## Complement component C1QA monoclonal antibody (C1QA/2953)

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

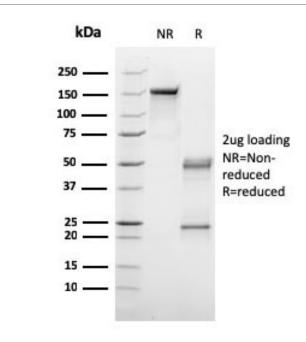
Order Online »

**ENZ-ABS872-0100** 

100µg

Manuals, SDS & CofA

**View Online »** 



SDS-PAGE analysis of purified, BSA-free C1QA antibody as confirmation of integrity and purity.

## **Handling & Storage**

Handling Store the C1QA antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder

(without azide).

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Complement C1q A-Chain

Application IHC (PS), WB

**Clone** C1QA/2953

Formulation Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA

(US sourced), 0.05% sodium azide.

**Host** Mouse

**Immunogen** Recombinant human C1QA (aa 104-237).

lgG2bk

Purity Detail Protein G affinity chromatography purified.

Recommendation Dilutions/Conditions Immunohistochemistry (FFPE) (1-2µg/ml)Western blot (1-

2μg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

UniProt ID P02745

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: July 12, 2024



info-

eu@enzolifesciences.com