

PRODUCT INFORMATION

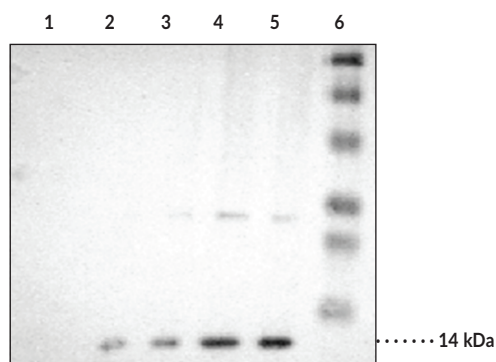


sPLA₂ (human Type IIA) Polyclonal Antiserum Item No. 160502

Overview and Properties

Contents:	This vial contains 100 µl of lyophilized antiserum
Synonyms:	GIIC sPLA ₂ , Group IIA Phospholipase A ₂ , NPS-PLA ₂ , Phosphatidylcholine 2-Acylhydrolase 2A, Phospholipase A ₂ Membrane Associated
Immunogen:	Human recombinant synovial sPLA ₂ (Type IIA)
Cross Reactivity:	(-) sPLA ₂ (Type V)
Species Reactivity:	(+) Human; (-) Bee venom
Uniprot No.:	P14555
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	Lyophilized from serum. Reconstitute in 100 µl of double distilled water.
Host:	Rabbit
Applications:	Western Blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Bee Venom PLA₂ (-) (2 µg)
Lane 2: sPLA₂ (IIA) (10 ng)
Lane 3: sPLA₂ (IIA) (20 ng)
Lane 4: sPLA₂ (IIA) (50 ng)
Lane 5: sPLA₂ (IIA) (100 ng)
Lane 6: Precision Plus Protein Standard

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.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/16/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

sPLA₂ 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 Human sPLA₂ (Type II_A) is secreted, 14 kDa protein which is dependent on Ca²⁺ for optimal activity. Human sPLA₂ (Type II_A) is 79% homologous to rat platelet and murine Type II sPLA₂s. Cayman's sPLA₂ antibody detects a band of approximately 15 kD by Western blot. For other applications, the working concentration of the antiserum must be determined empirically.

References

1. Dennis, E.A. Phospholipases in *The Enzymes*. Boyer, P., editor, XVI, Burlington:Elsevier (1983).
2. Dennis, E.A. Diversity of group types, regulation, and function of phospholipase A₂. *J. Biol. Chem.* **269**(18), 13057-13060 (1994).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM