



Promega

# Technical Bulletin

---

## MAO-Glo™ Assay

INSTRUCTIONS FOR USE OF PRODUCTS  
V1401, V1402, V1452 AND V1560.



# MAO-Glo™ Assay

All technical literature is available at: [www.promega.com/tbs/](http://www.promega.com/tbs/)  
Please visit the web site to verify that you are using the most current version of this  
Technical Bulletin. Please contact Promega Technical Services if you have questions on use  
of this system. E-mail: [techserv@promega.com](mailto:techserv@promega.com)

1. Description.....	1
2. Product Components and Storage Conditions .....	3
3. Performing the MAO-Glo™ Assay .....	4
A. General Considerations .....	4
B. Reagent Preparation.....	4
C. Protocol .....	7
4. Results.....	8
5. Composition of Buffers and Solutions .....	9
6. References .....	9
7. Appendix .....	10
A. Monoamine Oxidase Reaction Chemistry .....	10
B. Luciferin Detection Chemistry .....	14
C. Measuring the Affinity of Known Substrates and Inhibitors.....	15
D. Performing the MAO-Glo™ Assay with Promega MAO-A Enzyme .....	17
8. Related Products .....	18

## 1. Description

The MAO-Glo™ Assay<sup>(a-c)</sup> provides a homogeneous luminescent method for measuring monoamine oxidase (MAO) activity from recombinant and native sources and measuring the effects of test compounds on MAO activities (1). The MAO-Glo™ Assay is performed by incubating the MAO enzyme with a luminogenic MAO substrate. The substrate of the MAO-Glo™ Assay is a derivative of beetle luciferin ((4S)-4,5-dihydro-2-(6-hydroxybenzothiazolyl)-4-thiazolecarboxylic acid). MAO converts this luciferin derivative to methyl ester luciferin.

After the MAO reaction has been performed, the reconstituted Luciferin Detection Reagent is added to simultaneously stop the MAO reaction, convert the methyl ester derivative to luciferin and produce light (Figure 1). This addition initiates a stable glow-type luminescent signal with a half-life greater than 5 hours. The amount of light produced is directly proportional to the activity of MAO.



**Figure 1. Conversion of the MAO Substrate by MAO.** Monoamine oxidase acts on the luminogenic MAO Substrate to produce methyl ester luciferin, which is then converted into light by the esterase and luciferase enzymes in the reconstituted Luciferin Detection Reagent. Any effects of test compounds on luciferase or esterase activities have been minimized (Section 7.B).

The MAO-Glo™ Assay includes a luminogenic MAO Substrate, two MAO Reaction Buffers (one that can be used with either MAO-A or MAO-B and one that is designed specifically for MAO-B), a lyophilized Luciferin Detection Reagent and the Reconstitution Buffer with esterase. The user supplies the sample material containing MAO. MAO-A Enzyme is supplied with Cat.# V1560, or may be purchased separately. Protocols are configured for multiwell plate formats but can easily be adapted for single-tube applications.

The MAO-Glo™ Assay is ideal for many applications, including:

- Measuring native and recombinant MAO activities
- Screening drugs and chemical entities for their capacity to modulate MAO activities in native or recombinant fractions

Advantages of the MAO-Glo™ Assays include:

**Speed:** The luminescence format eliminates the need for time-consuming analyses such as HPLC.

**Simplified Method:** The simple “add and read” protocol makes the assay amenable to high-throughput screening in multiwell plates.

**Greater Sensitivity:** Less MAO enzyme is required in these assays than in typical HPLC or fluorometric methods because of the enhanced sensitivity.

**No Fluorescence Interference:** By using luminescence to monitor enzyme activity, any overlaps between the fluorescent excitation and emission wavelengths of reagents and test compounds are inconsequential. In fluorescent assays, such overlaps can confound analysis and present misleading or irrelevant data.

**Signal Stability:** Glow-type luminescence provides a stable signal with a half-life of greater than 5 hours. This eliminates the need for strictly timed luminescent detection.

**Greater Specificity:** These assays are more specific for MAO enzyme than assays that measure  $\text{H}_2\text{O}_2$ .

**Non-Radioactive Detection:** Scintillation cocktail or disposal of radioactive waste is not required.

## 2. Product Components and Storage Conditions

Product	Size	Cat.#
MAO-Glo™ Assay	200 assays	V1401

The kit contains sufficient reagents for 200 assays at 50µl per assay in 96-well plates or 330 assays at 30µl per assay in 384-well plates. Includes:

- 1 vial MAO Substrate
- 10ml MAO Reaction Buffer
- 10ml MAO B Reaction Buffer
- 10ml Reconstitution Buffer with esterase
- 1 vial Luciferin Detection Reagent

Product	Size	Cat.#
MAO-Glo™ Assay	1,000 assays	V1402

The kit contains sufficient reagents for 1,000 assays at 50µl per assay in 96-well plates or 1,660 assays at 30µl per assay in 384-well plates. Includes:

- 1 vial MAO Substrate
- 50ml MAO Reaction Buffer
- 50ml MAO B Reaction Buffer
- 50ml Reconstitution Buffer with esterase
- 1 vial Luciferin Detection Reagent

Product	Size	Cat.#
MAO-Glo™ Assay with MAO-A	1,000 assays	V1560

The kit contains sufficient reagents for 1,000 assays at 50µl per assay in 96-well plates or 1,660 assays at 30µl per assay in 384-well plates. Includes:

- 1 vial MAO Substrate
- 50ml MAO Reaction Buffer
- 50ml MAO B Reaction Buffer
- 50ml Reconstitution Buffer with esterase
- 1 vial Luciferin Detection Reagent
- 500µl MAO-A

Product	Size	Cat.#
MAO-A Enzyme	500µl	V1452

**Note:** MAO-A included in V1560 and V1452 is human recombinant enzyme expressed in yeast.

**Storage Conditions:** Store the MAO A Enzyme at -70°C. Store all other components at -20°C protected from light. See expiration date on the product label. Avoid multiple freeze-thaw cycles of all components.

