

**Purified Anti-Human Beta-Amyloid
Antibody *bmAbetaN***

Catalog number: V101140

Unit size: 100 µg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	Amyloid-beta (A β or Abeta, APP, AAA, ABETA, ABPP, AD1, APPI, CTFgamma, CVAP, PN-II, PN2, amyloid beta precursor protein, preA4)
Clone	bm-AbetaN

Biological Properties

Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS)

Applications
