

Flavivirus group antigen Recombinant monoclonal antibody (D1-4G2-4-15 (4G2)) (Mouse IgG2ak)

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

Ordering Information

[Order Online »](#)

ENZ-ABS492-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Protein E, Envelope Protein
Application	ELISA, Flow Cytometry, IF, IHC (PS), WB
Clone	D1-4G2-4-15 (4G2)
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Dengue Virus type 2 antigens.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Virus
Specificity	Recognizes flavivirus group specific antigens (Dengue virus, West Nile Virus, Japanese Encephalitis, Yellow Fever Virus, Zika virus etc). Binds to the fusion loop at the extremity of domain II of protein E.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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-
eu@enzolifesciences.com](mailto:info-eu@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)