

Data Sheet

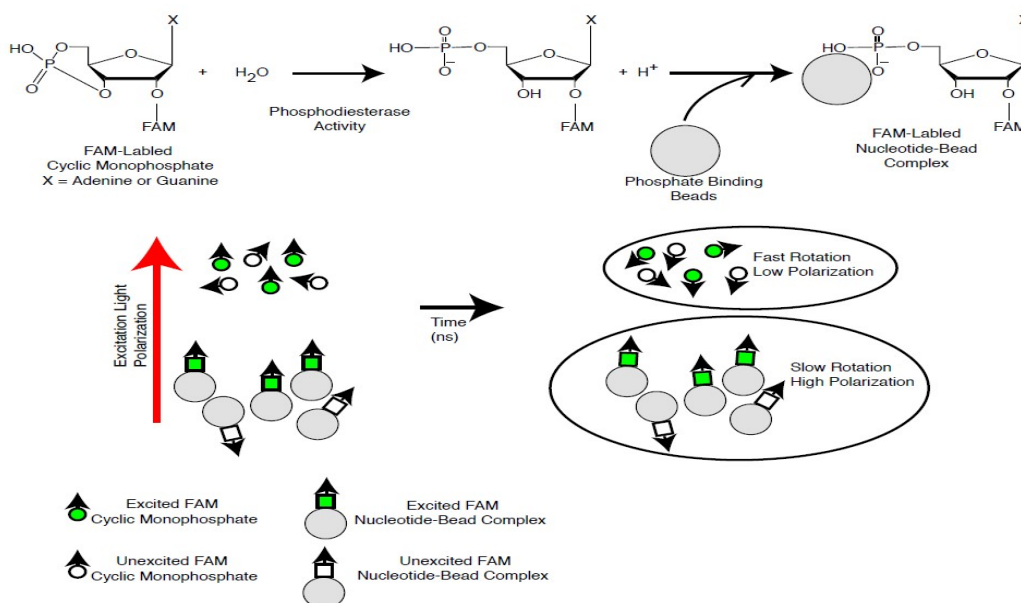
Mouse PDE1C Assay Kit

Catalog # 61312
 Size: 96 reactions

DESCRIPTION: Phosphodiesterases (PDEs) play an important role in the dynamic regulation of cAMP and cAMP signaling. Mouse PDE1C, also known as cAMP-inhibited phosphodiesterase, has been implicated in cardiovascular function and fertility.

The Mouse PDE1C Assay Kit is designed for identification of inhibitors of Mouse PDE1C using fluorescence polarization. The assay is based on the binding of a fluorescent nucleotide monophosphate generated by Mouse PDE1C to the binding agent.

Phosphodiesterases catalyze the hydrolysis of the phosphodiester bond in dye-labeled cyclic monophosphates. Beads selectively bind the phosphate group in the nucleotide product. This increases the size of the nucleotide relative to unreacted cyclic monophosphate. In the polarization assay, dye molecules with absorption transition vectors parallel to the linearly-polarized excitation light are selectively excited. Dyes attached to the rapidly-rotating cyclic monophosphates will obtain random orientations and emit light with low polarization. Dyes attached to the slowly-rotating nucleotide-bead complexes will not have time to reorient and therefore will emit highly polarized light.



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

The *Mouse PDE1C Assay Kit* comes in a convenient 96-well format, with purified Mouse PDE1C enzyme, fluorescently labeled substrate (cAMP), binding agent, and PDE assay buffer for 100 enzyme reactions. The key to the *Mouse PDE1C Assay Kit* is the specific binding agent. Using this kit, only two simple steps on a microtiter plate are required for Mouse PDE1C reactions. First, the fluorescently labeled cAMP is incubated with a sample containing Mouse PDE1C for 1 hour. Second, a binding agent is added to the reaction mix to produce a change in fluorescent polarization that can then be measured using a fluorescence reader equipped for the measurement of fluorescence polarization.

COMPONENTS:

| Catalog # | Component | Amount | Storage | |
|-----------|--------------------------------------|--------|------------|------------------------------------|
| 60012 | Mouse PDE1C recombinant enzyme | >1 µg | -80°C | (Avoid freeze/thaw cycles!) |
| 60200 | FAM-Cyclic-3', 5'-AMP (20 µM) | 50 µl | -80°C | |
| 60393 | PDE assay buffer | 25 ml | -20°C | |
| 60390 | Binding Agent | 100 µl | +4°C | |
| 60391 | Binding Agent Diluent (cAMP) | 10 ml | +4°C | |
| 79685 | Black, low binding, microtiter plate | 1 | Room temp. | |

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable to measure fluorescence polarization.
Adjustable micropipettor and sterile tips.
1,4-Dithiothreitol (DTT) 1 M in anhydrous DMSO.

APPLICATIONS: Great for studying enzyme kinetics and screening small molecular inhibitors for drug discovery and HTS applications.

STABILITY: 6 months from date of receipt when stored as directed.

REFERENCES:

1. Knight WE, Chen S, Zhang Y, *et al.* PDE1C deficiency antagonizes pathological cardiac remodeling and dysfunction. (2016). *Proc. Natl. Acad. Sci. USA* **113(45)**:E7116-E7125.
2. Vandeput, F., Wolda, S. L., Krall, J., Hambleton, R., Uher, L., McCaw, K. N., Radwanski, P. B., Florio, V., Movsesian, M. A. (2007). Cyclic nucleotide phosphodiesterase PDE1C1 in human cardiac myocytes. *J. Biol. Chem.* **282**: 32749-32757.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Dilute 20 μ M **FAM-Cyclic-3', 5'-AMP** stock 100-fold with **PDE assay buffer** to make a 200 nM solution. Make only sufficient quantity needed for the assay; store remaining 20 μ M stock solution in aliquots at -20°C .
- 2) Dilute 1M 1,4-Dithiothreitol (DTT) 1:500 into the diluted **FAM-Cyclic-3',5'-AMP**. For example, add 10 μ l DTT (1M) to 5 ml of diluted FAM-Cyclic-3', 5'-GMP (200 nM).
- 3) Add 25 μ l of **FAM-Cyclic-3',5'-AMP** (200 nM) to each well designated "Positive Control," "Test Inhibitor," and "Substrate Control."
- 4) Add 45 μ l of **PDE assay buffer** to each well designated "Blank" and add 20 μ l of **PDE assay buffer** to each well designated "Substrate Control."
- 5) Add 5 μ l of inhibitor solution to each well designated "Test Inhibitor." For the wells labeled "Positive Control," "Substrate Control," and "Blank," add 5 μ l of the same solution without inhibitor (inhibitor buffer).
- 6) Thaw **Mouse PDE1C** on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. Aliquot **Mouse PDE1C** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: Mouse PDE1C is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*

| | Positive Control | Test Inhibitor | Substrate Control | "Blank" Negative Control |
|---------------------------------|-----------------------------|-----------------------------|-----------------------------|-----------------------------|
| FAM-Cyclic-3',5'-AMP (200 nM) | 25 μ l | 25 μ l | 25 μ l | - |
| PDE assay buffer | - | - | 20 μ l | 45 μ l |
| Inhibitor (in PDE assay buffer) | - | 5 μ l | - | - |
| Inhibitor Buffer (no inhibitor) | 5 μ l | - | 5 μ l | 5 μ l |
| Mouse PDE1C (0.5 pg/ μ l) | 20 μ l | 20 μ l | - | - |
| Total | 50 μl | 50 μl | 50 μl | 50 μl |

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

7) Dilute **Mouse PDE1C** in **PDE assay buffer** to 0.5 pg/μl (0.01 ng/reaction)*. Initiate reaction by adding 20 μl of diluted **Mouse PDE1C** (0.5 pg/μl) to the wells designated "Positive Control" and "Test Inhibitor." Discard any remaining diluted enzyme after use.
**Note: Optimal enzyme concentration may vary with the specific activity of the enzyme.*

8) Incubate the plate at room temperature for 1 hour.

Step 2:

- 1) Mix **Binding Agent** thoroughly and dilute **Binding Agent** 1:100 with **Binding Agent Diluent**.
- 2) Add 100 μl of diluted **Binding Agent** to each microwell. Incubate at room temperature for 30 minutes with slow shaking.
- 3) Read the fluorescent polarization of the sample in a microtiter-plate reader equipped for the measurement of fluorescence polarization, capable of excitation at wavelengths ranging from 485 ± 5 nm and detection of emitted light ranging from 528 ± 10 nm. Blank value is subtracted from all other values.

