

Phospholipase A2 Assay Kit ab133089

[3 References](#) [1 Image](#)

Overview

Product name	Phospholipase A2 Assay Kit
Detection method	Colorimetric
Sample type	Plasma
Assay type	Enzyme activity (quantitative)
Species reactivity	Reacts with: Mammals, Other species
Product overview	Phospholipase A2 Assay Kit (ab133089) provides an accurate and convenient method for measurement of PLA2 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. 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)).

If you want to detect cytosolic PLA2, we recommend using **Cytosolic Phospholipase A2 Assay Kit (ab133090)**.

Notes

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate reader

Properties

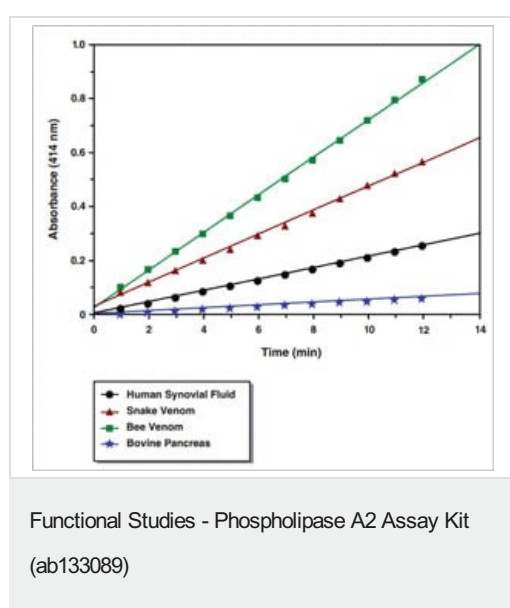
Storage instructions Store at -20°C. Please refer to protocols.

Components	96 tests
96-Well Plate (Colorimetric Assay)	1 x 96 tests
96-Well Plate Cover	1 unit
Bee venom PLA2 Control	1 vial
PLA2 Assay Buffer (10X)	1 vial

Components	96 tests
PLA2 Diheptanoyl Thio-PC (Substrate)	2 vials
PLA2 DTNB	4 vials

Function	Thought to participate in the regulation of the phospholipid metabolism in biomembranes including eicosanoid biosynthesis. Catalyzes the calcium-dependent hydrolysis of the 2-acyl groups in 3-sn-phosphoglycerides.
Sequence similarities	Belongs to the phospholipase A2 family.
Cellular localization	Membrane.

Images



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

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors