

## Data Sheet

### **Rat PDE4D Assay Kit**

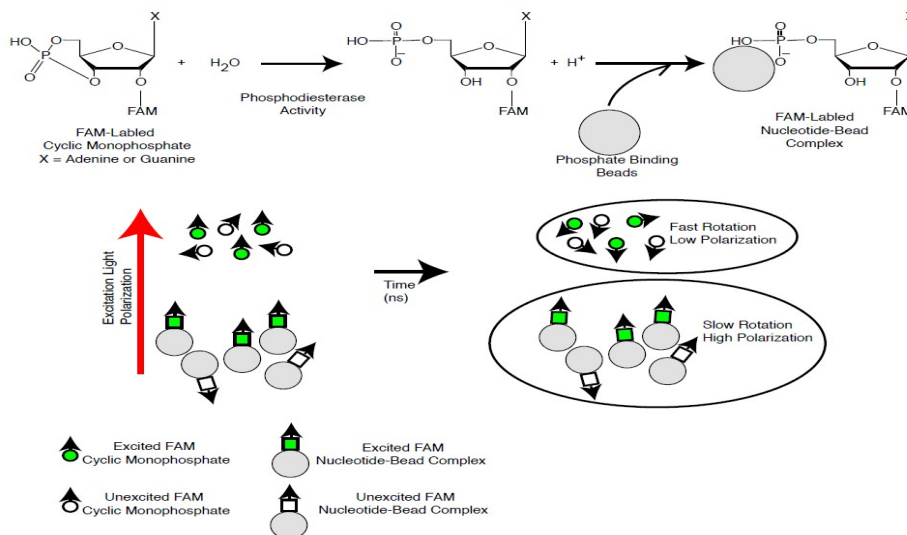
Catalog # 79629  
 Size: 96 reactions

#### DESCRIPTION:

Phosphodiesterase's (PDEs) play an important role in the dynamic regulation of cAMP and cAMP signaling. Rat PDE4D, also known as cAMP-inhibited phosphodiesterase, has been implicated in cardiovascular function and fertility.

The Rat PDE4D Assay Kit is designed for identification of inhibitors of Rat PDE4D using fluorescence polarization. The assay is based on the binding of a fluorescent nucleotide monophosphate generated by Rat PDE4D 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](mailto:info@bpsbioscience.com)

The Rat PDE4D Assay Kit comes in a convenient 96-well format, with purified Rat PDE4D enzyme, fluorescently labeled substrate (cAMP), binding agent, and PDE assay buffer for 100 enzyme reactions. The key to the *Rat PDE4D Assay Kit* is the specific binding agent. Using this kit, only two simple steps on a microtiter plate are required for Rat PDE4D reactions. First, the fluorescently labeled cAMP is incubated with a sample containing Rat PDE4D 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	
60054	Rat PDE4D 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	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)

#### APPLICATIONS:

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

#### STABILITY:

At least 6 months from date of receipt when stored as directed.

#### REFERENCES:

1. Clapcote, S.J. *Adv Neurobiol.* 2017;**17**:103-131
2. Braun, N.N., *et al.*, *NeuroReport.* 2007; Nov; **18**(17):1841-1844.

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)

## ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

### Step 1:

- 1) Dilute 20  $\mu$ 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  $\mu$ 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  $\mu$ l DTT (1M) to 5 ml of diluted FAM-Cyclic-3', 5'-GMP (200 nM).
- 3) Add 25  $\mu$ l of **FAM-Cyclic-3', 5'-AMP** (200 nM) to each well designated "Positive Control," "Test Inhibitor," and "Substrate Control."
- 4) Add 45  $\mu$ l of **PDE assay buffer** to each well designated "Blank" and add 20  $\mu$ l of **PDE assay buffer** to each well designated "Substrate Control."
- 5) Add 5  $\mu$ l of inhibitor solution to each well designated "Test Inhibitor." For the wells labeled "Positive Control," "Substrate Control," and "Blank," add 5  $\mu$ l of the same solution without inhibitor (inhibitor buffer).
- 6) Thaw **Rat PDE4D** on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. Aliquot **Rat PDE4D** enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: Rat PDE4D 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 $\mu$ l	25 $\mu$ l	25 $\mu$ l	-
PDE assay buffer	-	-	20 $\mu$ l	45 $\mu$ l
Inhibitor (in PDE assay buffer)	-	5 $\mu$ l	-	-
Inhibitor Buffer (no inhibitor)	5 $\mu$ l	-	5 $\mu$ l	5 $\mu$ l
Rat PDE4D (1 pg/ $\mu$ l)	20 $\mu$ l	20 $\mu$ l	-	-
<b>Total</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>

- 7) Dilute **Rat PDE4D** in **PDE assay buffer** to 1 pg/ $\mu$ l (0.02 ng/reaction)\*. Initiate reaction by adding 20  $\mu$ l of diluted **Rat PDE4D** (1 pg/ $\mu$ l) 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.*

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)

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  $\mu$ l 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  $485 \pm 5$  nm and detection of emitted light ranging from  $528 \pm 10$  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".

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

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)

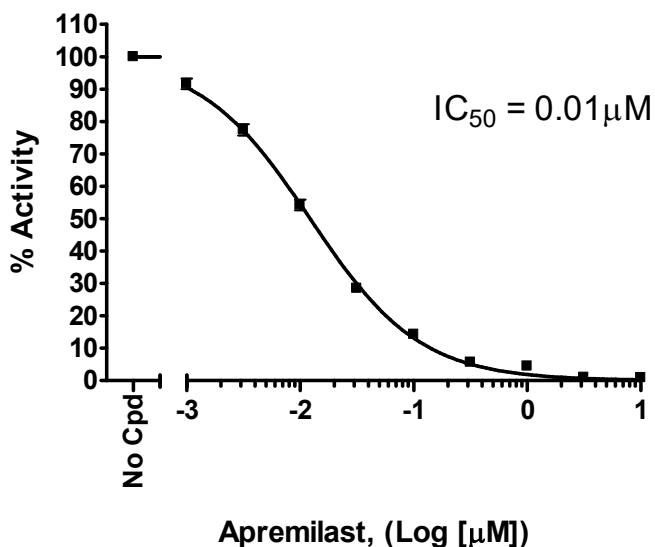


6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**EXAMPLE OF ASSAY RESULTS:**

**Rat PDE4D Activity**

Substrate Conc. = 100nM (cAMP)



Inhibition of Rat PDE4D by Apremilast measured using the *RAT PDE4D Assay Kit*, BPS Bioscience #79629. 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 Name</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
Rat PDE4B	60049	5 µg
Rat 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
PDE4C1	60044	5 µg
PDE4D2	60048	5 µg
PDE4D3	60046	5 µg
PDE4D7	60047	5 µg
PDE4A1A	60040	10 µg
PDE4B1	60041	10 µg
PDE4B2	60042	5 µg
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.
PDE4D2 Assay Kit	60345	96 rxns.
PDE4D 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](mailto:info@bpsbioscience.com)

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