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

C3 is a central component of the complement system, consisting of a betachain (75 kDa) and an alpha-chain (110 kDa) connected by cystein 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.

# **Ordering Information**

**Order Online** »

BPD-CAM-008-01-1	1mg
BPD-CAM-008-01-04	400μg

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 IgG1ĸ

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



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
infouk@enzolifesciences.com