

AEBSF . hydrochloride

Serine protease inhibitor

Specific irreversible inhibitor of serine proteases including trypsin, chymotrypsin, plasmin, plasma kallikrein, and thrombin. Stable, non-toxic alternative to PMSF (Prod. No. ALX-270-184). Inhibitor of NAD(P)H oxidase. Activator of MAP kinases JNK and p38.

Citations: 14

[View Online »](#)

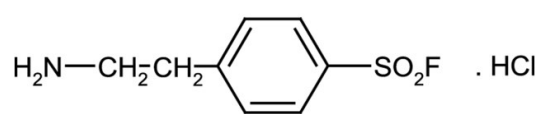
Ordering Information

[Order Online »](#)

ALX-270-022-M250	250mg
ALX-270-022-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from moisture and light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-(2-Aminoethyl)benzenesulfonylfluoride . HCl
Appearance	White to off-white powder.
CAS	30827-99-7
Couple Target	Jnk, NADPH oxidase, p38, Serine protease
Couple Type	Activator, Inhibitor
Formula	$C_8H_{10}FNO_2S \cdot HCl$
MW	239.7
Purity	≥98% (HPLC)
Solubility	Soluble in water or 100% ethanol.

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-
fr@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)