## Complement C3a des Arg (human), ELISA kit

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

**The Complement Grades 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 »

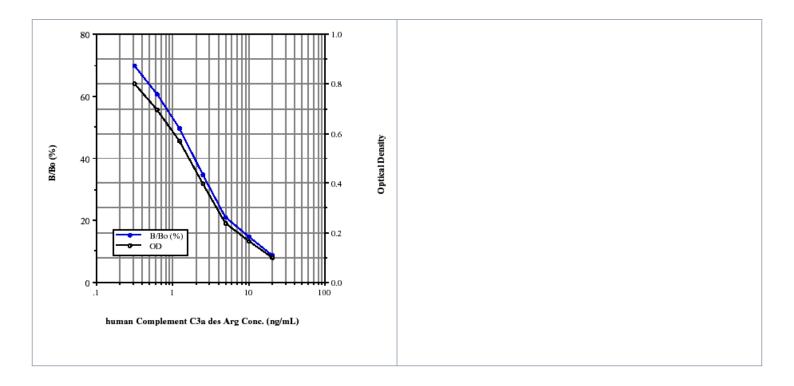
ADI-900-058

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com