

# sPLA<sub>2</sub> activity kit

The sPLA<sub>2</sub> activity kit is a higher throughput assay for the quantification of up to 42 samples in duplicate per 96 well kit. This kit provides safe, sensitive, and reliable results in under an hour.

Secretory Phospholipase A<sub>2</sub>s (sPLA<sub>2</sub>s) are a group of related enzymes that upon activation, cleave an acyl ester bond in the sn-2 position in glycerophospholipid. These are extracellular enzymes that have high disulfide bond content, low molecular mass (14KDa), and require mM levels of Ca<sup>+</sup> for catalysis. sPLA<sub>2</sub> enzymes play a crucial role in the generation of arachidonic acid, which is further metabolized to proinflammatory eicosanoids.

Citations: 2

[View Online »](#)

## Ordering Information

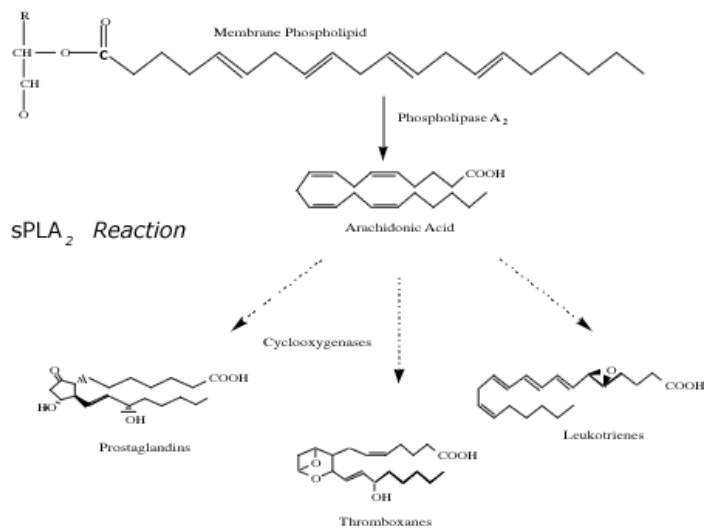
[Order Online »](#)

ADI-907-002	96 wells
-------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Higher throughput - assay up to 42 samples in duplicate per 96 well kit
- Sensitive - detect < 3 U/ml
- Fast - results in just 40 minutes!



sPLA<sub>2</sub> Reaction

## Handling & Storage

**Use/Stability** Store all components 4°C, except Substrate at -20°C.

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Secretory phospholipase A2

**Application** Activity assay, Colorimetric detection

**Application Notes** For the quantitative determination of sPLA<sub>2</sub> in culture supernatants, cell lysates, and buffers from any species.

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter plate, Standard, Reaction buffer concentrate, Substrate, Stop reagent, Color reagent

**Sensitivity** 2.34 Units/ml (range 5 – 80 Units/mL)

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)