

BEC . ammonium salt

Arginase inhibitor

Slow-binding competitive inhibitor of arginase I ($K_i=0.4-0.6\mu\text{M}$ & $K_D=2.22\mu\text{M}$ for recombinant rat liver arginase) and arginase II ($K_i=0.31\mu\text{M}$ at pH 7.5 and $0.03\mu\text{M}$ at pH 9.5 for human recombinant type II arginase). Enhances NO-dependent smooth muscle relaxation in human penile corpus cavernosum tissue. Does not inhibit nitric oxide synthase (NOS) and may serve as a valuable reagent to probe the physiological relationship between arginase and NOS. Inhibitor of rat aortic smooth muscle cell proliferation.

Citations: 10

[View Online »](#)

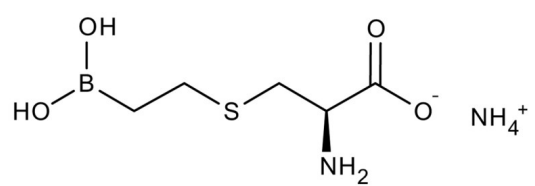
Ordering Information

[Order Online »](#)

| | |
|------------------|-----|
| ALX-270-345-M001 | 1mg |
| ALX-270-345-M005 | 5mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. |
| Handling | Very hygroscopic. Keep under inert gas. Keep cool and dry. |
| Long Term Storage | +4°C |
| Shipping | Ambient Temperature |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|------------------|---|
| Alternative Name | S-(2-Boronoethyl)-L-cysteine . NH ₄ |
| Appearance | White to off-white powder. |
| CAS | 63107-40-4 (non-salt) |
| Couple Target | Arginase |
| Couple Type | Inhibitor |
| Formula | C ₅ H ₁₁ BNO ₄ S . NH ₄ |
| MW | 192.0 . 18.0 |
| Purity | ≥95% (TLC) |
| Solubility | Soluble in water (10mg/ml). |



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)