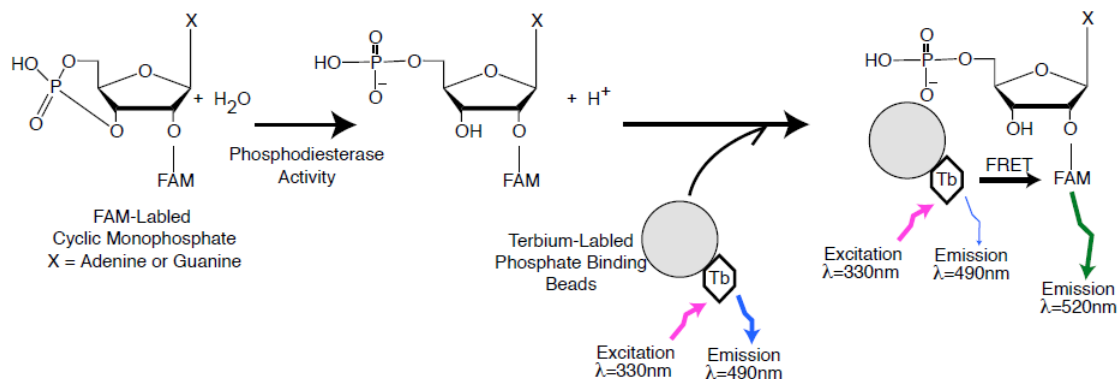


Data Sheet
PDE4D7 TR-FRET Assay Kit
Catalog # 60708

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 *PDE4D7 TR-FRET Assay Kit* is designed for identification of inhibitors of PDE4D7 using TR-FRET (Time Resolved Fluorescence Resonance Energy Transfer) technology. The assay is based on the generation of FAM-labeled nucleotide monophosphates by the phosphodiesterase. These phosphate groups bind to terbium-labeled nanoparticles, resulting in energy transfer from the terbium to the FAM, which emits a fluorescent signal at 520 nm. The change in fluorescent intensity can be easily measured using a fluorescence plate reader.



The *PDE4D7 TR-FRET Assay Kit* comes in a convenient 96-well format, with purified PDE4D7 enzyme, fluorescently labeled PDE substrate (cAMP), binding agent, and PDE assay buffer for 100 enzyme reactions. Using this kit, only two simple steps on a microtiter plate are required for the PDE4D7 activity assay. First, the fluorescent-labeled cAMP is incubated with a sample containing PDE4D7 for 1 hour. Second, a binding agent and a terbium donor are added to the reaction mix and incubated for 1 hour. Then, fluorescence intensity can be measured using a fluorescence reader.

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

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

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

COMPONENTS:

| Catalog # | Component | Amount | Storage | |
|-----------|-------------------------------------|--------|------------|-------------------------------------|
| 60047 | PDE4D7 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 | |
| 60394 | Tb donor | 50 µl | -80°C | |
| 60390 | Binding Agent | 200 µl | +4°C | |
| 78422 | Binding Buffer A | 20 ml | +4°C | |
| 78423 | Binding Buffer B | 20 ml | +4°C | |
| 79685 | Black, low binding microtiter plate | 1 | Room temp. | |

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable of measuring Time Resolved Fluorescence Resonance Energy Transfer (TR-FRET)

REFERENCE(S): Bender. A., *et al. Pharmacol. Rev.* 2006; **58**: 488-520.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Protocol for PDE4D7 assay

Step 1:

- 1) Dilute 20 µM FAM-Cyclic-3',5'-AMP substrate stock solution 100-fold with PDE buffer to make a 200 nM solution.
- 2) Add 25 µl of FAM-Cyclic-3',5'-AMP (200 nM) to each well designated "Substrate Control", "Positive Control", and "Test Inhibitor". Add 25 µl of PDE assay buffer to each well designated "Blank".
- 3) Add 5 µl of inhibitor solution to each well designated "Test Inhibitor". Add 5 µl of the same solution without inhibitor (inhibitor buffer) to the "Blank", "Substrate Control" and "Positive Control".
- 4) Thaw PDE4D7 on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full contents of the tube. *Note: PDE4D7 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 5) Dilute PDE4D7 in PDE buffer to 5 pg/µl (100 pg/reaction) in PDE buffer*. Add 20 µl of PDE assay buffer to the wells designated as the "Blank" and "Substrate Control". Initiate reaction by adding 20 µl of PDE4D7 (5 pg/µl) to the wells

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

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.*

| | Blank | Substrate Control | Positive Control | Test Inhibitor |
|---------------------------------|--------------|-------------------|------------------|----------------|
| FAM-Cyclic-3',5'-AMP (200 nM) | – | 25 µl | 25 µl | 25 µl |
| PDE assay buffer | 45 µl | 20 µl | – | – |
| Test Inhibitor | – | – | – | 5 µl |
| Inhibitor Buffer (no inhibitor) | 5 µl | 5 µl | 5 µl | – |
| PDE4D7 (5 pg/µl) | – | – | 20 µl | 20 µl |
| Total | 50 µl | 50 µl | 50 µl | 50 µl |

6) Incubate at room temperature for 1 hour.

Step 2:

- 1) Make binding dilution buffer by mixing equal volumes of Binding buffer A and Binding buffer B. For example, mix 1 ml Binding buffer A with 1 ml Binding buffer B.
- 2) Mix **binding agent** thoroughly and dilute **binding agent** 1:50 with binding dilution buffer made in Step 1.
- 3) Add Tb donor (1:200 dilution) to the mixture in Step 2.
- 4) Add 100 µl to each well. Incubate at room temperature for 1 hour with slow shaking.
- 5) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

Instrument Settings

| | |
|-----------------------|---------------|
| Reading Mode | Time Resolved |
| Excitation Wavelength | 330±20 |
| Emission Wavelength | 490±10 |
| Lag Time | 50 µs |
| Integration Time | 50 µs |
| Excitation Wavelength | 330±20 |
| Emission Wavelength | 520±10 |
| Lag Time | 50 µs |
| Integration Time | 50 µs |

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

CALCULATING RESULTS:

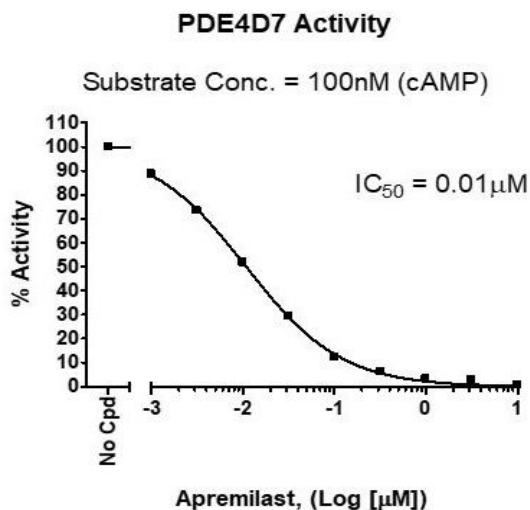
$$FRET = \frac{S_{520} - \left(\frac{Tb_{520}}{Tb_{490}} \times S_{490} \right)}{S_{490}} \times 1000$$

Where S_{520} = Sample 520 nm reading, S_{490} = Sample 490 nm reading, Tb_{520} = Tb only 520 nm reading, Tb_{490} = Tb only 490 nm reading. When percentage activity is calculated, the FRET value from substrate only control can be set as zero percent activity and the FRET value from positive control can be set as one hundred percent activity.

$$\% \text{ Activity} = \frac{FRET_s - FRET_{Sub}}{FRET_p - FRET_{Sub}} \times 100\%$$

Where $FRET_s$ = Sample FRET, $FRET_{Sub}$ = Substrate only control FRET, and $FRET_p$ = Positive control FRET.

EXAMPLE OF ASSAY RESULTS:



Inhibition of PDE4D7 by Apremilast, measured using the *PDE4D7 TR-FRET Assay Kit*, BPS Bioscience # 60708. 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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

RELATED PRODUCTS :

| <u>Product</u> | <u>Cat. #</u> | <u>Size</u> |
|--------------------------|---------------|-------------|
| PDE4D7 | 60047 | 5 µg |
| PDE4D3 | 60046 | 5 µg |
| PDE4D2 | 60048 | 5 µg |
| PDE4A1A | 60040 | 10 µg |
| PDE4A10 | 60038 | 10 µg |
| PDE4A4B | 60039 | 10 µg |
| PDE4B1 | 60041 | 10 µg |
| PDE4B2 | 60042 | 10 µg |
| PDE4C1 | 60044 | 5 µg |
| PDE4A FP Assay Kit | 60340 | 96 rxns. |
| PDE4B FP Assay Kit | 60343 | 96 rxns. |
| PDE4D3 FP Assay Kit | 60345 | 96 rxns. |
| PDE4D3 TR-FRET Assay Kit | 60701 | 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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com