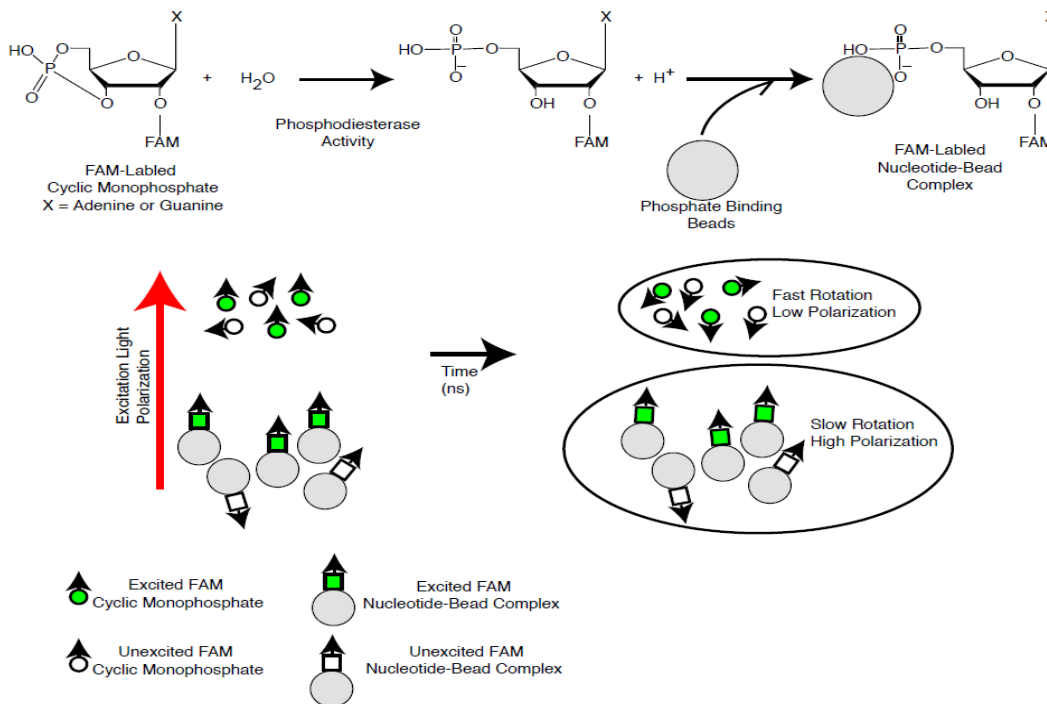


## Data Sheet **PDE10A1 Assay Kit** Catalog # 60385

**BACKGROUND:** Phosphodiesterases (PDEs) play an important role in dynamic regulation of cAMP and cGMP signaling. PDE10A is a dual substrate PDE highly expressed in striatal medium spiny neurons. PDE10A inhibitors can improve the cognitive symptoms of schizophrenia, and exhibit potential therapeutic value for Huntington's disease. PDE10A1 is located in the cytosol, whereas PDE10A2 is a membrane-associated protein.

**DESCRIPTION:** The PDE10A1 Assay Kit is designed for identification of PDE10A1 inhibitors using fluorescence polarization. The assay is based on the binding of a fluorescent nucleotide monophosphate generated by PDE10A1 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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://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 PDE10A1 inhibitor screening assay kit comes in a convenient 384-well format, including purified PDE10A1 enzyme, fluorescently labeled PDE10A1 substrate (cAMP), binding agent, and PDE assay buffer for 400 enzyme reactions. The key to the *PDE10A1 Assay Kit* is the specific binding agent. Using this kit, only two simple steps on a microtiter plate are required for PDE10A1 reactions. First, the fluorescently labeled cAMP is incubated with a sample containing PDE10A1 for 1 hour. Second, a binding agent is added to the reaction mix to produce a change in fluorescent polarization. The FP signal is measured using a fluorescent microplate reader *capable of measuring fluorescence polarization*.

#### COMPONENTS:

Catalog #	Component	Amount	Storage	
60099	PDE10A1 recombinant enzyme	1 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
60200	FAM-Cyclic-3', 5'-AMP: 20 µM	50 µl	-80°C	
60393	PDE assay buffer	25 ml	-20°C	
60390	Binding Agent	250 µl	+4°C	
60391	Binding Agent Diluent (cAMP)	25 ml	+4°C	
	Black, low binding, 384 microtiter plate	1	Room temp.	

#### MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable of measuring fluorescence polarization

**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) Nishi, A., and Snyder, G.L. *J. Pharmacol. Sci.*, 2010, **114**: 6-16.
- 2) Omori, K., and Kotera, J. *Circulation Res.* 2007; **100**: 309-327.

#### ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

##### Step 1:

- 1) Dilute **20 µM FAM-Cyclic-3',5'-AMP** substrate stock solution 100-fold with **PDE Assay Buffer** to make a 200 nM solution. Make only a sufficient quantity needed for the assay; store remaining stock solution in aliquots at -20°C.
- 2) Add 12.5 µl of **FAM-Cyclic-3',5'-AMP** (200 nM) to each well designated "Substrate Control", "Positive Control", and "Test Inhibitor". Add 12.5 µl of PDE assay buffer to each well designated "Blank".

