# **BEC**. ammonium salt

#### **Arginase inhibitor**

Slow-binding competitive inhibitor of arginase I ( $K_i$ =0.4-0.6µM &  $K_D$ =2.22µM for recombinant rat liver arginase) and arginase II ( $K_i$ =0.31µM at pH 7.5 and 0.03µM 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 »

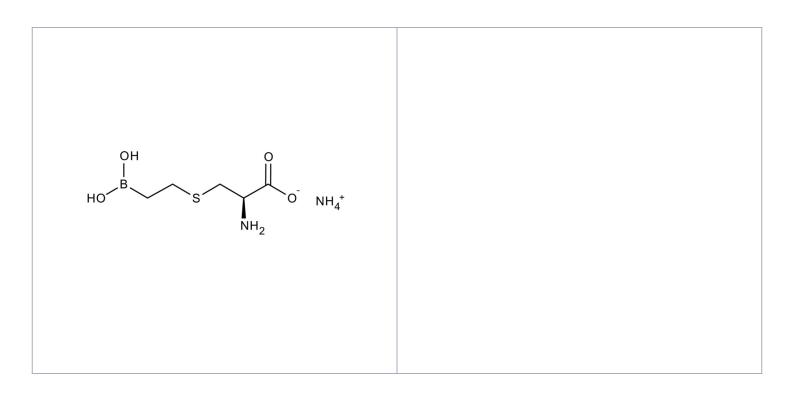
## **Ordering Information**

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 . NH4

**Appearance** White to off-white powder.

**CAS** 63107-40-4 (non-salt)

Couple Target Arginase

Couple Type Inhibitor

Formula  $C_5H_{11}BNO_4S \cdot NH_4$ 

**MW** 192.0 . 18.0

**Purity** ≥95% (TLC)

**Solubility** Soluble in water (10mg/ml).

Last modified: May 29, 2024