The resuspended MAO Substrate can be stored at -20°C for up to 3 months. The reconstituted Luciferin Detection Reagent can be stored at room temperature (approximately 22°C) for 1 week or -20°C for 2 months with no change in activity. The activity of the reconstituted Luciferin Detection Reagent decreased approximately 10% when the reagent was stored at 4°C for 2 months.

### 3. Performing the MAO-Glo™ Assay


#### 3.A. General Considerations


MAO-Glo™ Assays are performed in two steps (Figure 2).

**Step 1. The MAO Reaction:** The first step of the MAO-Glo™ Assay combines the MAO Substrate with MAO enzyme to generate methyl ester luciferin. The recommended final substrate concentrations are the apparent  $K_m$  values for the respective reactions, and the recommended amount of MAO should give strong signals and high signal-to-background ratios. However, the magnitude of the MAO-Glo™ Assay signal will vary with the specific activity of the MAO preparation. MAO-Glo™ Assays are performed in MAO Reaction Buffer; MAO-B reactions may also be performed in MAO-B Reaction Buffer, which contains 10% DMSO to increase the activity of MAO-B by approximately fourfold (Section 7.A).

**Step 2. Luciferin Detection:** The methyl ester luciferin, which is produced in Step 1 by the action of MAO on the MAO Substrate, reacts with esterase and luciferase to produce light in Step 2. The MAO-Glo™ Assays use a proprietary stabilized luciferase (Ultra-Glo™ Luciferase) and a proprietary detection reagent formulation to produce a stable “glow-type” luminescent signal with a half-life greater than 5 hours (extrapolated from Figure 9, Section 7.B).

#### 3.B. Reagent Preparation

 The resuspended MAO Substrate can be stored at  $-20^{\circ}\text{C}$  for up to 3 months. The reconstituted Luciferin Detection Reagent can be stored at room temperature (approximately  $22^{\circ}\text{C}$ ) for 1 week or  $-20^{\circ}\text{C}$  for 2 months with no change in activity. The activity of the reconstituted Luciferin Detection Reagent decreased approximately 10% when the reagent was stored at  $4^{\circ}\text{C}$  for 2 months. Avoid multiple freeze-thaw cycles of all components.

 Use the same reaction buffer, either MAO Reaction Buffer or MAO B Reaction Buffer, for each dilution within a single experiment.

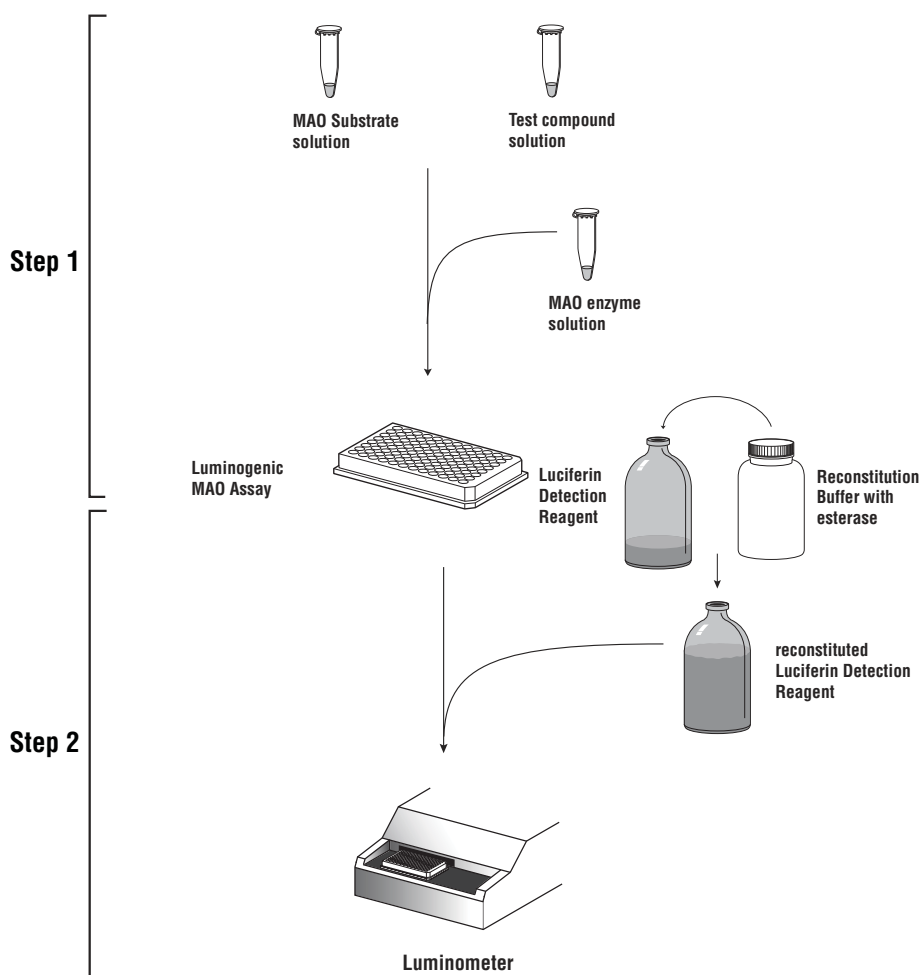
**Note:** The volumes given here are intended for 96-well plates. For 384-well plates, we recommend using 60% of the volumes indicated.

