

# Complement C3a des Arg (human), ELISA kit

## The Complement C3a des Arg (human), ELISA kit is a colorimetric competitive enzyme immunoassay kit

The Complement C3a des Arg (human), EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit comes ready-to-use with sample preparation reagents included. The Complement C3a des Arg (human), EIA kit is specific, offering low cross-reactivity to structurally related molecules, and sensitive with quantification of as little as 120 pg/mL Complement C3a des Arg.

Activation of the complement cascade is initiated by multiple pathways, including C1q binding of antigen-antibody complexes on pathogens (classical pathway), binding of serum lectin to pathogen-unique mannose residue patterns (MBL pathway), or the spontaneous hydrolysis of C3 and subsequent processing to C3 convertase (alternative pathway). C3a and C4a are two complement cleavage products of 77- and 76-amino acid residues, respectively, that are liberated into the bloodstream and processed to their more stable metabolites C3a des Arg and C4a des Arg. Although both fragments have anaphylactic properties, C3a is considered a more potent stimulator of smooth muscle cell contraction, histamine release, and inflammation than C4a. Complement activation is associated with sepsis, respiratory distress syndrome, multi-organ failure, and transplant organ rejection.

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

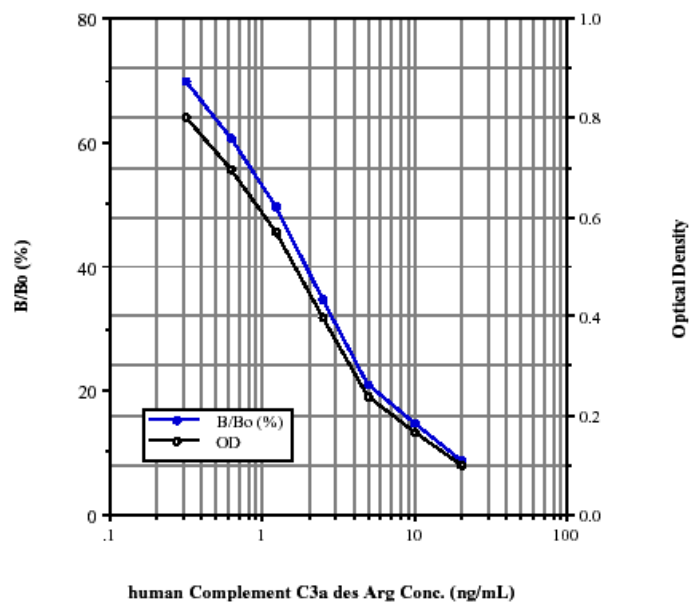
ADI-900-058

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Ready-to-use - sample preparation reagents provided
- Specific - low cross-reactivity to structurally related molecules
- Sensitive - detect as little as 120 pg/ml
- Higher throughput - results from up to 37 samples in duplicate in just 3 hours



## Handling & Storage

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

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of Complement C3a des arg in human EDTA plasma.

**Assay Time** 3 hours

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

**Contents** GxR IgG Microtiter plate, Conjugate, Antibody, Assay buffer 10 concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution, Complement reagent A, Complement reagent B

**Crossreactivity** C3a des Arg (100%), C3 (1.28%), C4a des Arg (0.40%), C4 (0.31%), C5a des Arg (0.10%), C5 (0.05%)

**Sensitivity** 0.120 ng/ml (range 0.313 – 20 ng/ml)

**Species Reactivity** Human

**UniProt ID** P01024

**Wavelength** 405 nm



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-  
fr@enzolifesciences.com](mailto:info-fr@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)