

Phospholipase A2 Assay Kit ab133089

3 References 画像数 1

製品の概要

製品名	Phospholipase A2 Assay Kit
検出方法	Colorimetric
サンプルの種類	Plasma
アッセイタイプ	Enzyme activity (quantitative)
種交差性	交差種: Mammals, Other species
製品の概要	<p>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)).</p> <p>If you want to detect cytosolic PLA2, we recommend using <u>Cytosolic Phospholipase A2 Assay Kit (ab133090)</u>.</p>
特記事項	<p>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.</p> <p>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.</p>
試験プラットフォーム	Microplate reader

製品の特性

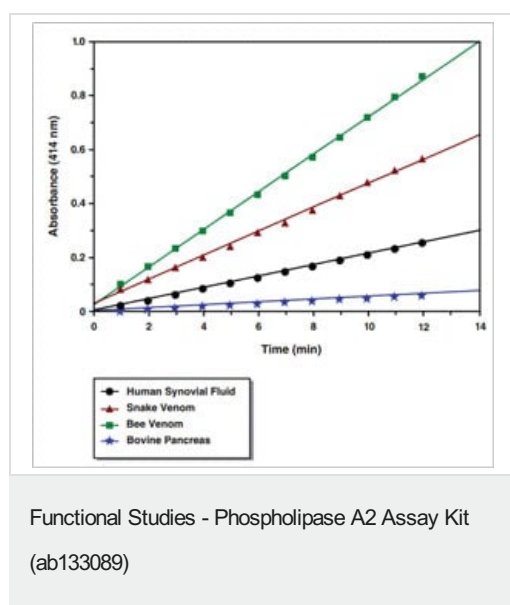
保存方法 Store at -20°C. Please refer to protocols.

内容	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

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

機能	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.
配列類似性	Belongs to the phospholipase A2 family.
細胞内局在	Membrane.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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