#### Materials to Be Supplied by the User

- MAO enzyme preparation
- test compound
- dimethyl sulfoxide (DMSO)

#### Preparation of the 2X MAO Enzyme Solution

For each reaction, prepare  $26.25\mu\text{l}$  of a 2X MAO enzyme preparation. This volume allows 5% extra material to compensate for pipetting error. Dilute the MAO-containing material with MAO Reaction Buffer to achieve the desired concentration.



5-6080A

**Figure 2. Schematic diagram showing reagent preparation and use of the MAO-Glo™ Assays.**

### 3.B. Reagent Preparation (continued)

#### Preparation of the 4X Test Compound

Prepare a 4X solution of the compound to be tested. Dilute or dissolve the test compound in MAO Reaction Buffer.

#### Preparation of the MAO Substrate Solution

1. Dissolve the dried MAO Substrate in DMSO. For Cat.# V1401, dissolve the substrate in 105 $\mu$ l of DMSO, for Cat.# V1402 and V1560, dissolve the substrate in 525 $\mu$ l of DMSO. This will create a 4mM stock solution.
2. For each reaction, prepare 13.13 $\mu$ l of a 4X MAO Substrate solution. This volume allows 5% extra material to compensate for pipetting error. Dilute the resuspended MAO Substrate from Step 1 with MAO Reaction Buffer to a concentration of 160 $\mu$ M (25-fold dilution) to assay MAO-A or 16 $\mu$ M (250-fold dilution) to assay MAO-B.

For example, to perform a 96-well plate of reactions for MAO-A, use 50 $\mu$ l of MAO Substrate diluted with MAO Reaction Buffer to a final volume of 1,250 $\mu$ l; if assaying MAO-B, use 5.0 $\mu$ l of MAO Substrate diluted to a final volume of 1,250 $\mu$ l.

#### Preparation of the Reconstituted Luciferin Detection Reagent

1. Equilibrate the Reconstitution Buffer with esterase and lyophilized Luciferin Detection Reagent to room temperature.
2. Transfer the entire contents of the bottle of Reconstitution Buffer with esterase to the amber bottle of lyophilized Luciferin Detection Reagent. Mix by swirling or inverting several times to obtain a homogeneous solution. Do not vortex.

---

### 3.C. Protocol

1. Add 12.5µl of 4X MAO Substrate solution per well.
2. Add 12.5µl of 4X test compound per well. If the assay is to be performed without a test compound, add 12.5µl of MAO Reaction Buffer containing the appropriate carrier or solvent.
3. To initiate the MAO reaction, add 25µl of 2X MAO enzyme solution per well, and mix briefly. For negative control reactions, add 25µl of MAO Reaction Buffer.

**Note:** If you are using the MAO-A enzyme provided by Promega (Cat.# V1452), use 120µU of enzyme in a volume of 25µl.

4. Incubate the plate at room temperature for 1 hour.  
**Note:** The net luminescent signal of the MAO-Glo™ Assay depends on both time and temperature (Section 7.A).
5. Add 50µl of reconstituted Luciferin Detection Reagent per well, and mix briefly.
6. Incubate the plate at room temperature for 20 minutes to generate and stabilize the luminescent signal.
7. Measure and record the luminescent signal using a plate-reading luminometer or CCD camera. Values are displayed as relative light units (RLU).

**Note:** Luminometer settings will depend on the manufacturer. Use an integration time of 0.25–1 second per well as a guideline. Relative light units are arbitrary units that vary, depending on the instrument manufacturer and model. Absolute readings from one luminometer may not match those from a different model.

#### 4. Results

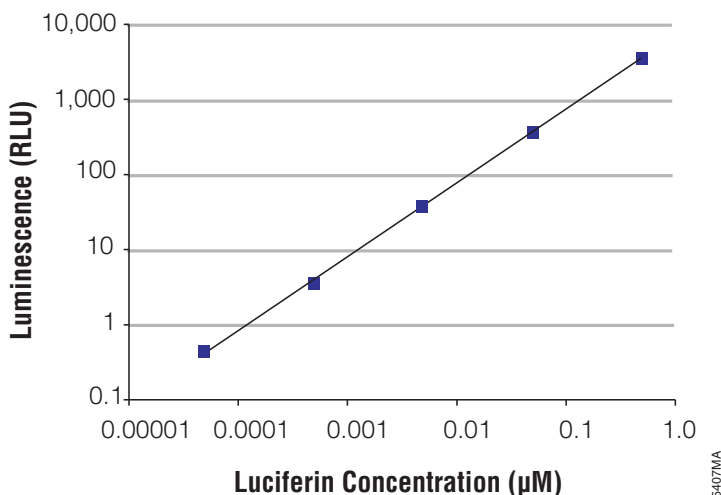
Calculate net MAO-dependent luminescence (net RLU) by subtracting the average luminescence of the negative control reactions without MAO enzyme from that of the MAO-containing reactions. The net signal from MAO reactions in the absence of test compound represents the total MAO activity. Changes from the average net signal for total MAO activity to the net signals for reactions with test compound reflect the effect of the compound on MAO activity. Changes in luminescent signal will typically be seen as decreases because of the inhibition of MAO or use of alternative substrates. However, some compounds may cause signals to increase if they activate the MAO enzyme.

The product of MAO and esterase enzyme activities is D-luciferin. D-luciferin reacts with luciferase to produce light that is directly proportional to the D-luciferin concentration (Figure 3). For example, an inhibitor concentration that causes a 50% drop in luminescence has also caused a 50% decrease in D-luciferin accumulation. Data may be expressed directly in relative light units (RLU) or percent activity, or the values may be converted to reaction rates by dividing RLU by reaction time and amount of enzyme. For example, obtaining 100,000 RLU from a 20-minute reaction with 0.5pmol of MAO-A corresponds to a specific activity of 10,000 RLU/pmol MAO A/minute. Alternatively, RLU may be converted to a corresponding D-luciferin concentration for assays that include a D-luciferin standard curve. However, to detect MAO inhibition at a single concentration of a test compound or to measure the  $IC_{50}$ ,  $K_i$  or  $K_m$  value for an inhibitor or substrate, it is not necessary to convert luminescence to D-luciferin concentration.

