



sPLA₂ Assay Kit

Item No. 765001

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

GENERAL INFORMATION	3	Materials Supplied
	4	Precautions
	4	If You Have Problems
	4	Storage and Stability
	4	Materials Needed but Not Supplied
INTRODUCTION	5	Background
	5	About This Assay
PRE-ASSAY PREPARATION	7	Reagent Preparation
	8	Sample Preparation
ASSAY PROTOCOL	9	Plate Set Up
	11	Performing the Assay
ANALYSIS	12	Calculations
	13	Performance Characteristics
RESOURCES	15	Interferences
	16	Troubleshooting
	17	References
	18	Plate Template
	19	Notes
	19	Warranty and Limitation of Remedy

GENERAL INFORMATION

Materials Supplied

Item Number	Item	Quantity
765010	sPLA ₂ Assay Buffer (10X)	1 vial
765012	sPLA ₂ DTNB	4 vials
765015	sPLA ₂ Diheptanoyl Thio-PC (Substrate)	2 vials
765016	Bee venom PLA ₂ Control	1 vial
400014	96-Well Plate (Colorimetric Assay)	1 plate
400012	96-Well Cover Sheet	1 cover

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



WARNING: THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

Safety Data

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent *via* email to your institution.

Precautions

Please read these instructions carefully before beginning this assay.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888
Fax: 734-971-3641
Email: techserv@caymanchem.com
Hours: M-F 8:00 AM to 5:30 PM EST

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored at -20°C and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. A plate reader capable of measuring absorbances at 405 or 414 nm
2. Adjustable pipettes and a repeating pipettor
3. A source of pure water; glass distilled water or HPLC-grade water is acceptable

INTRODUCTION

Background

Phospholipase A₂ (PLA₂) catalyzes the hydrolysis of phospholipids at the *sn*-2 position yielding a free fatty acid and a lysophospholipid.¹ The release of arachidonic acid from membrane phospholipids by PLA₂ is believed to be a key step in the control of eicosanoid production within the cell.²

About This Assay

Cayman's secretory PLA₂ (sPLA₂) Assay Kit provides an accurate and convenient method for measurement of sPLA₂ activity. This assay uses the 1,2-dithio analog of diheptanoyl phosphatidylcholine which serves as a substrate for most PLA₂s (e.g., bee and cobra venoms, pancreatic, etc.) with the exception of cytosolic PLA₂.^{3,4} Upon hydrolysis of the thio ester bond at the *sn*-2 position by PLA₂, free thiols are detected using DTNB (5,5'-dithio-bis-(2-nitrobenzoic acid)) (see Figure 1, on page 6).

Reagent Preparation

Some of the kit components are in lyophilized form and need to be reconstituted prior to use. Follow the directions carefully to ensure proper volumes of water or Assay Buffer are used to reconstitute the vial components.

1. sPLA₂ Assay Buffer - (Item No. 765010)

Dilute 3 ml of Assay Buffer concentrate with 27 ml of HPLC-grade water. This final Assay Buffer (25 mM Tris-HCl, pH 7.5, containing 10 mM CaCl₂, 100 mM KCl, and 0.3 mM Triton X-100) should be used for reconstitution of substrate and dilution of samples prior to assaying.

2. sPLA₂ DTNB - (Item No. 765012)

Reconstitute the contents of one vial with 1.0 ml of HPLC-grade water to yield 10 mM DTNB in 0.4 M Tris-HCl, pH 8.0. Store the reconstituted reagent on ice in the dark and use within eight hours.

3. sPLA₂ Diheptanoyl Thio-PC (Substrate) - (Item No. 765015)

Evaporate the ethanolic solution of diheptanoyl thio-PC to dryness under a gentle stream of inert gas (e.g., nitrogen, argon). Reconstitute the contents of each vial with 12 ml of diluted Assay Buffer to achieve a final concentration of 1.66 mM. Make sure to vortex until the substrate solution becomes clear (high background absorbance may result if the substrate is not completely dissolved). The substrate, when stored at -20°C in diluted Assay Buffer, is stable for at least two weeks. *NOTE: If not using the entire plate, then reconstitute only one of the substrate vials.*

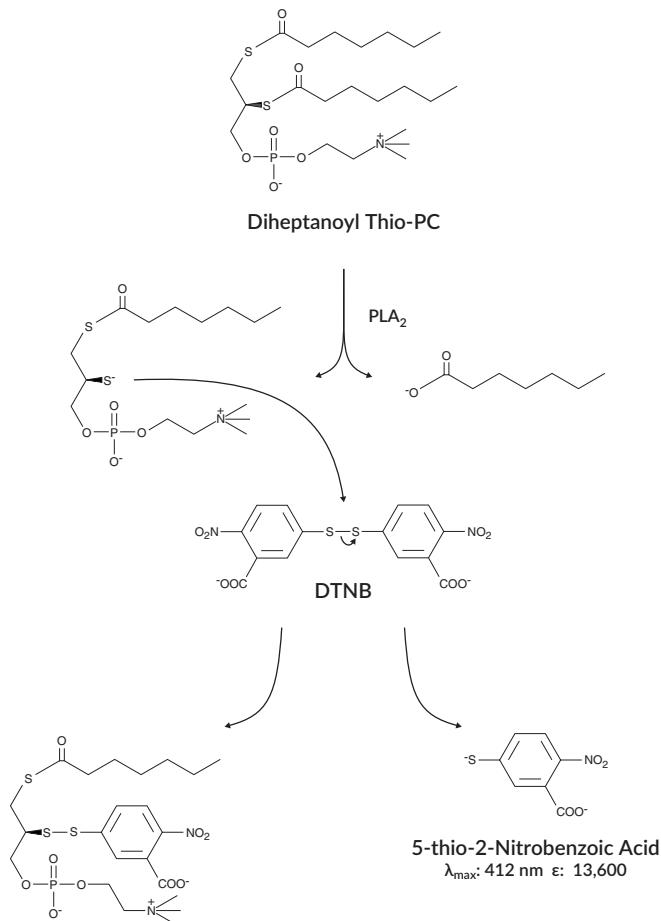


Figure 1. Scheme

4. Bee Venom PLA₂ Control - (Item No. 765016)

A solution of bee venom PLA₂ is supplied as a positive control. To avoid repeated freezing and thawing, the PLA₂ can be aliquoted into several small vials. Bee venom PLA₂, when stored at -20°C, is stable for one year. Transfer 10 µl of the supplied enzyme to another vial and dilute with 990 µl of diluted Assay Buffer prior to use. Store the enzyme on ice and use within one hour. A 10 µl aliquot of this diluted enzyme per well causes an increase of approximately 0.1 absorbance unit/minute under the standard assay conditions described below. The PLA₂ can be further diluted with Assay Buffer if a slower reaction rate is desired.

Sample Preparation

In general, any sPLA₂ sample can be measured by this assay. The sample must be free of particulates to avoid interferences in the absorbance measurement. Thiols, thiol 'scavengers', and PLA₂ inhibitors must be removed from the samples before performing the assay (extensive dialysis will eliminate most of the interfering substances of small molecular size). If the samples are too dilute, they can be concentrated using an Amicon centrifuge concentrator with a molecular weight cut-off of 3,000 Da.

Plate Set Up

There is no specific pattern for using the wells on the plate. However, it is necessary to have some wells (at least two) designated as non-enzymatic controls. The absorbance rate of these wells must then be subtracted from the absorbance rate measured in the sample wells. We suggest that you have at least two wells designated as positive controls, and that you record the contents of each well on the template sheet provided (see page 18).

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blk	Blk	S7	S7	S15	S15	S23	S23	S31	S31	S39	S39
B	+	+	S8	S8	S16	S16	S24	S24	S32	S32	S40	S40
C	S1	S1	S9	S9	S17	S17	S25	S25	S33	S33	S41	S41
D	S2	S2	S10	S10	S18	S18	S26	S26	S34	S34	S42	S42
E	S3	S3	S11	S11	S19	S19	S27	S27	S35	S35	S43	S43
F	S4	S4	S12	S12	S20	S20	S28	S28	S36	S36	S44	S44
G	S5	S5	S13	S13	S21	S21	S29	S29	S37	S37	S45	S45
H	S6	S6	S14	S14	S22	S22	S30	S30	S38	S38	S46	S46

Blk = Blank Wells

+ = Positive Control Wells

S1-S46 = Sample Wells

Figure 2. Sample plate format

Pipetting Hints

- It is recommended that an adjustable pipette be used to deliver Substrate, DTNB, and Buffer to the wells. This saves time and helps to maintain more precise times of incubation.
- Use different tips to pipette Substrate, DTNB, and sample.
- Before pipetting each reagent, equilibrate the pipette tip in that reagent (*i.e.*, slowly fill the tip and gently expel the contents, repeat several times).
- Do not expose the pipette tip to the reagent(s) already in the well.

