

# Procathepsin K (human), (recombinant)

Member of the papain cysteine proteinase family identified as the predominant proteinase responsible for the resorption of the bone matrix.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-201-239-C010	10µg
------------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable for several months when stored at -80°C. After activation, cathepsin is highly autoproteolytic. When not used immediately, add MMTS (Methyl methanethiosulfonate) to 1mM final concentration for stabilization and shock freeze in liquid nitrogen. Do not store dilutions of the enzyme.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Procathepsin O, Procathepsin O2, Procathepsin X
<b>Concentration</b>	150µg/ml
<b>Formulation</b>	Liquid. In 25mM TRIS, pH 8.0, containing 500mM sodium chloride.
<b>MW</b>	~35kDa
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Source</b>	Produced in <i>E. coli</i> (aa 19-329). Methionine was introduced at aa 18 to create a new N-terminal sequence (MEEIL).
<b>Specific Activity</b>	≥1'000mU/mg protein. One unit is defined as the amount of enzyme that hydrolyzes 1µmol of substrate per min. at 37°C, pH 5.5.
<b>Technical Info / Product Notes</b>	<b>Inhibitors:</b> Cathepsin K is inhibited by leupeptin (Prod. No. ALX-260-009) ( $IC_{50}$ =70nM), E-64 ( $IC_{50}$ =5nM) or cystatine. Minimal effects exhibit pepstatin, phenylmethylsulfonyl fluoride (Prod. No. ALX-270-184), inhibitors of aspartyl and serine proteases. No inhibition was observed by addition of EDTA or phenanthroline, classical inhibitors of metalloproteases.
<b>UniProt ID</b>	P43235



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-be@enzolifesciences.com](mailto:info-be@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)