**Note:**  $IC_{50}$  refers to the concentration of a compound that inhibits a reaction by 50%. In the case of competitive inhibition,  $IC_{50} = 2K_i$  when the substrate is present at the  $K_m$  concentration, as per the relationship:

$$K_i = IC_{50} / [1 + (\text{substrate concentration} / K_m)].$$

If  $IC_{50}$  values obtained with the MAO-Glo™ Assays will be compared to  $IC_{50}$  or  $K_i$  values obtained with other MAO assays, you must consider this relationship and note that direct comparisons should only be made when the respective substrates are present at their  $K_m$  concentrations.



**Figure 3. The correlation between luminescent signal and D-luciferin concentration.** Luciferin was titrated in MAO Reaction Buffer and mixed with an equal volume of reconstituted Luciferin Detection Reagent. The luminescent signal was measured 20 minutes after reagent addition. For all samples  $n = 4$ , and the standard deviation is smaller than the data points shown.

## 5. Composition of Buffers and Solutions

### MAO Reaction Buffer

100mM HEPES (pH 7.5)  
5% glycerol

### MAO B Reaction Buffer

100mM HEPES (pH 7.5)  
5% glycerol  
10% dimethyl sulfoxide

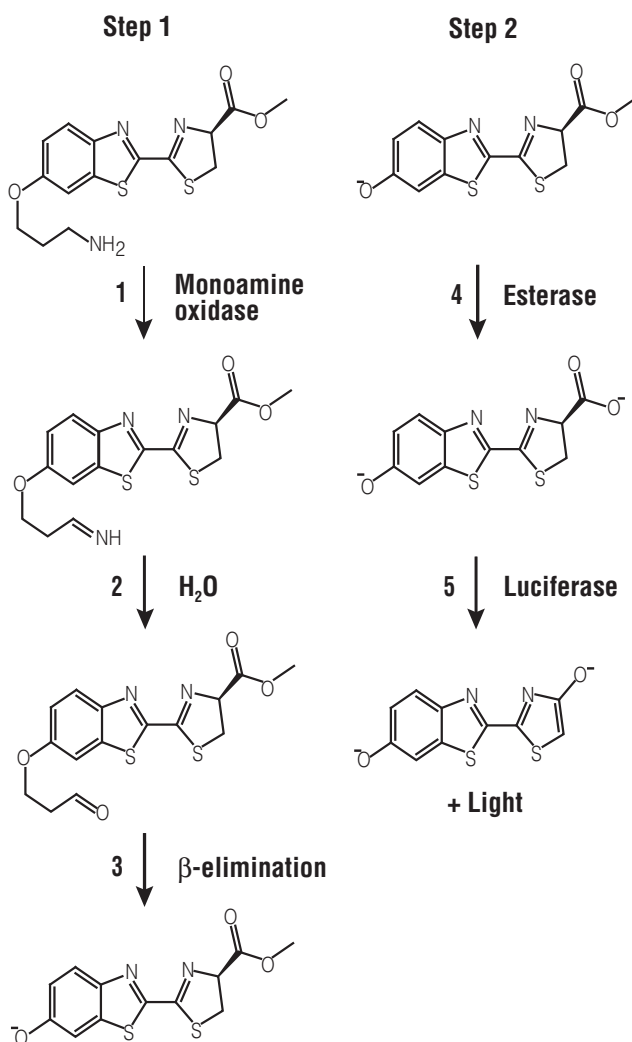
## 6. References

1. Valley, M., *et al.* (2006) A bioluminescent assay for monoamine oxidase activity. *Anal Biochem.* **359**, 238–46.
2. Johnston, J.P. (1968) Some observations upon a new inhibitor of monoamine oxidase in brain tissue. *Biochem. Pharmacol.* **17**, 1285–97.
3. Tsugeno, Y. *et al.* (1995) Regions of the molecule responsible for substrate specificity of monoamine oxidase A and B: A chimeric enzyme analysis. *J. Biochem.* **118**, 974–80.
4. Schoepp, D.D. and Azzaro, A.J. (1981) Specificity of endogenous substrates for types A and B monoamine oxidase in rat striatum. *J. Neurochem.* **36**, 2025–31.
5. Tipton, K.F., Fowler, C.J. and Houslay, M.D. (1982) In: *Monoamine oxidase: Basic and clinical frontiers*, Kamijo, K., Usdin, E. and Nagatsu, T., eds., *Excerpta Medica*, Amsterdam, 87–99.
6. Geha, R.M. *et al.* (2001) Substrate and inhibitor specificities for human monoamine oxidase A and B are influenced by a single amino acid. *J. Biol. Chem.* **276**, 9877–82.

## 7. Appendix

### 7.A. Monoamine Oxidase Reaction Chemistry

Monoamine oxidase (MAO) catalyzes the oxidative deamination of amine-containing substrates. In the MAO-Glo™ Assay, an aminopropylether analog of methyl ester luciferin is provided as a substrate for the MAO reaction. In Step 1 (Figure 4, reactions 1-3), the MAO enzyme oxidizes the amine to an imine, which is subsequently hydrolyzed by water to the corresponding aldehyde. The aldehyde then spontaneously undergoes a  $\beta$ -elimination reaction to generate methyl ester luciferin. Since the latter two reactions are not rate-limiting, the amount of methyl ester luciferin produced is proportional to the activity of MAO.

