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

Last modified: May 29, 2024