CALCULATING RESULTS:

Definition of Fluorescence Polarization

$$P = \frac{I_{\parallel} - I_{\perp}}{I_{\parallel} + I_{\perp}}$$

where I_{\parallel} = Intensity with polarizers parallel and I_{\perp} = Intensity with polarizers perpendicular.

The equation above assumes that light is transmitted equally well through both parallel and perpendicular oriented polarizers. In practice, this is generally not true and a correction must be made to measure the absolute polarization state of the molecule. This correction factor is called the "G Factor".

$$FP(\text{measured}) = \frac{([I_{\parallel}] - G * [I_{\perp}])}{([I_{\parallel}] + G * [I_{\perp}])} * 1000$$

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



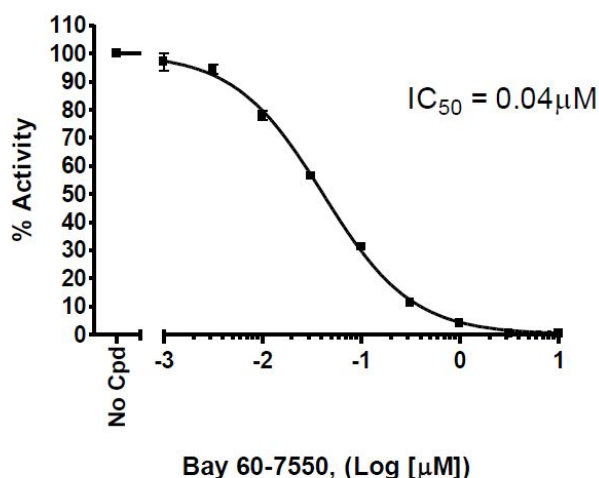
6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

The G-factor is instrument-dependent and may vary slightly depending upon instrument and conditions. Please check the manual of your instrument to obtain the information about the establishment of the G-factor.

EXAMPLE OF ASSAY RESULTS:

Mouse PDE1C Activity

Substrate Conc. = 100nM (cAMP)



Inhibition of murine PDE1C by Bay 60-7550 measured using the *Mouse PDE1C Assay Kit*, BPS Bioscience #61312. Fluorescence polarization was measured at 528 nm using a Tecan M1000 fluorescent microplate reader. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS :

| <u>Product Name</u> | <u>Catalog #</u> | <u>Size</u> |
|-------------------------------------|-------------------------|--------------------|
| Mouse PDE1C | 60012 | 10 µg |
| Mouse PDE3a | 60036 | 5 µg |
| Mouse PDE2a V1 | 60017 | 5 µg |
| Mouse PDE2a V2 | 60018 | 5 µg |
| Mouse PDE2a V4 | 60019 | 5 µg |
| Mouse PDE6C | 60065 | 5 µg |
| Mouse PDE5A | 60051 | 10 µg |
| Mouse PDE7A | 60072 | 10 µg |
| Mouse PDE7B | 60073 | 10 µg |
| Mouse PDE10C | 60101 | 10 µg |
| Mouse PDE11A | 60064 | 5 µg |
| Rat PDE4B | 60049 | 5 µg |
| rRat PDE4D | 60054 | 5 µg |
| Rat PDE1B | 60009 | 10 µg |
| Rat PDE2A | 60022 | 5 µg |
| Rat PDE7A | 60074 | 10 µg |
| Rat PDE7B | 60075 | 10 µg |
| Rat PDE10A | 60102 | 5 µg |
| PDE4B2 | 60042 | 5 µg |
| Mouse PDE3A1 Assay Kit | 79606 | 96 rxns. |
| Mouse PDE5A1 Assay Kit | 79602 | 96 rxns. |
| Mouse PDE5A1 Assay Kit | 79602 | 96 rxns. |
| Dog PDE4B | 60055 | 5 µg |
| Rat PDE4B Assay Kit | 79571 | 96 rxns. |
| PDE Assay Kit | 60300 | 96 rxns. |
| PDE4B2 Assay Kit | 60343 | 96 rxns. |
| PDE4A1A Assay Kit | 60340 | 96 rxns. |
| PDE7A2 Assay Kit | 60345 | 96 rxns. |
| PDE7A Cell-Based Activity Assay Kit | 60505 | 500 rxns. |

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com