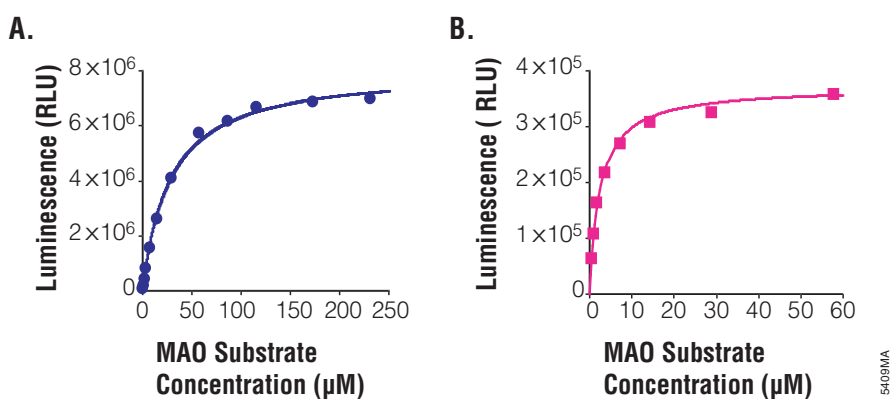


**Figure 4. The chemical and enzymatic reactions required to generate light in the MAO-Glo™ Assay.**

540BMA

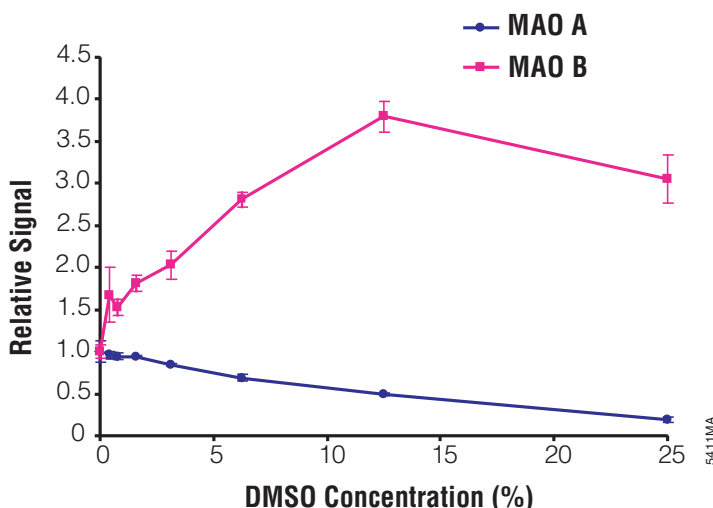
The MAO-A and MAO-B enzymes both use the MAO Substrate, but the two enzymes have different  $K_m$  and  $RLU_{max}$  values (Figure 5). We recommend performing the MAO reactions with the MAO Substrate at their individual  $K_m$  values of  $40\mu M$  and  $4\mu M$  (MAO-A and MAO-B, respectively).

The amount of MAO used and the reaction time will affect the level of luminescence (Figure 6). Increasing the length of the MAO reaction permits the use of less MAO enzyme per reaction. Conversely, increasing the amount of MAO enzyme may permit reaction times as short as 10–20 minutes. We recommend running MAO reactions with  $1\mu g$  of protein from microsomes containing MAO for 1 hour, but the amount of enzyme and time should be optimized for each application. For example, reactions with  $0.2\mu g$ ,  $1\mu g$  or  $5\mu g$  of protein incubated for 3 hours, 40 minutes or 10 minutes, respectively, all give signal-to-noise ratios greater than 100-fold.



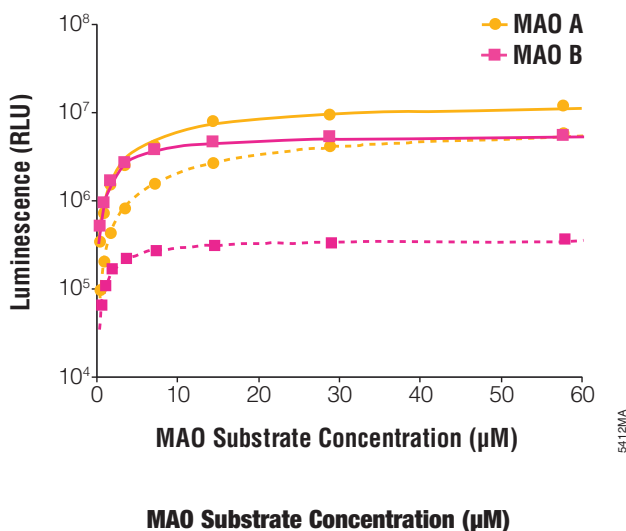
**Figure 5.  $K_m$  values for the MAO Substrate.** MAO-A reactions were performed in MAO Reaction Buffer (**Panel A**), and MAO-B reactions were performed in MAO-B Reaction Buffer (**Panel B**) for 1 hour at  $25^\circ C$ . Microsomes containing the MAO-A or MAO-B enzymes (Sigma-Aldrich, Inc., Cat.# M7316 and M7441, respectively) were used at  $1\mu g$  of protein per  $50\mu l$  reaction. Addition of an equal volume of reconstituted Luciferin Detection Reagent terminated the MAO reaction and generated a luminescent signal, which was read 20 minutes after addition. The data were fit (solid lines) with the program TableCurve (Systat Software, Inc.).

MAO-A and MAO-B are affected differently by environmental conditions. Although solvents such as DMSO are detrimental to the MAO-A reaction, low concentrations of solvent increase the reaction rate of MAO-B (Figure 7) without significantly affecting its  $K_m$  value for the MAO Substrate (data not shown). For this reason, the MAO Reaction Buffer is a general formulation that can be used with either MAO enzyme, whereas the MAO-B Reaction Buffer, which contains 10% DMSO, has been optimized to maximize the luminescent signal in MAO B reactions. Temperature also has a much greater effect on the MAO-B reaction compared to the MAO-A reaction, such that increasing the temperature of the MAO-B reaction can yield a significant increase in the net luminescent signal (Figure 8).



