

(S)-ESBA . HCl

Kynurenine aminotransferase inhibitor

Potent and selective mouse/rat kynurenine aminotransferase II (KAT II) inhibitor. $IC_{50}=6.1\mu M$.

Citations: 5

[View Online »](#)

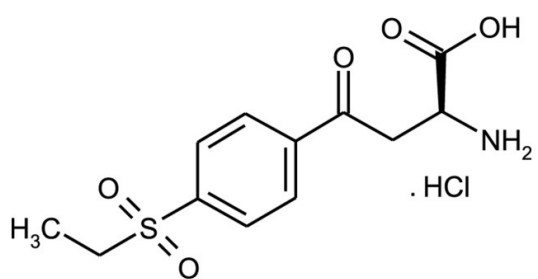
Ordering Information

[Order Online »](#)

ALX-550-529-M005	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 (S)-(4-Ethylsulfonyl)benzoylalanine . hydrochloride

Appearance White to off-white powder.

Couple Target Kynurenine aminotransferase

Couple Type Inhibitor

Formula $C_{12}H_{15}NO_5S \cdot HCl$

MW 285.3 . 36.5

Purity ≥97% (HPLC)

Solubility Soluble in water or DMSO.

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)