

# Complement component C9 (human) monoclonal antibody (C9 53)

Complement factor C9 is a 71kDa single-chain ellipsoid  $\alpha$ -globular molecule of 558 amino acids including a leader sequence of 20 amino acids. C9 binds the C5b8 complex on membranes and has the unique function of forming tubular structures containing between 12 and 18 monomers, closely resembling the membrane attack complex (MAC). The primary site of synthesis is hepatocytes; secondary sites are monocytes, fibroblasts and glial cells. The normal level of C9 in plasma is 60  $\mu$ g/mL, and the level is up regulated from hepatocytes by IFN- $\gamma$ .

This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

|                      |   |
|----------------------|---|
| Alternative Name     | C9 esterase   |
| Application          | ELISA   |
| Clone                | C9 53   |
| Formulation          | Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide. |
| GenBank ID           | 735   |
| Host                 | Mouse   |
| Immunogen            | Human complement factor C9.   |
| Isotype              | IgG1κ   |
| Purity Detail        | Protein A/G-purified.   |
| Source               | From culture supernatant.   |
| Species Reactivity   | Human   |
| UniProt ID           | P02748  |
| Worry-free Guarantee | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>                              |

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)