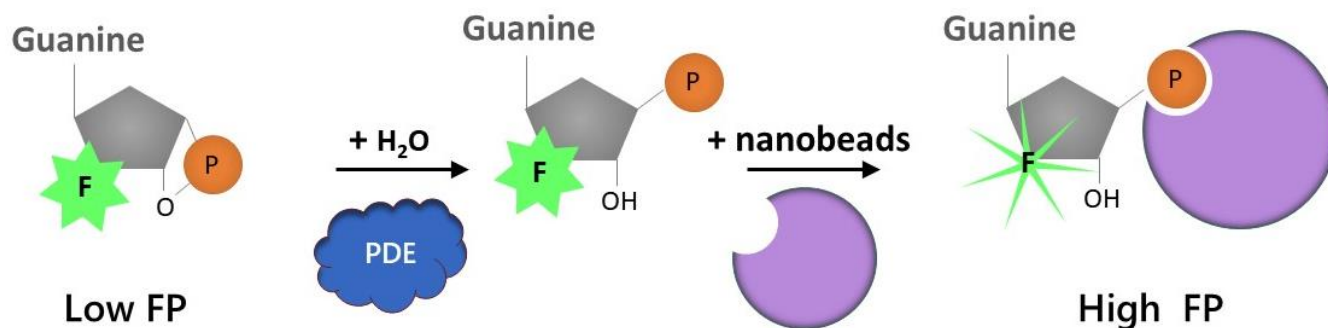


**Description**

The Phosphodiesterase 5A1 (PDE5A1) Assay Kit is a fluorescence polarization (FP), homogeneous 384-well assay kit designed for the screening & profiling of PDE5A1 inhibitors. This assay takes advantage of a specific fluorescent phosphate-binding nanoparticle. The kit contains enough purified recombinant PDE5A1, fluorescent probe, PDE assay buffer, Binding Agent, and diluent for 400 reactions.



*Figure 1: Illustration of the PDE5A1 assay principle.*

The assay uses a fluorescein-labeled cyclic guanine monophosphate (cGMP-FAM), in which the phosphate group is engaged within the cyclic nucleotide. This is a very small molecular that can rotate fast (low FP). PDE5A1 catalyzes the hydrolysis of the phosphodiester bond in the cyclic nucleotide and frees the phosphate group. In a second step the free phosphate group is recognized by a specific phosphate-binding nanobead (Binding Agent) leading to the formation of large complex, with restricted movement (high FP). FP is proportional to PDE activity.

This assay requires a fluorescent microplate reader **capable of measuring fluorescence polarization (FP)** to read the FP signal. For more information FP technology, visit our Tech Note: [FP, assay principles and applications](#).

**Background**

Phosphodiesterases (PDEs) play an important role in the dynamic regulation of the second messengers cAMP and cGMP signaling, by hydrolyzing them. PDE5A1, also known as cGMP-inhibited phosphodiesterase, is found in cerebellum, kidney, pancreas, lung, and vascular smooth muscle, and it has been implicated in cardiovascular function and fertility. Several inhibitors have been identified, such as sildenafil, vardenafil and tadalafil, for the treatment of pulmonary hypertension and erectile dysfunction. The reported side effects on vision may reveal new roles for PDE5 in the retina, so further studies into the functions of PDE5A1 and development of new inhibitors will be advantageous.

**Applications**

Study enzyme kinetics and screen small molecule inhibitors for drug discovery in high throughput screening (HTS) applications.

**Supplied Materials**

Catalog #	Name	Amount	Storage
60050	PDE5A1, GST-Tag*	>1 µg	-80°C
60201	20 µM FAM-Cyclic-3', 5'-GMP	50 µl	-80°C
60393	PDE Assay Buffer	25 ml	-20°C
60390	Binding Agent	250 µl	4°C
60392	cGMP Binding Agent Diluent	25 ml	4°C
79961	Low binding, black 384-well plate	1	Room Temp.

