

Zika Virus VLP monoclonal antibody (ID5)

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

Ordering Information

[Order Online »](#)

ENZ-ABS838-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, WB
Clone	IG3-1-A12-G9
Crossreactivity	It also shows cross-reactivity with VLPs and envelope proteins from Dengue virus (DENV), West Nile virus, Japanese Encephalitis virus and Chikungunya virus. It recognises recombinant His-tagged DENV envelope proteins produced in mammalian and insect cells. Fc-tagged envelope proteins may not be detected in ELISA. Antibody detects ZIKV Env and VLP in non-reduced western blot, but not in reduced conditions.
Formulation	Liquid. In PBS, pH 7.4.
Host	Mouse
Isotype	IgG
Purity Detail	Protein G affinity purified.
Source	Purified from hybridoma tissue culture supernatant.
Specificity	Recognizes Zika virus VLPs and envelope protein.
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee