

# Granzyme B assay kit for drug discovery

Granzyme B is a serine protease found in the lymphocyte granules of cytotoxic T-lymphocytes (CTLs) and natural killer (NK) cells. Activated granzyme B prefers an aspartic acid residue at the P1 site of its substrates; a preference shared by only one other group of mammalian proteases, the caspases. Granzyme B is required for the rapid apoptotic signal delivered by CTLs and NK cells, an effect mediated by its ability to cleave and activate various procaspases. Granzyme B plays a crucial role in the development of acute graft-vs.-host disease. Thus, the inhibition of this enzyme may prove to be an important tool in controlling immune responses. Conversely, the failure of granzyme B-induced apoptosis may contribute to certain pathologies. Such apoptosis is important for the elimination of virus-infected and malignant cells and resistance to granzyme B action, conferred by the serpin PI-9, may aid in the escape of some tumors from immune surveillance. The Granzyme B Assay Kit for Drug Discovery is a complete assay system designed to measure the protease activity of granzyme B. It contains a colorimetric substrate (IEPD-pNA) to measure this activity; cleavage of the p-nitroanilide (pNA) group from the colorimetric substrate increases absorption at 405nm. The assays are performed in a convenient, 96-well microplate format. The kit is useful to screen for inhibitors of granzyme B, a potential therapeutic target. An inhibitor, 3, 4-Dichloroisocoumarin, is also included for use as a control.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-AK711-0001

96 wells

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Activity assay, Colorimetric detection, HTS

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**UniProt ID** P10144

**Last modified: January 9, 2025**



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)