

## Technical Bulletin

# Granzyme B Activity Assay Kit

**Catalog Number MAK176**

## Product Description

cytotoxic T lymphocyte (CTL)-mediated DNA fragmentation and apoptosis in target cells. Granzyme B, a serine protease, cleaves proteins after aspartic acid residues. Granzyme B-mediated release of mitochondrial cell death factors are required for the activation of caspase-3.<sup>1</sup> Furthermore, the degradation of granzyme B is known to decrease the susceptibility of cancer cells to NK-cell-mediated lysis during hypoxia.<sup>2</sup> Granzyme B activity assays can be used for studying T cell immunity, CTL responses, cell-mediated cytotoxicity, and apoptosis.

The Granzyme B Activity Assay Kit provides a rapid, reliable, and sensitive procedure for measuring granzyme B activity in a variety of samples. Granzyme B activity is determined using an enzyme-catalyzed reaction in which the non-fluorescent substrate, Ac-IEPD-AFC, is hydrolyzed resulting in the release of AFC ( $\lambda_{\text{Ex}} = 380 \text{ nm}/\lambda_{\text{Em}} = 500 \text{ nm}$ ) proportional to the enzymatic activity present. One unit of granzyme B is the amount of enzyme that hydrolyzes 1 pmol of the substrate per minute at 37 °C.

## Components

The kit is sufficient for 100 assays in 96-well plates.

- |   |        |
|---|--------|
| • Granzyme B Assay Buffer<br>Catalog Number MAK176A     | 25 mL  |
| • Granzyme B Substrate<br>Catalog Number MAK176B        | 0.5 mL |
| • Granzyme B Positive Control<br>Catalog Number MAK176C | 1 vial |
| • AFC Standard, 1 mM<br>Catalog Number MAK176D          | 0.1 mL |

## Reagents and Equipment Required but Not Provided

- 96-well flat-bottom plate – It is recommended to use black plates with clear bottoms for fluorescence assays. Cell culture or tissue culture treated plates are **not** recommended.
- Fluorescence multiwell plate reader

## Precautions and Disclaimer

For Research Use Only. Not for use in diagnostic procedures. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

## Storage/Stability

The kit is shipped on wet ice. Store components at 2–8 °C, protected from light.

## Preparation Instructions

Briefly centrifuge vials prior to opening. To maintain reagent integrity, avoid repeated freeze/thaw cycles.

**Granzyme B Assay Buffer:** Allow buffer to come to room temperature before use.

**Granzyme B Positive Control:** Reconstitute with 20  $\mu\text{L}$  of Granzyme B Assay Buffer. Mix well by pipetting (Do Not Vortex), then store in single-use aliquots, protected from light, at  $-20\text{ }^{\circ}\text{C}$ . Use within 1 month of reconstitution.

## Procedure

All samples and standards should be run in duplicate.

### Sample Preparation

Tissue (50 mg) or cells ( $1 \times 10^6$ ) can be homogenized in 200  $\mu\text{L}$  of ice-cold Granzyme B Assay Buffer. Centrifuge the samples at  $13,000 \times g$  for 10 minutes to remove insoluble material.

Serum samples can be directly added to wells.

**Note:** For unknown samples, it is suggested to test several sample dilutions to ensure the readings are within the linear range of the standard curve.

For all samples, bring samples to a final volume of 50  $\mu\text{L}$  per well with Granzyme B Assay Buffer.

### Positive Control

Add 2  $\mu\text{L}$  of the Positive Control into the wells and adjust to a final volume of 50  $\mu\text{L}$  per well with Granzyme B Assay Buffer.

### Standard Curve Preparation

1. Prepare a 10  $\mu\text{M}$  (10 pmol/ $\mu\text{L}$ ) AFC Standard by diluting 10  $\mu\text{L}$  of the 1 mM AFC Standard with 990  $\mu\text{L}$  of the Granzyme B Assay Buffer.
2. Prepare AFC Standards according to Table 1. Mix well.

**Table 1.**

Preparation of Resorufin Standards

Well	10 $\mu\text{M}$ AFC Standard	Granzyme B Assay Buffer	AFC (pmol/well)
1	0 $\mu\text{L}$	100 $\mu\text{L}$	0
2	5 $\mu\text{L}$	95 $\mu\text{L}$	50
3	10 $\mu\text{L}$	90 $\mu\text{L}$	100
4	15 $\mu\text{L}$	85 $\mu\text{L}$	150
5	20 $\mu\text{L}$	80 $\mu\text{L}$	200
6	25 $\mu\text{L}$	75 $\mu\text{L}$	250

### Assay Reaction

1. Set up the Reaction Mix according to the scheme in Table 2. 50  $\mu\text{L}$  of the Reaction Mix is required for each Sample and Positive Control well.

**Table 2.**

Reaction Mix

Reagent	Volume
Granzyme B Assay Buffer	45 $\mu\text{L}$
Granzyme B Substrate	5 $\mu\text{L}$

2. Add 50  $\mu\text{L}$  of the Reaction Mix to each of the Sample and Positive Control wells. Mix well using a horizontal shaker or by pipetting.
3. Incubate the plate at  $37\text{ }^{\circ}\text{C}$ . After 2-3 minutes, take the initial measurement ( $T_1$ ). Measure the fluorescence intensity ( $\text{RFU}_1$ ) using  $\lambda_{\text{Ex}} = 380\text{ nm}$  /  $\lambda_{\text{Em}} = 500\text{ nm}$ .

**Note:** It is essential the initial reading ( $\text{RFU}_1$ ) is within the linear range of the Standard curve.



4. Continue to incubate the plate at 37 °C taking measurements (RFU) every 5 minutes. Protect the plate from light during the incubation.
5. Continue taking measurements until the value of the most active Sample is greater than the value of the highest Standard (250 pmol/well). At this time the most active Sample is near or exceeds the end of the linear range of the Standard curve.
6. The final measurement (RFU<sub>2</sub>) for calculating the enzyme activity would be the penultimate reading or the value before the most active Sample is near or exceeds the end of the linear range of the Standard curve (See Step 5). The time at which this reading is taken is designated T<sub>2</sub>.

Note: It is essential the final measurement (RFU<sub>2</sub>) falls within the linear range of the Standard curve.

## Results

### Calculations

1. Correct for the background by subtracting the final measurement (RFU<sub>2</sub>) obtained for the 0 (blank) AFC Standard (Well 1) from the final measurement (RFU<sub>2</sub>) of each of the Standards and Samples. Background values can be significant and must be subtracted from all readings.
2. Plot the AFC Standard curve using the corrected (RFU<sub>2</sub>) values.

Note: A new Standard curve must be set up each time the assay is run.

3. Calculate the change in Fluorescent measurement (ΔRFU) from T<sub>1</sub> to T<sub>2</sub> for each of the Samples.

$$\Delta\text{RFU} = \text{RFU}_2 - \text{RFU}_1$$

4. Compare the ΔRFU of each Sample to the Standard curve to determine the amount (B) of AFC generated by the Granzyme B assay between T<sub>1</sub> and T<sub>2</sub>.
5. Calculate the Granzyme B activity of the Sample by the following equation:

$$\text{Granzyme B Activity (pmol/min/mL)} = \frac{\mu\text{U/mL}}{\Delta T \times V}$$

$$\frac{B \times DF}{\Delta T \times V}$$

where

B = Amount (pmol) of AFC generated between T<sub>1</sub> and T<sub>2</sub>.

ΔT = Reaction time T<sub>2</sub> – T<sub>1</sub> (minutes)

V = Sample volume (mL) added to well

DF = Dilution factor of the Sample.

DF = 1 for undiluted Samples

One unit of Granzyme B is the amount of enzyme that will generate 1.0 pmol of AFC from the substrate per minute at 37 °C.

### Example

AFC amount (B) = 5,840 pmol  
First reading (T<sub>1</sub>) = 3 minutes  
Second reading (T<sub>2</sub>) = 32 minutes  
Sample volume (V) = 0.05 mL  
Sample dilution is 1

For this example, Granzyme B activity is:

$$\frac{5,840 \times 1}{(32 - 3) \times 0.05} = 4028 \mu\text{U/mL}$$

## References

1. Lord, S.J. et al., Granzyme B: a natural born killer. *Immunol. Rev.*, **193**, 31-38 (2003).
2. Lindner, S., et al., Interleukin 21-induced granzyme B-expressing B cells infiltrate tumors and regulate T cells. *Cancer Res.*, **73(8)**, 2468-2479 (2013).



## Troubleshooting Guide

Problem	Possible Cause	Suggested Solution
Assay not working	Cold assay buffer	Assay Buffer must be at room temperature
	Omission of step in procedure	Refer and follow Technical Bulletin precisely
	Plate reader at incorrect wavelength	Check filter settings of instrument
	Type of 96 well plate used	For Fluorometric assays, use black plates with clear bottoms
Samples with erratic readings	Samples prepared in different buffer	Use the Assay Buffer provided or refer to Technical Bulletin for instructions
	Cell/Tissue culture samples were incompletely homogenized	Repeat the sample homogenization, increasing the length and extent of homogenization step.
	Samples used after multiple freeze-thaw cycles	Aliquot and freeze samples if needed to use multiple times
	Presence of interfering substance in the sample	If possible, dilute sample further
	Use of old or inappropriately stored samples	Use fresh samples and store correctly until use
Lower/higher readings in samples and standards	Improperly thawed components	Thaw all components completely and mix gently before use
	Use of improperly stored reagents	Check the storage temperature prior to using the kit and store the components appropriately
	Allowing the reagents to sit for extended times on ice	Prepare fresh Master Reaction Mix before each use
	Incorrect incubation times or temperatures	Refer to Technical Bulletin and verify correct incubation times and temperatures
	Incorrect volumes used	Use calibrated pipettes and aliquot correctly
Non-linear standard curve	Use of partially thawed components	Thaw and resuspend all components before preparing the reaction mix
	Pipetting errors in preparation of standards	Avoid pipetting small volumes
	Pipetting errors in the Reaction Mix	Prepare a Master Reaction Mix whenever possible
	Air bubbles formed in well	Pipette gently against the wall of the plate well
	Standard stock is at incorrect concentration	Refer to the standard dilution instructions in the Technical Bulletin
	Calculation errors	Recheck calculations after referring to Technical Bulletin
	Substituting reagents from older kits/lots	Use fresh components from the same kit
Unanticipated results	Samples measured at incorrect wavelength	Check the equipment and filter settings
	Samples contain interfering substances	If possible, dilute sample further
	Sample readings above/below the linear range	Concentrate or dilute samples so readings are in the linear range

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

## Technical Service

Visit the tech service page on our web site at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

## Standard Warranty

The applicable warranty for the products listed in this publication may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).