

Complement component C3 (human) monoclonal antibody (F1-8)

C3 is a central component of the complement system, consisting of a beta-chain (75 kDa) and an alpha-chain (110 kDa) connected by cysteine bridges. Cleavage of C3 into C3b (177 kDa) and C3a (9 kDa) is a pivotal step in the complement activation cascade, which can be initiated by three distinct pathways—the classical, the lectin and the alternative complement pathway.

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

Application	ELISA, Immunoaffinity chromatography, IP
Clone	F1-8
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Complement component 3 (C3) purified from human plasma.
Isotype	IgG1k
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
Specificity	Specific for human complement component 3 (C3/C3b/C3c).
UniProt ID	P01024
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com