

## ab283404 – PLTP Inhibitor Drug Screening Kit (Fluorometric)

For the screening of potential PLTP inhibitors.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab283404>

### Storage and Stability

On receipt aliquot and store Positive Control at -20°C. Store rest of the kit components at 4°C, protected from light. Upon opening, use kit within 6 months.

### Materials Supplied

Item	Quantity	Storage Condition
PLTP Acceptor Molecule	1 mL	4°C
PLTP Assay Buffer (10X)	5 mL	4°C
PLTP Donor Molecule	1 mL	4°C
Positive Control (Rabbit Serum)	0.3 mL	-20°C

### Reagent Preparation

Before using the kit, spin the tubes prior to opening. Warm Assay Buffer to room temperature (RT) before use. All kit components are supplied as ready to be used. Keep on ice while in use.

### Assay Protocol

- **General Consideration for Using Fluorometer and Plate Reader:**
    - We recommend using a microtiter plate for the assay. The microtiter plates should be sealed as tightly as possible and incubate in a sealed, humidified chamber to prevent evaporation.
    - If using a regular fluorometer for sample reading, the samples should be diluted to 500 µL with 1X PLTP Assay Buffer before read.
  - **Preparation of Standard Curve:**
    - Standard curve is prepared by making serial dilutions of the donor molecule in isopropanol and subsequently recording the fluorescence intensity of each dilution, using isopropanol alone as a blank.
1. Prepare 6 test tubes labeled T0 to T5, each contains 0.2 mL of isopropanol; the tube labeled T5 should contain an additional 0.2 mL of isopropanol.
  2. Add 2 µL Donor Molecule to T5, vortex to mix well.
  3. Transfer 0.2 mL from T5 to T4. Mix and then transfer 0.2 mL from T4 to T3. Mix and then transfer 0.2 mL from T3 to T2. Mix and then transfer 0.2 mL from T2 to T1. The Donor Molecule solution contains 0.1 mM labeled lipids and thus the standard curve samples contain 0, 6.25, 12.5, 25, 50, 100 pmol donor molecules.
  4. Read the fluorescence intensity (Ex/Em: 465/535) of the standard samples from T0 to T5.
  5. Apply the fluorescence intensity values of the standard curve directly to your results to express specific activity of the plasma sample (moles/µL plasma/hr).

### Assay Procedure:

1. Prepare each testing sample in 160 µL of dH<sub>2</sub>O. Add 3 µL of Rabbit Serum.

- Prepare a blank that contains no rabbit serum as background. Prepare a positive control assay containing 3 µL of Rabbit Serum, but no testing inhibitors.
2. Prepare a Master Mix for each assay containing the follows. Mix well.

Donor Molecule	10 µL
Acceptor Molecule	10 µL
PLTP Assay Buffer (10X)	20 µL

3. Add 40 µL of the Master Mix into each sample including the blank and positive control also as prepared in Step 1. Mix well and incubate at 37°C for 30-60 minutes.
4. Measure the fluorescence intensity of the blank, test samples, and positive control using a fluorescence plate reader or fluorometer (Ex/Em: 465/535).

**Δ Note:** Due to the nature of the self-quenched probe, background fluorescence is usually high and therefore fluorescence intensity from each sample should be corrected by subtracting the background fluorescence intensity (without rabbit serum). The increase in fluorescence intensity with PLTP (Rabbit Serum) is usually 1.5 - 2.0-fold over blank (Organic solvent may increase background readings and therefore proper control may be needed if the testing inhibitor is prepared in organic solvent).

5. Comparison of the fluorescence intensity of the testing inhibitor with positive control to determine the inhibition efficiency of the testing inhibitors.

### Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.  
For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)  
[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)  
[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)