

# PRODUCT INFORMATION

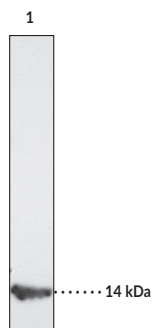


## sPLA<sub>2</sub> (human Type IIA) Monoclonal Antibody (Clone SCACC353) Item No. 160500

### Overview and Properties

<b>Contents:</b>	This vial contains 350 µg of protein G-purified monoclonal antibody.
<b>Synonyms:</b>	Secretory Phospholipase A <sub>2</sub> (human synovial)
<b>Immunogen:</b>	Human recombinant synovial sPLA <sub>2</sub> Type IIA
<b>Cross Reactivity:</b>	(-) Human sPLA <sub>2</sub> (type V), Bee venom sPLA <sub>2</sub> (Type III)
<b>Species Reactivity:</b>	(+) Human sPLA <sub>2</sub> (type IIA)
<b>Uniprot No.:</b>	P14555
<b>Form:</b>	Solid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol, 0.1% BSA and 0.02% NaN <sub>3</sub>
<b>Clone:</b>	SCACC353
<b>Host:</b>	Mouse
<b>Isotype:</b>	IgG1
<b>Applications:</b>	Immunoprecipitation (IP) and Western blot (WB); the recommended starting dilution is 1:150 and 1:200, respectively. <sup>1</sup> Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. <sup>2,3</sup>

### Image



Lane 1: Human sPLA<sub>2</sub> type II (250 ng)

### Description

PLA<sub>2</sub> catalyzes the hydrolysis of phospholipids at the *sn*-2 position yielding a free fatty acid and a lysophospholipid.<sup>4</sup> The release of arachidonic acid from membrane phospholipids by PLA<sub>2</sub> is believed to be a key step in the control of eicosanoid production within the cell.<sup>5</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 mouse type II sPLA<sub>2</sub>s.

### References

1. Lousse, J.-C., Defrère, S., Colette, S., et al. *Hum. Reprod.* **25**(3), 734-741 (2010).
2. Hulthe, J., Wiklund, O., Hurt-Camejo, E., et al. *Arterioscler. Thromb. Vasc. Biol.* **21**, 269-274 (2001).
3. Peilot, H., Rosengren, B., Bondjers, G., et al. *J. Biol. Chem.* **275** (30), 22895-22904 (2000).
4. Dennis, E.A. *Enzymes XVI*, 307-353 (1983).
5. Dennis, E.A. *J. Biol. Chem.* **269**, 13057-13060 (1994).

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, 11/15/2023

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