.

CAS 329-98-6

Couple Target FAAH, Serine protease

Couple Type Inhibitor

Formula $C_7H_7FO_2S$

MW 174.2

Purity $\geq 99\%$ (HPLC)

RTECS XT8040000

Solubility Soluble in 100% ethanol, methanol or isopropanol; insoluble in water.

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)