**Figure 7. The effect of DMSO concentration on MAO activity.** The MAO Substrate (40 $\mu$ M or 4 $\mu$ M) was incubated for 1 hour in MAO Reaction Buffer with microsomes containing MAO-A or MAO-B (1 $\mu$ g of protein per 50 $\mu$ l reaction), respectively, and various concentrations of DMSO. The luminescent signal was measured 20 minutes after adding the reconstituted Luciferin Detection Reagent. The relative signal values were normalized to the luminescent signal in the absence of DMSO. For all samples, n = 3, and the standard deviation is illustrated by the error bars shown.

## 7.A. Monoamine Oxidase Reaction Chemistry (continued)

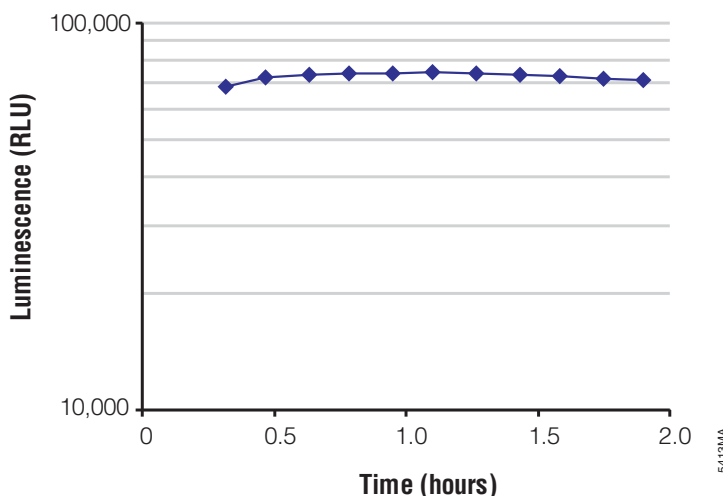


**Figure 8. The effect of temperature on MAO activity.** Reaction conditions are the same as those in Figure 5, except the MAO-A and MAO-B reactions were performed at 25°C (dotted lines) and 35°C (solid lines).

## 7.B. Luciferin Detection Chemistry

In Step 2, the reconstituted Luciferin Detection Reagent inactivates MAO and provides both the enzymes and cosubstrates necessary to convert the product of Step 1 into a luminescent signal. As shown in Figure 4 (reactions 4–5), an esterase converts the methyl ester luciferin into luciferin, which is subsequently consumed by luciferase to yield oxyluciferin and light. Since time is required for the esterase reaction and stabilization of the luciferase signal, the reconstituted Luciferin Detection Reagent should be added to the plates at least 20 minutes before quantifying luminescence (Figure 9). After this period, the luminescent signal is extremely stable, allowing the quantitation of numerous plates of samples at any time over many hours. For the most consistent data, we recommend incorporating a common control on each plate to account for any small variations in luminescence that can occur over time, such as those due to changes in temperature or the instrument signal.

The potential for the luciferase and esterase in the MAO-Glo™ Assays to affect the outcome of an assay has been minimized by maintaining high enzyme concentrations and creating specific chemistries to reduce the effects of potential inhibitors. For example, 10µM of the esterase substrates ethyl butyrate, ethyl acetate and 4-nitrophenyl acetate had little to no effect on the luminescent output of a MAO-B reaction (98.5% + 2.1%; 98.8% + 1.1%; and 98.4% + 1.6% of control samples, respectively; n = 3).



**Figure 9. The luminescent signal is stable in Step 2.** In a 50 $\mu$ l reaction, 40 $\mu$ M MAO Substrate was incubated with 2 $\mu$ g of protein from microsomes containing MAO-A for 45 minutes. After addition of an equal volume of reconstituted Luciferin Detection Reagent, the luminescent signal was read repeatedly over 2 hours. Similar trends were noted in experiments with MAO-B.

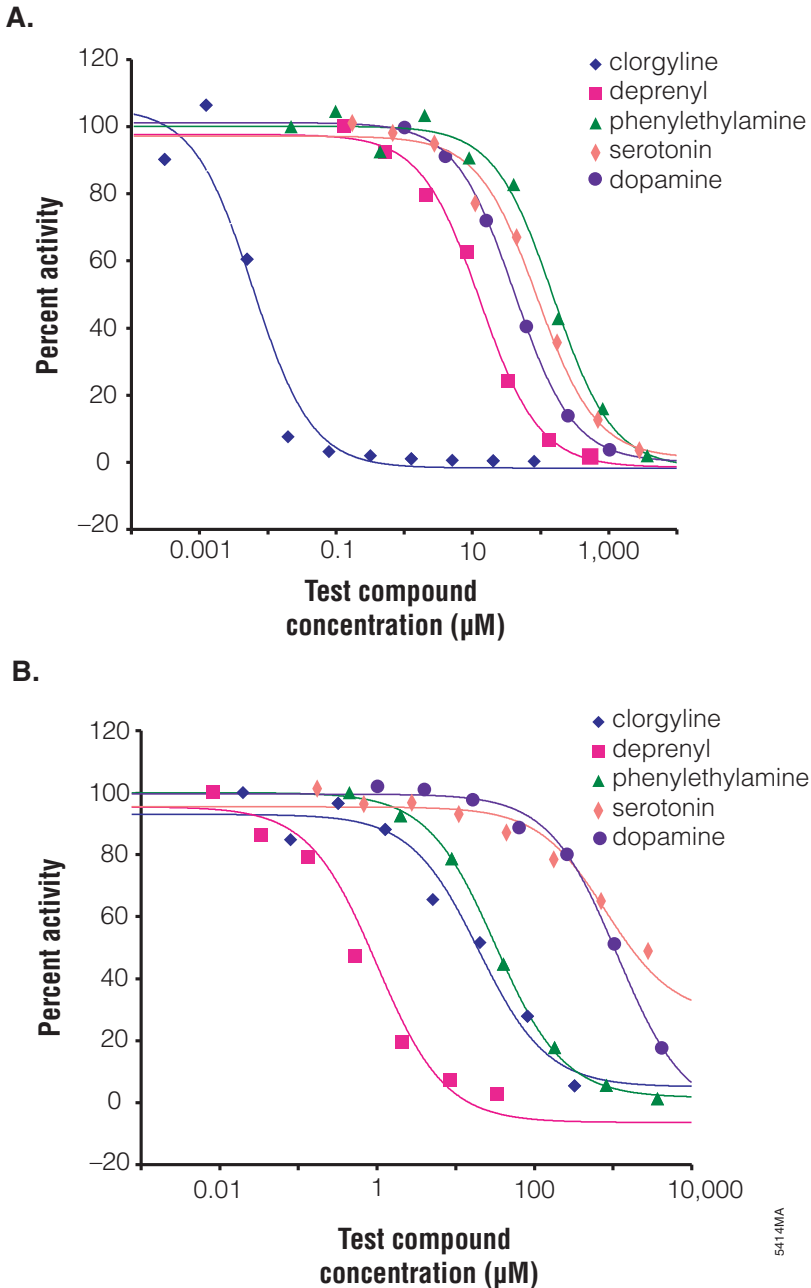
### 7.C. Measuring the Affinity of Known Substrates and Inhibitors

