

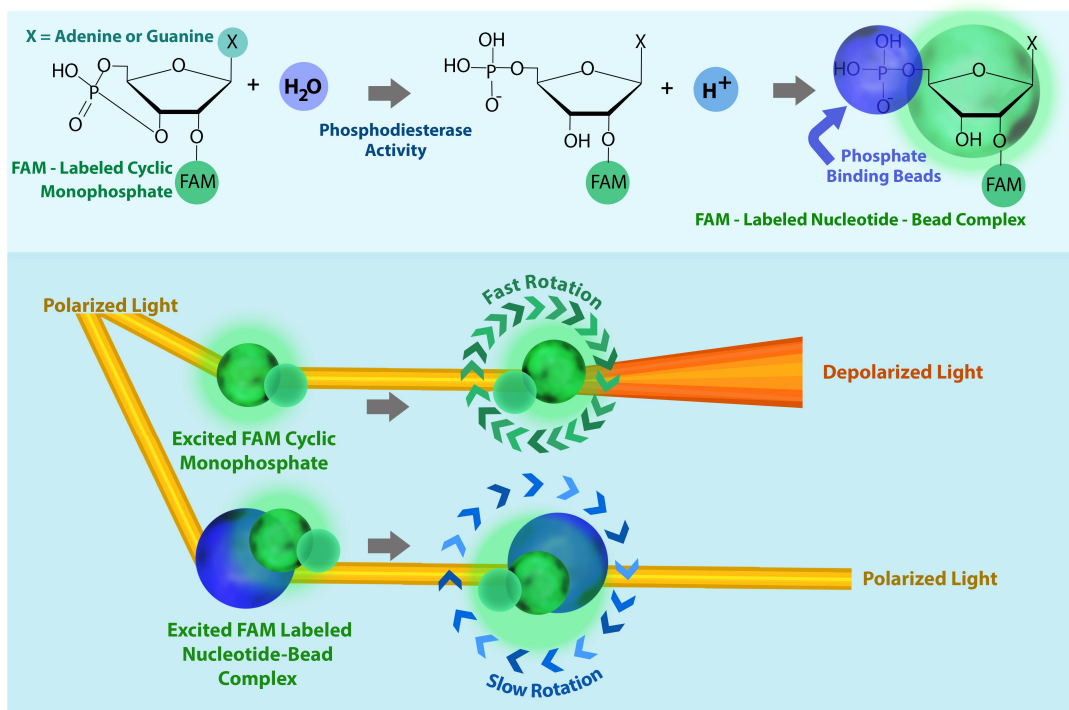
Data Sheet

PDE4D2 Assay Kit

Catalog # 60345
Size: 96 reactions

DESCRIPTION: Phosphodiesterases (PDEs) play an important role in dynamic regulation of cAMP and cGMP signaling. PDE4D is a regulator of airway smooth-muscle contractility, and has been identified as a potential risk predictor for ischemic stroke. Additionally, PDE4D has been associated with asthma pathophysiology and bone formation. The PDE4D gene encodes at least 9 different isoforms. The *PDE4D2 Assay Kit* is designed for identification of inhibitors of PDE4D2 using fluorescence polarization. The assay is based on the binding of a fluorescent nucleotide monophosphate generated by PDE4D2 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 *PDE4D2 Assay Kit* comes in a convenient 96-well format, with purified PDE4D2 enzyme, fluorescently labeled PDE4D substrate (cAMP), binding agent, and PDE assay buffer for 100 enzyme reactions. The key to the *PDE4D2 Assay Kit* is the specific binding agent. Using this kit, only two simple steps on a microtiter plate are required for PDE4D2 reactions. First, the fluorescently labeled cAMP is incubated with a sample containing PDE4D2 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	
60048	PDE4D2 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

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.

REFERENCE: Chandrasekaran A, *et al.*, *Cell Signal*. 2008; **20(1)**: 139-53.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Dilute 20 µM FAM-Cyclic-3', 5'-AMP stock 100-fold with PDE buffer to make a 200 nM solution. Make only sufficient quantity needed for the assay; store remaining 20 µM stock solution in aliquots at -20°C.
- 2) Add 25 µl of FAM-Cyclic-3',5'-AMP (200 nM) to each well designated "Positive Control", "Test Inhibitor", and "Substrate Control".

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

- 3) Add 20 μ l of PDE assay buffer to each well designated "Substrate Control" and 45 μ l of PDE assay buffer to each well designated "Blank".
- 4) 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).
- 5) Thaw PDE4D2 on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. Aliquot PDE4D2 enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -70°C immediately. *Note: PDE4D2 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 6) Dilute PDE4D2 in PDE buffer to 2.5 pg/ μ l (0.05 ng/reaction)*. Initiate reaction by adding 20 μ l of diluted PDE4D2 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.*
- 7) Incubate at room temperature for 1 hour.

	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
PDE4D2 (2.5 pg/ μ l)	20 μ l	20 μ l	–	–
Total	50 μl	50 μl	50 μl	50 μl

Step 2:

- 1) Mix **binding agent** thoroughly and dilute **binding agent** 1:100 with binding agent diluent.
- 2) Add 100 μ l diluted binding agent to each microwell. Incubate at room temperature for 1 hour 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.

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

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**

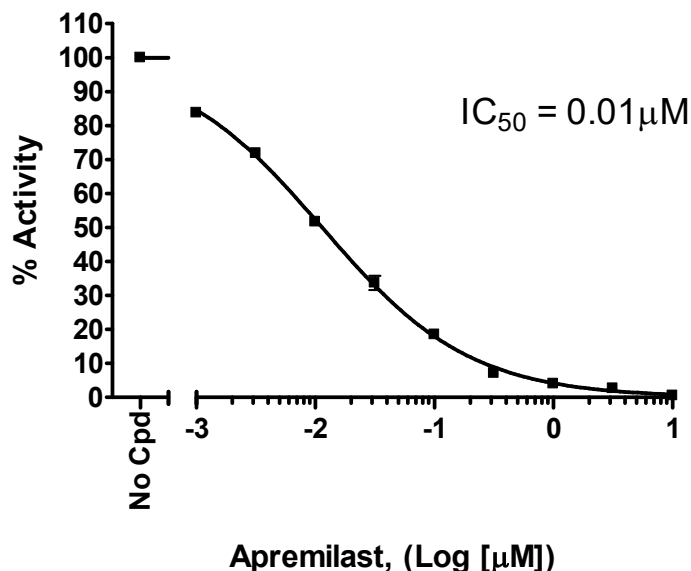
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

EXAMPLE OF ASSAY RESULTS:

PDE4D2 Activity

Substrate Conc. = 100nM (cAMP)



Inhibition of PDE4D2 by Apremilast, measured using the *PDE4D2 Assay Kit*, BPS Bioscience # 60345. 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*

RELATED PRODUCTS :

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
PDE4D2	60048	5 μg
PDE4D3	60046	5 μg
PDE4D7	60047	5 μg
PDE Assay Kit	60300	96 rxns.
PDE4D2 Assay Kit	60345	96 rxns.
PDE4D Cell-Based Activity Assay Kit	60505	500 rxns.
PDE4D2 TR-FRET Assay Kit	60707	96 rxns.
PDE4D3 TR-FRET Assay Kit	60701	96 rxns.
PDE4D7 TR-FRET Assay Kit	60708	96 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