General Information

- The final volume of the assay is 225 μl in all wells.
- It is not necessary to use all the wells on the plate at one time.
- If the appropriate enzyme dilution is not known, it may be necessary to assay at several dilutions.
- Use the Assay Buffer (dilute) in the assay.

Performing the Assay

1. **Blank Wells (Non-enzymatic controls)** - add 10 μl DTNB and 15 μl Assay Buffer to at least two wells (if performing inhibitor studies,* add 5 μl DMSO and 10 μl Assay Buffer instead of 15 μl Assay Buffer).
2. **Positive Control Wells (Bee Venom PLA₂)** - add 10 μl DTNB, 10 μl Bee Venom PLA₂, and 5 μl Assay Buffer to at least two wells (if performing inhibitor studies,* add 5 μl DMSO instead of 5 μl Assay Buffer).
3. **Sample Wells** - add 10 μl DTNB, 10 μl sample, and 5 μl Assay Buffer to at least two wells (if performing inhibitor studies,* add 5 μl of inhibitor dissolved in DMSO instead of 5 μl Assay Buffer). To obtain reproducible results, the amount of PLA₂ added to the well should cause an absorbance increase between 0.01 and 0.1/min. When necessary, samples should be concentrated or diluted with Assay Buffer to bring the enzymatic activity to this level. *NOTE: The amount of sample added to the well should always be 10 μl .*
4. Initiate the reactions by adding 200 μl Substrate Solution to all the wells. Make sure to note the precise time you started and add the Substrate Solution as quickly as possible.
5. Carefully shake the plate to mix.
6. Read the absorbance every minute at 414 (or 405) nm using a plate reader to obtain at least five time points.

*Inhibitors should be dissolved in DMSO and should be added to the assay in a final volume of 5 μl . In the event that the appropriate concentration of inhibitor is completely unknown, we recommend that several different dilutions of the inhibitor in DMSO be made.

Calculations

- Determine change in absorbance (ΔA_{414}) per minute by:
 - Plotting the absorbance values as a function of time to obtain the slope (rate) of the linear portion of the curve (an example is shown on page 14 using bee venom, snake venom, human synovial, and bovine pancreas PLA₂s).

OR

- Select two point on the linear portion of the curve and determine the change in absorbance during that time using the following equation:

$$\Delta A_{414} = \frac{A_{414}(\text{Time 2}) - A_{414}(\text{Time 1})}{\text{Time 2 (min.)} - \text{Time 1 (min.)}}$$

- Determine the rate of $\Delta A_{414}/\text{min}$ for the non-enzymatic controls (Blanks) and subtract this rate from that of the sample wells.
- Use the following formula to calculate the sPLA₂ activity. The reaction rate at 414 nm can be determined using the DTNB extinction coefficient of 10.66 mM⁻¹. **One unit of enzyme hydrolyzes one μmol of diheptanoyl Thio-PC per minute at 25°C.

$$\text{sPLA}_2 \text{ Activity } (\mu\text{mol}/\text{min}/\text{ml}) = \frac{\Delta A_{414}/\text{min.}}{10.66 \text{ mM}^{-1}} \times \frac{0.225 \text{ ml}}{0.01 \text{ ml}} \times \text{Sample dilution}$$

**The actual extinction coefficient for DTNB at 414 nm is 13.6 mM⁻¹cm⁻¹. This value has been adjusted for the pathlength of the solution in the well (0.784 cm). The extinction coefficient for DTNB at 405 nm is 12.8 mM⁻¹cm⁻¹. The adjusted value is 10.0 mM⁻¹.

Performance Characteristics

Sensitivity:

The detection range of this assay is from 0.02 to 0.2 $\mu\text{mol}/\text{min}/\text{ml}$ of sPLA₂ activity which is equivalent to an absorbance increase of 0.01 to 0.1 per minute.

Precision:

When a series of 16 bee venom PLA₂ measurements were performed on the same day, the intra-assay coefficient of variation was 2.5%. When a series of 16 bee venom PLA₂ measurements were performed on five different days under the same experimental conditions, the inter-assay coefficient of variation was 4.2%.