The ability of the MAO-Glo™ Assay to detect the effect of test compounds on MAO activity was demonstrated by performing the assay in the presence of known substrates or inhibitors of MAO enzymes (Figure 10). Clorgyline and deprenyl are MAO-A- and MAO-B-specific inhibitors, respectively, while phenylethylamine, serotonin and dopamine are substrates with varying specificity for each MAO isozyme. In Table 1, the  $K_m$  or  $K_i$  values obtained with the MAO-Glo™ Assay are compared to those previously published. The values for known substrates or inhibitors shown in Table 1 were calculated from the data shown in Figure 10.

	MAO-A		MAO-B	
	$K_i$ or $K_m$ value ( $\mu$ M)	published value ( $\mu$ M)	$K_i$ or $K_m$ value ( $\mu$ M)	published value ( $\mu$ M)
clorgyline	0.003 $\pm$ 0.001	0.0039 <sup>1</sup>	10 $\pm$ 4	4 <sup>1</sup>
deprenyl	7 $\pm$ 1	5 <sup>2</sup>	0.5 $\pm$ 0.2	0.13 <sup>2</sup>
phenylethylamine	78 $\pm$ 16	78 <sup>3</sup>	16 $\pm$ 1	20 <sup>4</sup>
serotonin	45 $\pm$ 8	80 <sup>5</sup>	410 $\pm$ 140	2032 <sup>4</sup>
dopamine	21 $\pm$ 1	120 <sup>3</sup>	570 $\pm$ 120	301 <sup>3</sup>

**Table 1.  $K_i$  or  $K_m$  Values for Known MAO Substrates or Inhibitors.**

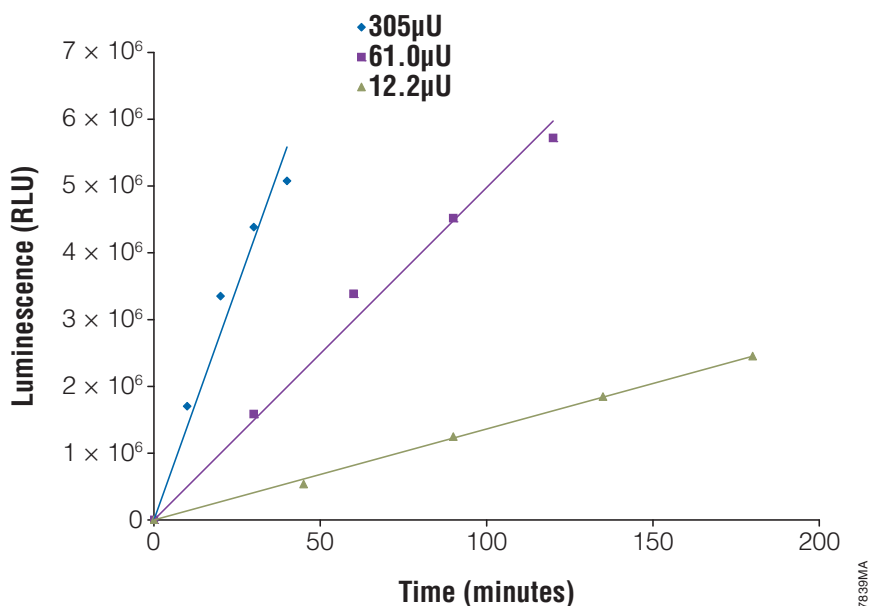
<sup>1</sup>As published by Johnston (2). <sup>2</sup>As published by Tsugeno *et al.* (3). <sup>3</sup>As published by Schoepp and Azzaro. (4). <sup>4</sup>As published by Tipton, Fowler and Houslay (5). <sup>5</sup>As published by Geha *et al.* (6).



**Figure 10. The MAO-Glo™ Assay can measure the binding of known substrates and inhibitors.** MAO-A (Panel A) and MAO-B (Panel B) reactions were performed in the presence of various inhibitors for 1 hour at 25°C, and the luminescent signal was measured 20 minutes after adding the reconstituted Luciferin Detection Reagent. The data was fit (solid lines) with the program TableCurve (Systat Software, Inc.). Test compounds were purchased from Sigma-Aldrich, Inc., and dissolved in MAO Reaction Buffer prior to use.

