

sPLA₂ activity kit

The sPLA₂ 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 A2s (sPLA₂s) 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⁺ for catalysis. sPLA₂ enzymes play a crucial role in the generation of arachidonic acid, which is further metabolized to proinflammatory eicosanoids.

Citations: 2

[View Online »](#)

Ordering Information

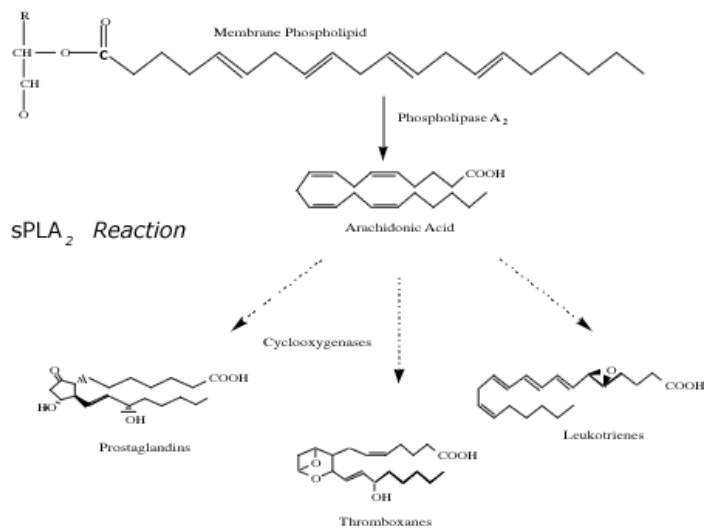
[Order Online »](#)

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!



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 sPLA₂ 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)



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)