

Ficolin H (human) monoclonal antibody (FCN334) (biotin conjugate)

H-Ficolin (Hakata antigen, Ficolin-3) is an innate immunity pattern recognition molecule found in human serum. H-Ficolin binds to distinct pathogen-associated molecular patterns (PAMP) formed by carbohydrates (GlcNAc, GalNAc and fucose) activating complement through the lectin pathway by means of the associated MASPs ((MBL)-associated serine proteases). H-Ficolin consists of 35kDa protein chains associated in triplet subunits, which are again associated in high molecular weight oligomers by means of disulfide bonds at the N-terminal region.

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

Ordering Information

[Order Online »](#)

BPD-RIG-334-01B-015	150µg
---------------------	-------

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	Hakata antigen, Ficolin-3
Application	ELISA
Clone	FCN334
Crossreactivity	No cross-reaction has been observed with L- or M-ficolin.
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
GenBank ID	8547
Host	Mouse
Immunogen	Recombinant human H-ficolin.
Isotype	IgG1k
Species Reactivity	Human
UniProt ID	O75636
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](mailto:info-usa@enzolifesciences.com)

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