

Elafin (human), (recombinant)

Belongs to the chelonianin family of protease inhibitors and is involved in the physiological control of neutrophil serine proteinases, where it targets NE and proteinase 3. Shows antiprotease, antibacterial and antiinflammatory activity and primes innate immunity.

Citations: 20

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-2400	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ESI, Elastase-specific inhibitor, SKALP, Skin-derived antileukoprotease, Protease inhibitor WAP3
Appearance	White to faint yellow powder.
Biological Activity	Inhibits human leukocyte elastase ($K_i=0.17\text{nM}$) and human proteinase 3 ($K_i=0.42\text{nM}$).
Formulation	Lyophilized.
MW	~5.9kDa
Purity	≥90% (HPLC, SDS-PAGE, MS)
Reconstitution	Reconstitute in sterile distilled water. Further dilutions should be made in aqueous buffers, pH 4.0-8.0.
Source	Produced in yeast.
Specific Activity	Inhibition of human leukocyte elastase: ≥90%
UniProt ID	P19957

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com