Procathepsin K (human), (recombinant)

Cathepsin K, a member of the papain family of cysteine proteases, acts upon proteins such as collagen, AL amyloid, kinin and elastin. It is a lysosomal protease expressed primarily in osteoclasts, but also in other cell types such as macrophages. It functions in bone remodeling and is implicated in disease states such as atherosclerosis, arthritis and pycnodysostosis.

Citations: 1

View Online »

Ordering Information

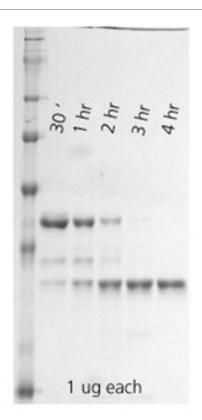
Order Online »

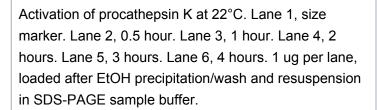
BML-SE367-0010

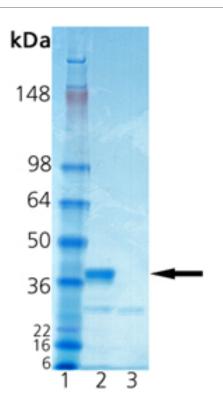
10µg

Manuals, SDS & CofA

View Online »







SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0 μg of purified Procathepsin K; Lane 3: 1.0 μg Cathepsin K after activation at low pH for 3 hours.

Handling & Storage

Use/Stability -80°C. After initial defrost, aliquot product into individual tubes and refreeze at -80°C.

Avoid repeated freeze/defrost cycles and extended periods unfrozen. This enzyme is stable when stored as received under the above conditions. Procedures such as dil

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity To activate proenzyme, add 32.5mM NaOAc, pH 3.5, to

20% of procathepsin K volume (e.g., 5µl procathepsin K +

1µI 32.5mM NaOAc, pH 3.5), and incubate at room temperature for 25-30 minutes, or at +4°C for

approximately seven hours.

Alternative Name Procathepsin O, Procathepsin O2, Procathepsin X

Application Notes Useful tool to study enzyme kinetics, cleave target

substrates and screen for inhibitors.

Formulation Liquid. In 27.8µl 20mM sodium phosphate, pH 6.9, 500mM

NaCl, 1 mM EDTA.

MW 35.3kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in insect cells. Full length glycosylated human

procathepsin K (also known as cathepsin O or cathepsin O2) cloned from human cDNA. Produced in a baculovirus

expression system.

UniProt ID P43235

Last modified: May 29, 2024



info-

eu@enzolifesciences.com