α-Synuclein (human) monoclonal antibody (LB509)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS218-0500	0.5ml
ENZ-ABS218-0200	200μΙ

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, IHC, WB

Clone LB509

Crossreactivity Does not cross-react with rodent α -synuclein.

Formulation Liquid. In PBS. Does not contain any preservatives or

carrier proteins.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 115-122 of human

α-synuclein.

Isotype IgG1ĸ

Purity Detail Purified.

Species Reactivity Human

UniProt ID P37840

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