\* The concentration of protein is lot-specific and will be indicated on the tube containing the protein.

**Materials Required but Not Supplied**

Fluorescent plate reader capable of measuring fluorescence polarization.

**Stability**

This assay kit will perform optimally for up to **6 months** from date of receipt when the materials are stored as directed.

**Safety**

This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous and is harmful by inhalation, in contact with skin, eyes, clothing, and if swallowed. If contact occurs, wash thoroughly.

**Assay Protocol**

- All samples and controls should be tested in duplicate.
  - The assay should include “Blank”, “Negative Control”, “Positive Control” and “Test Inhibitor”.
  - It is recommended to use inhibitor Vardefanil at 0.1 µM as an internal control for the assay.
  - We recommend maintaining the diluted protein on ice during use.
  - For detailed information on protein handling please refer to [Protein FAQs \(bpsbioscience.com\)](https://www.bpsbioscience.com/protein-faqs).
1. Dilute 20 µM FAM-Cyclic-3', 5'-GMP stock solution 100-fold using PDE Assay Buffer to make a 200 nM dilution (you will need 12.5 µl/well).
  2. Add 12.5 µl of diluted FAM-Cyclic-3', 5'-GMP to the “Negative Control”, “Positive Control”, and “Test Inhibitor” wells.
  3. Add 22.5 µl of PDE Assay Buffer to the “Blank” wells.
  4. Add 10 µl of PDE Assay Buffer to the “Negative Control” wells.
  5. Prepare the Test Inhibitor (2.5 µl/well): for a titration prepare serial dilutions at concentrations 10-fold higher than the desired final concentrations. The final volume of the reaction is 25 µl.

5.1 If the Test Inhibitor is water-soluble, prepare serial dilutions 10-fold more concentrated than the desired final concentrations using PDE Assay Buffer.

For the positive and negative controls, use PDE Assay Buffer as Diluent Solution.

**OR**

5.2 If the Test inhibitor is soluble in DMSO, prepare the inhibitor in 100% DMSO at a concentration 100-fold higher than the highest desired concentration, then dilute the inhibitor 10-fold in PDE Assay Buffer to prepare the highest concentration of the 10-fold intermediate dilutions. The concentration of DMSO is now 10%.

Use 10% DMSO in PDE Assay Buffer (vol/vol) for the serial dilution to keep the concentration of DMSO constant.

For positive and negative controls, prepare 10% DMSO in PDE Assay Buffer (vol/vol) so that all wells contain the same amount of DMSO (Diluent Solution).

*Note: The final concentration of DMSO should not exceed 1%.*

6. Add 2.5 µl of the inhibitor serial dilution to the “Test Inhibitor” wells.
7. Add 2.5 µl of the Diluent Solution to the “Blank”, “Negative Control”, and “Positive Control” wells.
8. Thaw **PDE5A1** on ice. Briefly spin the tube containing the enzyme to recover its full content.

*Note: The concentration of protein is lot-specific and will be indicated on the tube containing the protein.*

9. Dilute PDE5A1 with PDE Assay Buffer to 10 pg/µl (10 µl/well).

*The diluted protein is not stable upon freeze/thaw. Do not freeze and re-use the diluted enzyme.*

10. Initiate the reaction by adding 10 µl of diluted PDE5A1 to the “Positive Control” and “Test Inhibitor” wells.
11. Incubate at Room Temperature (RT) for 1 hour.

