

# Procathepsin V (human), (recombinant) (His- tag)

Full-length proenzyme. Can be activated in one step.

Cathepsin V, a member of the papain family of cysteine proteases, has 78% identity to cathepsin L, but unlike cathepsin L is not widely expressed, being localized to thymus, testis, corneal epithelium, and macrophages. It is a strong elastase, also cleaving proteins such as the invariant chain (Ii), plasminogen, and kininogen. It is implicated in disease states such as cancer, angiogenesis, atherosclerosis, and myasthenia gravis.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-SE554-0010	10µg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Avoid extended periods unfrozen. This enzyme is stable for 6 months when stored as received under the above conditions.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Cathepsin L2
<b>Application Notes</b>	Useful tool to study enzyme kinetics, cleave target substrates and screen for inhibitors.
<b>Formulation</b>	Liquid. In 25mM TRIS-HCl, pH 8.0, containing 100mM sodium chloride, 0.05% Tween-20 and 10% glycerol.
<b>MW</b>	~37kDa (SDS-PAGE)
<b>Purity</b>	≥90%
<b>Source</b>	Produced in insect cells. Recombinant glycosylated procathepsin V cloned from human cDNA and purified as full-length proenzyme. Produced in a baculovirus expression system.
<b>Specific Activity</b>	436 U/μg protein. One unit will hydrolyze one pmole Z-Leu-Arg-AMC substrate (Prod. No. BML-P229, 25 μM) per minute at 25°C, in 25mM NaOAc pH 5.5, 100mM NaCl, 1.0mM DTT.
<b>Technical Info / Product Notes</b>	The proenzyme can be activated as in Adachi et al. (Reference year 2008): Dilute proenzyme into 200mM NaAcetate, pH 6.0, with 0.05% SDS, 2.5mM DTT, and incubate at 37°C for 5-30 minutes. Alternatively, pre-incubate proenzyme 5-30 minutes in assay buffer (25mM NaOAc pH 5.5, 100mM sodium chloride, 1.0mM DTT), then add substrate to begin assay. Incubation times must be determined empirically; activation is dependent on factors such as buffer, temperature, and enzyme concentration, and cathepsin V will autodegrade once activated.



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)