

Cathepsin L1 (human), (recombinant)

Cathepsin L1 plays an important role in the overall degradation of proteins in lysosomes. Its substrates include collagen and elastin, as well as α -1 protease inhibitor, a major controlling element of neutrophil elastase activity. It has been implicated in several pathologic processes, including myofibril necrosis in myopathies and in myocardial ischemia, and in the renal tubular response to proteinuria.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CTSL1, MEP, CATL, CTSL

Formulation Lyophilized from 0.22µm filtered solution in 50mM sodium acetate, pH 5.0, containing 100mM NaCl. Generally 5-8% mannitol or trehalose is added as a protectant before lyophilization.

MW ~36.7 kDa

Purity ≥98% (SDS-PAGE)

Source Produced in HEK 293 cells.

UniProt ID P07711

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com