

Cathepsin K drug discovery kit

The Cathepsin K Drug Discovery Kit is a complete assay system designed to screen inhibitors of cathepsin K. Using a convenient microplate format, cathepsin K activity is monitored fluorimetrically using the substrate provided in the kit.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK430-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The enzyme component should be thawed as close to the time of use as is possible, and held on ice while in use. The enzyme can be frozen and thawed up to 5 times, but care should be taken to minimize both the time the enzyme spends on ice, and the number of times it is frozen and thawed. Do not hold the diluted enzyme at assay temperature (e.g 37°C) for an extended time before starting the assay.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Activity assay, Fluorescent detection, HTS
Application Notes	Screening potential activators or inhibitors of cathepsin K using the recombinant enzyme supplied in the kit. This kit is NOT suitable for measuring cathepsin K activity in biological samples
Contents	Cathepsin K (human) (recombinant), AMC substrate, AMC calibration standard, inhibitor, assay buffer, 1/2-volume white microplate.
UniProt ID	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)