### 7.D. Performing the MAO-Glo™ Assay with Promega MAO-A Enzyme

If you are using the MAO-A enzyme provided by Promega (Cat.# V1452), use 120 $\mu$ U of enzyme in a volume of 25 $\mu$ l in Step 3 of our suggested protocol (Section 3.C, Step 3). Figure 11 shows the correlation between net signal, amount of enzyme, and the time of the MAO reaction.



**Figure 11. The luminescent signal varies with the amount of enzyme and time in Step 1.** The MAO Substrate (40 $\mu$ M) was incubated for various lengths of time with microsomes containing the Promega MAO-A enzyme. Each 50 $\mu$ l reaction contained 305, 61.0 or 12.2  $\mu$ U of enzyme. The luminescent signal was measured 20 minutes after adding the reconstituted Luciferin Detection Reagent. For all samples, n = 3, and the standard deviation is smaller than the data points shown.

## 8. Related Products

### Cytochrome P450 Assays

Product	Size	Cat.#
P450-Glo™ CYP1A1 Assay	10ml	V8751
	50ml	V8752
P450-Glo™ CYP1B1 Assay	10ml	V8761
	50ml	V8762
P450-Glo™ CYP1A2 Assay	10ml	V8771
	50ml	V8772
P450-Glo™ CYP2C8 Assay	10ml	V8781
	50ml	V8782
P450-Glo™ CYP2C9 Assay	10ml	V8791
	50ml	V8792
P450-Glo™ CYP3A4 Assay	10ml	V8801
	50ml	V8802
P450-Glo™ CYP3A7 Assay	10ml	V8811
	50ml	V8812
P450-Glo™ CYP2C19 Assay	10ml	V8881
	50ml	V8882
P450-Glo™ CYP2D6 Assay	10ml	V8891
	50ml	V8892
P450-Glo™ CYP1A2 Screening System	1,000 assays	V9770
P450-Glo™ CYP2C9 Screening System	1,000 assays	V9790
P450-Glo™ CYP3A4 Screening System	1,000 assays	V9800
P450-Glo™ CYP2C19 Screening System	1,000 assays	V9880
P450-Glo™ CYP2D6 Screening System	1,000 assays	V9890

### Luminogenic Enzyme Substrates

Product	Size	Cat.#
Luciferin-NAT2	3mg	P1721
Luciferin-3A7	3mg	P1741
Luciferin-4A11	3mg	P1621
Luciferin-4F2/3	3mg	P1651
Luciferin-4F12	3mg	P1661
Luciferin-2J2/4F12 ester	3mg	P1671
Luciferin-MultiCYP ester	3mg	P1731
Luciferin Detection Reagent	50ml	V8921
Luciferase Detection Reagent with Esterase	50ml	V8931

Additional sizes available. **Note:** Use Cat.# V8921 with Cat.# P1621, P1651, P1661, P1721, and P1741. Use Cat.# V8931 with Cat.# P1671 and P1731.

<b>Product</b>	<b>Size</b>	<b>Cat.#</b>
Beetle Luciferin, Potassium Salt	5mg	E1601
	50mg	E1602
	250mg	E1603
	1g	E1605

### P-Glycoprotein Assays

<b>Product</b>	<b>Size</b>	<b>Cat.#</b>
Pgp-Glo™ Assay System	10ml	V3591
Pgp-Glo™ Assay System with P-glycoprotein	10ml	V3601

### Cell Viability and Apoptosis Assays

<b>Product</b>	<b>Size</b>	<b>Cat.#</b>
CellTiter-Glo® Luminescent Cell Viability Assay	10ml	G7570
	10 × 10ml	G7571
	100ml	G7572
	10 × 100ml	G7573
CellTiter 96® AQueous One Solution Cell Proliferation Assay	200 assays	G3582
	1,000 assays	G3580
	5,000 assays	G3581
CellTiter-Blue® Cell Viability Assay	20ml	G8080
	100ml	G8081
	10 × 100ml	G8082
Caspase-Glo® 3/7 Assay	2.5ml	G8090
	10ml	G8091
	100ml	G8092
	10 × 100ml	G8093
Caspase-Glo® 8 Assay	2.5ml	G8200
	10ml	G8201
	100ml	G8202
Caspase-Glo® 9 Assay	2.5ml	G8210
	10ml	G8211
	100ml	G8212
Apo-ONE® Homogeneous Caspase-3/7 Assay	10ml	G7790
	100ml	G7791

---

## Luminometers

<b>Product</b>	<b>Size</b>	<b>Cat.#</b>
GloMax <sup>®</sup> Multi Base Instrument	1 each	E7031
GloMax <sup>®</sup> Multi Luminescence Module	1 each	E7041
GloMax <sup>®</sup> Multi Fluorescence Module	1 each	E7051
GloMax <sup>®</sup> Multi Absorbance Module	1 each	E7061
GloMax <sup>®</sup> 96 Microplate Luminometer	1 each	E6501
GloMax <sup>®</sup> 20/20 Luminometer	1 each	E5311

<sup>(a)</sup>U.S. Pat. Nos. 6,602,677 and 7,241,584, Australian Pat. Nos. 754312 and 785294, European Pat. No. 1131441 and other patents pending.

<sup>(b)</sup>Patent Pending.

<sup>(c)</sup>The method of recombinant expression of *Coleoptera* luciferase is covered by U.S. Pat. Nos. 5,583,024, 5,674,713 and 5,700,673.

© 2005-2011 Promega Corporation. All Rights Reserved.

Apo-ONE, Caspase-Glo, CellTiter 96, CellTiter-Blue, CellTiter-Glo and GloMax are registered trademarks of Promega Corporation. MAO-Glo, P450-Glo, Pgp-Glo and Ultra-Glo are trademarks of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.