## Cathepsin H (human liver), (purified)

Cathepsin H is a lysosomal protease (monopeptidase and aminopeptidase) that can also be secreted from the cell and is involved in protein degradation and cancer.

Citations: 2

View Online »

**Ordering Information** 

Order Online »

BML-SE200-0025

25µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -80°C

**Shipping** Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

**Application Notes** Useful for inhibitor screening, kinetic studies, and cleavage

of substrates in vitro or in tissue culture.

Source Isolated from human liver.

**UniProt ID** P09668

Last modified: May 29, 2024



info-

eu@enzolifesciences.com