

# Amyloid $\beta$ Recombinant monoclonal antibody (WO2) (Mouse IgG2ak)

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

## Ordering Information

[Order Online »](#)

ENZ-ABS386-0100	100µg
-----------------	-------

## 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

Application	ELISA, IHC (PS), WB
Clone	WO2
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Synthetic peptides, corresponding to aa 1-42 of human beta amyloid, conjugated to KLH.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Specificity	Recognizes aa 4-10 of human Amyloid $\beta$ .
UniProt ID	P05067
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>