

# Product Information



## sPLA<sub>2</sub> (human Type IIA) Western Ready Control

Catalog No. 10011382

<b>Synonyms:</b>	Secretory Phospholipase A <sub>2</sub> (human Type IIA), sPLA <sub>2</sub> (human synovial)
<b>Source:</b>	Human recombinant protein expressed in <i>E. coli</i>
<b>M<sub>r</sub>:</b>	16 kDa
<b>Specificity:</b>	Single band on a western blot when probed with Cayman's sPLA <sub>2</sub> (human Type IIA) Monoclonal Antibody (Catalog No. 160500).
<b>Supplied in:</b>	Laemmli buffer (2% SDS, 10% glycerol, 62.5 mM Tris-HCl, pH 6.8, 5% 2-mercaptoethanol, bromophenol blue)
<b>Stability:</b>	≥2 years at -20°C
<b>Applications:</b>	Positive control in western blotting

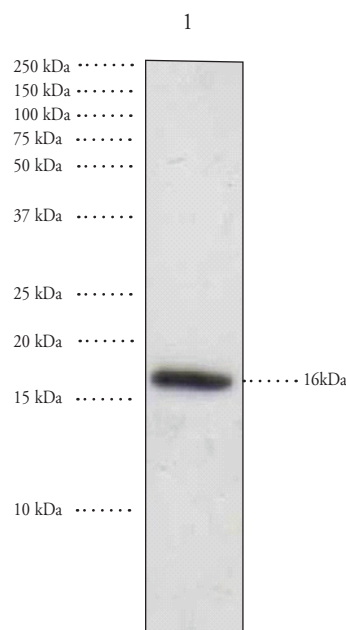
Phospholipase A<sub>2</sub> (PLA<sub>2</sub>) catalyzes the hydrolysis of phospholipids at the *sn*-2 position yielding a free fatty acid and a lysophospholipid.<sup>1</sup> The release of arachidonic acid from the membrane phospholipids by PLA<sub>2</sub> is believed to be a key step in the control of eicosanoid production within the cell.<sup>2</sup> Human synovial sPLA<sub>2</sub> is a secreted 14 kDa protein which is dependent on Ca<sup>2+</sup> for optimal activity. Human synovial sPLA<sub>2</sub> is 79% homologous to rat platelet and murine type II sPLA<sub>2</sub>s.

### Laboratory Procedures

Western Ready Controls are designed to serve as positive controls on western blots. This sPLA<sub>2</sub> (human Type IIA) Western Ready Control is supplied pre-boiled in 55 µl of Laemmli buffer. To use, thaw and load 10 µl of sPLA<sub>2</sub> (human Type IIA) Western Ready Control per well and follow a standard western blotting protocol. The amount needed per well will depend on the antibody and detection system used.

### References

1. Dennis, E.A. Phospholipases. *Enzymes XVI*, 307-353 (1983).
2. Dennis, E.A. Diversity of group types, regulation, and function of phospholipase A<sub>2</sub>. *J. Biol. Chem.* **269**, 13057-13060 (1994).



### Related Products

sPLA<sub>2</sub> (human Type IIA) Monoclonal Antibody - Cat. No. 160500 • sPLA<sub>2</sub> (human Type V) Monoclonal Antibody - Cat. No. 160510 • sPLA<sub>2</sub> (human Type IIA) EIA Kit - Cat. No. 585000 • sPLA<sub>2</sub> Assay Kit - Cat. No. 765001 • sPLA<sub>2</sub> (human recombinant Type V) - Cat. No. 10009563

**WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/26/2008

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com