

Product Information

Thrombin Activity Fluorometric Assay Kit

Catalog Number **MAK242**
Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

Thrombin (EC 3.4.21.5, Factor IIa), a serine protease, is an important clotting factor in the coagulation cascade that involves the conversion of soluble fibrinogen to insoluble active fibrin strands. In this pathway, prothrombin is proteolytically converted into active thrombin.

Thrombin is also a potent vasoconstrictor and mitogen implicated as a major factor in vasospasm following subarachnoid hemorrhage. Ruptured cerebral aneurysm blood clots around a cerebral artery releases thrombin, which in turn induces acute and prolonged narrowing of the blood vessel, potentially resulting in cerebral ischemia and infarction (stroke). In addition, it is a pro-inflammatory enzyme that may influence the onset and progression of atherosclerosis.

This Thrombin Activity Fluorometric Assay Kit utilizes the ability of thrombin to proteolytically cleave a synthetic substrate and release a fluorophore, AMC, which can be easily quantified by fluorescence. This assay kit is simple, rapid, and can detect thrombin activity as low as 1 ng in samples.

Components

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

Thrombin Dilution Buffer Catalog Number MAK242A	1 mL
Thrombin Assay Buffer Catalog Number MAK242B	15 mL
Thrombin Enzyme Standard Catalog Number MAK242C	5 μL
Thrombin Substrate Catalog Number MAK242D	0.5 mL

Reagents and Equipment Required but Not Provided.

- 96 well flat-bottom plate – white plates are recommended for this assay.
- Fluorescence multiwell plate reader

Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Preparation Instructions

Briefly centrifuge small vials at low speed prior to opening.

Thrombin Assay Buffer – Bring to room temperature before use.

Thrombin Enzyme Standard – Prepare a stock solution of Thrombin Enzyme (50 ng/ μL) by adding 12 μL of Thrombin Dilution buffer to 4 μL of Thrombin Enzyme Standard. Mix, aliquot, and store at $-80\text{ }^{\circ}\text{C}$. Avoid repeated freeze/thaw cycles.

Storage/Stability

The kit is shipped on wet ice and storage at $-20\text{ }^{\circ}\text{C}$, protected from light, is recommended. Briefly centrifuge all small vials prior to opening.

Procedure

Read entire protocol before performing the assay.

Sample Preparation

Add 2–50 μL of sample containing thrombin per well of 96 well plate and adjust the volume to 50 μL with Thrombin Assay Buffer.

Standard Curve Preparation

Dilute Thrombin Enzyme Standard to 2.5 $\text{ng}/\mu\text{L}$ by adding 38 μL of Thrombin Dilution Buffer to 2 μL of Thrombin Enzyme stock solution (50 $\text{ng}/\mu\text{L}$). Mix and add 0, 2, 4, 6, 8 and 10 μL of diluted Thrombin Enzyme Standard (2.5 $\text{ng}/\mu\text{L}$) into a series of wells in a 96 well plate. Adjust the volume to 50 μL with Thrombin Assay Buffer to prepare 0, 5, 10, 15, 20 and 25 ng/well of Thrombin Enzyme Standard.

Note: Store the diluted Thrombin Enzyme Standard solution at $-80\text{ }^{\circ}\text{C}$.

Substrate Mix

Prepare enough reagents for the number of assays to be performed. Prepare 50 μL of Substrate Mix for each standard and sample well, see Table 1.

Table 1.

Preparation of Substrate Mix

Reagents	Volume
Thrombin Assay Buffer	45 μL
Thrombin Substrate	5 μL

Mix and add 50 μL of Thrombin Substrate Mix into each standard and sample well, and mix well.

Measurement

Measure fluorescence in kinetic mode for 30–60 minutes at $37\text{ }^{\circ}\text{C}$ ($\lambda_{\text{ex}} = 350\text{ nm}/\lambda_{\text{em}} = 450\text{ nm}$). Choose two time points (T_1 & T_2) in the linear range of the plot and obtain the corresponding values for the fluorescence (RFU_1 and RFU_2).

ResultsCalculations

Apply sample's ΔRFU to to obtain corresponding

Subtract 0 Standard reading from all readings. Plot the Thrombin Standard Curve. Apply the ΔRFU of the sample to the Thrombin Standard Curve to obtain corresponding Thrombin (B, in ng) and calculate the activity of Thrombin in the sample as:

$$\text{Fxa Activity} = \frac{\text{B}}{\text{V}} \times \text{dilution factor}$$

(ng/mL) V
($\mu\text{g}/\text{L}$)

B = Thrombin amount from Standard Curve (ng)

V = sample volume added into the reaction well (mL)

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	White plates are recommended for this assay.
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 samples will be used 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 expired kit or improperly stored reagents	Check the expiration date and store the components appropriately
	Allowing the reagents to sit for extended times on ice	Prepare fresh 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 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

SJ,MAM 04/16-1