

Product Information

Monoamine Oxidase Assay Kit

Catalog Number **MAK136**
Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

Monoamine oxidases (MAO, EC 1.4.3.4) are a family of mitochondrial enzymes that catalyze the oxidative deamination of monoamines. In mammals, MAO consists of two isoenzymes, MAO-A and MAO-B. MAO dysfunction is thought to be responsible for a number of neurological disorders. Unusually high or low levels of MAO in the body have been associated with depression, schizophrenia, substance abuse, attention deficit disorder, migraines, and irregular sexual maturation. MAO inhibitors are one of the major classes of drug prescribed for the treatment of depression.

The Monoamine Oxidase Assay Kit provides a convenient fluorimetric means to measure MAO enzyme activity in biological samples. In the assay, MAO reacts with *p*-tyramine, a substrate for both MAO-A and MAO-B, resulting in the formation of H_2O_2 , which is determined by a fluorimetric method ($\lambda_{\text{ex}} = 530/\lambda_{\text{em}} = 585\text{ nm}$). The assay is simple, sensitive, stable, and high-throughput adaptable. Unit definition: one unit of MAO catalyzes the formation of $1\text{ }\mu\text{mole}$ of H_2O_2 per minute under the assay conditions.

Components

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

| | |
|--|------------------|
| Assay Buffer, pH 7.4 Catalog Number MAK136A | 12 mL |
| Pargyline, 20 mM Catalog Number MAK136B | 50 μL |
| Clorgyline, 20 mM Catalog Number MAK136C | 50 μL |

| | |
|--|-------------------|
| Hydrogen Peroxide, 3% H_2O_2 Catalog Number MAK136D | 100 μL |
| <i>p</i> -Tyramine Catalog Number MAK136E | 120 μL |
| HRP Enzyme Catalog Number MAK136F | 120 μL |
| Dye Reagent Catalog Number MAK136G | 120 μL |

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.
- 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

Use ultrapure water for the preparation of reagents. Equilibrate all components to room temperature before use. Briefly centrifuge vials before opening. Keep thawed tubes on ice during assay.

Storage/Stability

The kit is shipped on dry ice. If desired, store Assay Buffer and Hydrogen Peroxide at $2\text{--}8\text{ }^{\circ}\text{C}$ and all other components at $-20\text{ }^{\circ}\text{C}$.

Procedure

Notes: Thiols (β -mercaptoethanol, dithioerythritol, etc.) $>10 \mu\text{M}$ will interfere with this assay and should be avoided in sample preparation.

Samples should be free of particle or precipitates. MAO can be extracted from a tissue by homogenization and differential centrifugation.¹ Store sample at $-80 \text{ }^\circ\text{C}$.

Prior to assay, concentrations of protein, inhibitor, substrate, and incubation time may need to be established for a given sample.

Inhibitor Preparation

Add $5 \mu\text{L}$ of 20 mM inhibitor to 10 mL water to generate a $10 \mu\text{M}$ solution. Clorgyline is an MAO-A inhibitor and pargyline is an MAO-B inhibitor.

Sample Preparation

1. Dilute samples as needed in Assay Buffer. Aliquot $45 \mu\text{L}$ of each sample into two separate wells, one well will serve as a sample well and one will serve as a sample control well.
2. Add $5 \mu\text{L}$ of water to the sample well and add $5 \mu\text{L}$ of the $10 \mu\text{M}$ inhibitor to the control well. Use clorgyline when assaying for MAO-A and use pargyline when assaying for MAO-B.
3. Mix well using horizontal shaker or by pipetting, and incubate at room temperature for 10 minutes.

For unknown samples, it is suggested to test several sample dilutions.

Hydrogen Peroxide Standards Preparation

Add $5 \mu\text{L}$ of $3\% \text{ H}_2\text{O}_2$ to $1,375 \mu\text{L}$ of water. Further dilute $5 \mu\text{L}$ of the diluted H_2O_2 solution with $795 \mu\text{L}$ of water to generate a $20 \mu\text{M} \text{ H}_2\text{O}_2$ standard solution. Add $0, 12.5, 25,$ and $50 \mu\text{L}$ of $20 \mu\text{M} \text{ H}_2\text{O}_2$ standard solution into separate wells of the 96 well plate and bring volume to $50 \mu\text{L}$ with water generating $0, 5, 10,$ and $20 \mu\text{M} \text{ H}_2\text{O}_2$ standards.

Assay Reaction

1. Set up the Master Reaction Mix according to the scheme in Table 1. $50 \mu\text{L}$ of the Master Reaction Mix is required for each sample and standard reaction (well).

Table 1.
Master Reaction Mix

| Reagent | Sample and Standards |
|--------------------|----------------------|
| Assay Buffer | $50 \mu\text{L}$ |
| <i>p</i> -Tyramine | $1 \mu\text{L}$ |
| HRP Enzyme | $1 \mu\text{L}$ |
| Dye Reagent | $1 \mu\text{L}$ |

2. Add $50 \mu\text{L}$ of the Master Reaction Mix to each of the sample, sample control, and standard wells. Mix well using a horizontal shaker or by pipetting, and incubate the reaction 20 minutes at room temperature. Protect the plate from light during the incubation.
3. Measure the fluorescence (FLU) of the samples and standards at $\lambda_{\text{ex}} = 530/\lambda_{\text{em}} = 585$.

Note: To screen for MAO inhibitors or characterize inhibitor potency (IC_{50}), mix $5 \mu\text{L}$ of $10 \mu\text{M}$ inhibitor solution with $45 \mu\text{L}$ of sample and incubate for at least 10 minutes to allow the inhibitor to interact with the enzyme, prior to adding the Working Reagent.

Results

Plot H₂O₂ standard curve and determine the slope (μM⁻¹).

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

MAO enzyme activity in the sample is calculated as:

$$\text{MAO Activity (units/L)} = \frac{\text{FLU}_{\text{sample}} - \text{FLU}_{\text{control}}}{\text{Slope} \times t}$$

FLU_{sample} = Absorbance measured in unknown sample

FLU_{control} = Absorbance measured in sample control
(sample in presence of inhibitor)

Slope = Determined from calibration curve (μM⁻¹)

t = incubation time (i.e., 20 minutes)

Unit definition: one unit of MAO catalyzes the formation of 1 μmole of H₂O₂ per minute under the assay conditions.

Reference

Biochem. J., **108**, 95 (1968).

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 fluorescence 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 |
| | 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 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 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 calibrator curve | Pipetting errors in preparation of calibrator | 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 |
| | Calibrator stock is at incorrect concentration | Refer to the calibrator 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 |

KVG,MF,LS,MAM 05/16-1