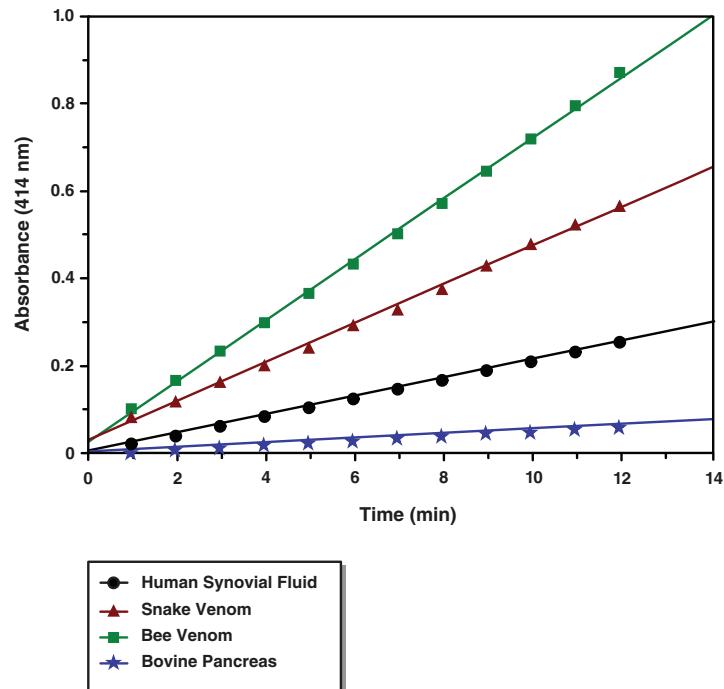


Figure 3. Absorbance times for bee venom, snake venom, human synovial, and bovine pancreas PLA₂s

RESOURCES

Interferences

1. Solvents

A dramatic decrease in enzymatic activity was observed when ethanol is added to the assay. Methanol also affects enzymatic activity but to a lesser extent. The addition of DMSO has little effect on enzymatic activity.

2. Culture Media and Buffers

All buffers and media should be tested for background absorbance before doing any experiments. If the initial background absorbance is higher than 0.3 absorbance units then the samples should be diluted in Assay Buffer or water before performing the assay. Tris, HEPES, and phosphate buffers (25 mM) work in the assay but imidazole buffers have high background absorbances. The following media also work in the assay: Grace's, RPMI 1640, DMEM, HAM's F12, and IMDM.

3. Thiols and Thiol-Scavengers

Samples containing thiols (e.g., glutathione, cysteine, dithiothreitol, or 2-mercaptoethanol) will exhibit high background absorbances and interfere with PLA₂ activity determination. Samples containing thiol-scavengers (e.g., N-ethylmaleimide) will inhibit color development. Extensive dialysis will eliminate most of the interfering substances of small molecular size.

Troubleshooting

Problem	Possible Causes	Recommended Solutions
Erratic values; dispersion of duplicates/triplicates	A. Poor pipetting/ technique B. Bubble in the well(s)	A. Be careful not to splash the contents of the wells B. Carefully tap the side of the plate with your finger to remove bubbles
No color development	A. DTNB or sample was not added to well(s) B. The enzymatic activity was too low	A. Make sure to add all components to the wells B. Standardize the assay with Bee Venom PLA ₂
The color development was too fast	Too much enzyme added to well(s)	Dilute your samples with diluted Assay Buffer and re-assay
High background absorbance	A. Substrate not in solution B. Thiols present in sample	A. Make sure to vortex the substrate until a clear solution is made B. Remove thiols or thiol reagents from sample
The reaction rate is not linear at high absorbance	Plate reader not sensitive enough at high absorbance	A. Use only the points at lower concentrations in the linear portion for making the curve B. Dilute your sample with diluted Assay Buffer and re-assay

References

1. Dennis, E.A. Phospholipases. *Enzymes XVI*, 307-353 (1983).
2. Dennis, E.A. Diversity of group types, regulation, and function of phospholipase A₂. *J. Biol. Chem.* **269**, 13057-13060 (1994).
3. Hendrickson, H.S., Kendrickson, E.K., and Dybvig, R.H. Chiral synthesis of a dithiolester analog of phosphatidylcholine as a substrate for the assay of phospholipase A₂. *J. Lipid Res.* **24**, 1532-1537 (1983).
4. Reynolds, L.J., Hughes, L.L., and Dennis, E.A. Analysis of human synovial fluid phospholipase A₂ on short chain phosphatidylcholine-mixed micelles: Development of a spectrophotometric assay suitable for a microtiterplate reader. *Anal. Biochem.* **204**, 190-197 (1992).