Component	Blank	Negative Control	Positive Control	Test Inhibitor
FAM-Cyclic-3',5'-GMP (200 nM)	-	12.5 µl	12.5 µl	12.5 µl
PDE Assay Buffer	22.5 µl	10 µl	-	-
Test Inhibitor	-	-	-	2.5 µl
Diluent Solution	2.5 µl	2.5 µl	2.5 µl	-
Diluted PDE5A1 (10 pg/µl)	-	-	10 µl	10 µl
<b>Total</b>	<b>25 µl</b>	<b>25 µl</b>	<b>25 µl</b>	<b>25 µl</b>

12. Gently mix the tube containing the **Binding Agent** and dilute 1:100 with the **cGMP Binding Agent Diluent**.
13. Add 50 µl of diluted Binding Agent to each well.
14. Incubate at RT for 20 minutes with gentle agitation.
15. Read FP in a fluorescence plate reader capable of measuring fluorescence polarization ( $\lambda_{exc} = 485 \text{ nm}$ ;  $\lambda_{em} = 528 \text{ nm}$ ) and set to FP.
16. Subtract the “Blank” value from all other values.

**Calculating Results; Definition of Fluorescence Polarization:**

Users may ignore the G-factor when all experiments are performed using the same instrument since the G-factor is instrument-dependent.

If desired, the G-factor is set before measurements are performed. It needs to be determined by the investigator when not clearly indicated by the manufacturer. The instrument manual will contain information about how to establish the **G-factor**. For example, BPS Bioscience’s scientists use a Tecan M1000 fluorescent plate reader which has a G-factor set to 22 mP.

Instruments provide measurement in milli-Polarization = mP.

Results are calculated as follows.

1. Subtract the “Blank” value from all other values.
2. Calculate  $\Delta\text{mP}$  for all samples:

$$\Delta\text{mP} = (\text{mP value of the sample}) - (\text{mP of the Reference control})$$

Where mP refers to milli-Polarization values provided by the instrument and Reference control is the mP value obtained in the condition containing only the fluorescent probe (a condition in which the probe is in free state).

## Example of Assay Results

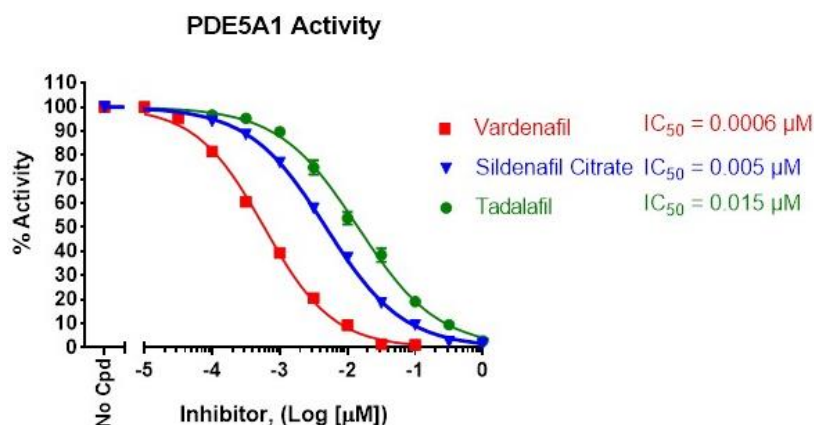


Figure 2: Titration of various inhibitors of PDE5A1.

PDE5A1 was incubated with increasing concentrations of sildenafil citrate (Axxora #LKT-S3313), vardenafil (Sigma #SML2103), and tadalafil (SelleckChem #S1512) in the presence of 100 nM FAM-Cyclic-3', 5'-GMP substrate. Fluorescence Polarization was measured using a Tecan M1000 fluorescent microplate reader. Results are expressed in % of activity, in which FP in the absence of inhibitor (positive control) is set to 100%.

Data shown is representative. For lot-specific information, contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## Troubleshooting Guide

Visit [bpsbioscience.com/assay-kits-faq](https://bpsbioscience.com/assay-kits-faq) for detailed troubleshooting instructions. For further questions, email [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## References

Maurice DH. *Front. Biosci.* 2005; 10: 1221-1228

## Related Products

Products	Catalog #	Size
PDE5A1 (Mouse), GST-Tag Recombinant	60051	5 μg
PDE5A1 (Mouse) Assay Kit	79602	96 reactions
PDE5A1 Assay Kit	60350	96 reactions

Version 110923