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-201-240-C100

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

antileukoproteinase, Protease inhibitor WAP3

Appearance White to faint yellow powder.

Biological Activity Inhibits human leukocyte elastase (K_i=0.17nM) and human

proteinase 3 (K_i=0.42nM) .

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

