

Complement factor H (human) monoclonal antibody (L20/3) (biotin conjugate)

Factor H is a regulatory factor of the alternative pathway of complement activation. Factor H is a glycoprotein consisting of a single polypeptide chain with a molecular mass of 150kDa.

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	Factor H
Application	ELISA
Clone	L20/3
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
GenBank ID	3075
Host	Mouse
Immunogen	Human complement factor H.
Isotype	IgG1κ
Species Reactivity	Human
Specificity	Specific for complement factor H in human serum and plasma.
UniProt ID	P08603
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