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.

Ordering Information

Order Online »

BPD-GAU-020-03B-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 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 IgG1ĸ

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 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com