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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

- 3) Add 2.5  $\mu$ l of inhibitor solution to each well designated "Test Inhibitor". Add 2.5  $\mu$ l of the same solution without inhibitor (inhibitor buffer) to the wells labeled "Blank", "Substrate Control" and "Positive Control".
- 4) Add 10  $\mu$ l of **PDE Assay Buffer** to the wells designated as the "Blank" and "Substrate Control".
- 5) Thaw **PDE10A1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. Aliquot **PDE10A1** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: PDE10A1 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 6) Dilute **PDE10A1** in **PDE Assay Buffer** to 1 pg/ $\mu$ l (10 pg/reaction)\*. Initiate reaction by adding 10  $\mu$ l of diluted **PDE10A1** to the wells designated for the "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.*
- 7) Incubate at room temperature for 1 hour.

	Blank	Substrate Control	Positive Control	Test Inhibitor
FAM-Cyclic-3',5'-AMP (200 nM)	-	12.5 $\mu$ l	12.5 $\mu$ l	12.5 $\mu$ l
PDE assay buffer	22.5 $\mu$ l	10 $\mu$ l	-	-
Test Inhibitor	-	-	-	2.5 $\mu$ l
Inhibitor Buffer (no inhibitor)	2.5 $\mu$ l	2.5 $\mu$ l	2.5 $\mu$ l	-
PDE10A1 (1 pg/ $\mu$ l)	-	-	10 $\mu$ l	10 $\mu$ l
<b>Total</b>	<b>25 <math>\mu</math>l</b>	<b>25 <math>\mu</math>l</b>	<b>25 <math>\mu</math>l</b>	<b>25 <math>\mu</math>l</b>

## Step 2:

- 1) Shake the tube containing the **Binding Agent** to ensure it is thoroughly mixed. Mix **binding agent** thoroughly and dilute **binding agent** 1:100 with the cAMP **Binding Agent Diluent**.
- 2) Add 50  $\mu$ l diluted **Binding Agent** to each well. Incubate at room temperature for 20 minutes with slow shaking.
- 3) Read the fluorescent polarization of the sample in a microtiter-plate reader capable of excitation at wavelengths ranging from 485  $\pm$  5 nm and detection of emitted light ranging from 528  $\pm$  5 nm. Blank value is subtracted from all other values.

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

## 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. Most instruments display fluorescence polarization in units of mP.

$$mP = \left( \frac{I_{\parallel} - I_{\perp}}{I_{\parallel} + I_{\perp}} \right) \times 1000$$

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".

$$mP = \left( \frac{I_{\parallel} - G(I_{\perp})}{I_{\parallel} + G(I_{\perp})} \right) \times 1000 \quad \text{OR} \quad mP = \left( \frac{G(I_{\parallel}) - I_{\perp}}{G(I_{\parallel}) + I_{\perp}} \right) \times 1000$$

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.

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

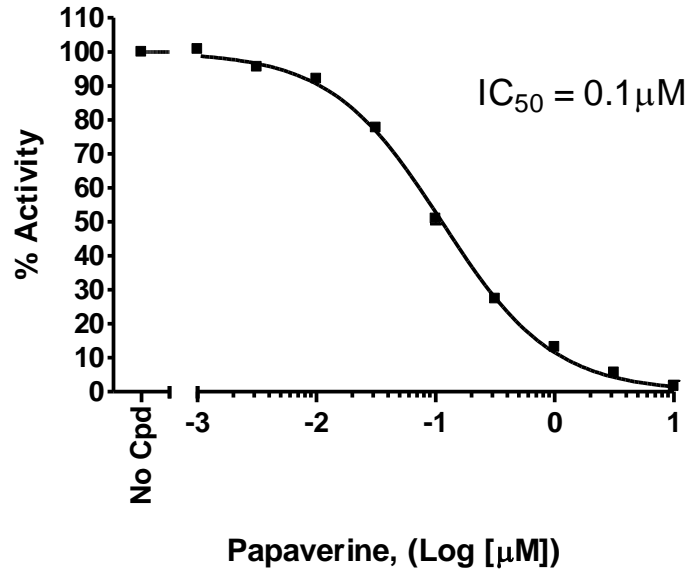


6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**EXAMPLE OF ASSAY RESULTS:**

**PDE10A1 Activity**

Substrate Conc. = 100nM (cAMP)



Inhibition of PDE10A1 by Papaverine, measured using the *PDE10A1 Assay Kit*, BPS Bioscience #60385. 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](mailto: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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS :**

<b><u>Product</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
PDE10A1	60099	10 µg
PDE10A2	60100	10 µg
Mouse PDE10A	60101	10 µg
PDE1A1	60010	10 µg
PDE1B	60011	10 µg
PDE3A	60030	10 µg
hPDE10A-HEK293 Cell Line	60410	2 vials
PDE10A1 TR-FRET Assay Kit	60709	96 rxns.
PDE10A2 TR-FRET Assay Kit	60710	96 rxns.
PDE1A Assay Kit	60310	96 rxns.
PDE1B Assay Kit	60311	96 rxns.
PDE3A Assay Kit	60330	96 rxns.
PDE4A Assay Kit	60340	96 rxns.
PDE5A Assay Kit	60350	96 rxns.
PDE10A Assay Kit	60400	96 rxns.
FAM-cAMP Substrate	60200	100 nmole.

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\*\*](mailto:info@bpsbioscience.com)

Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)