

L-Kynurenine

Key intermediate in the breakdown pathway of tryptophan.

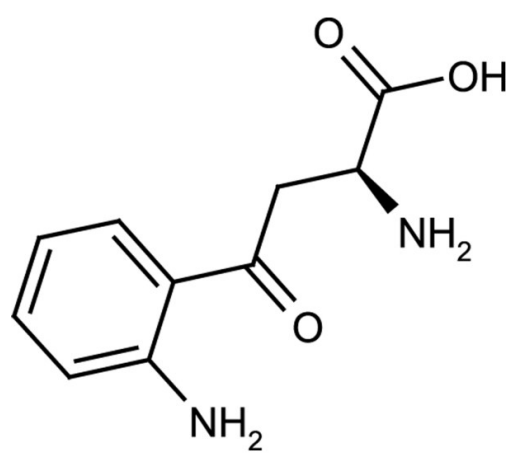
Ordering Information

[Order Online »](#)

ALX-550-408-M050	50mg
------------------	------

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 β -Anthraniloyl-L-alanine, L-2-Amino-4-(2-aminophenyl)-4-oxobutanoic acid

Appearance Light yellow solid.

CAS 2922-83-0

Formula $C_{10}H_{12}N_2O_3$

MI 14: 5328

MW 208.2

Purity $\geq 95\%$ (HPLC)

Solubility Soluble in 0.5 M hydrochloric acid (50mg/ml); very poorly soluble in water (0.4g/100ml, 20°C).

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)