

## Data Sheet

### Anti-β-Amyloid, Plaque Neutralizing Antibody

Human, Recombinant, Monoclonal

**Catalog #:** 71223

**Lot #:** 191121

**Conc.:** 5.54 mg/ml

**Host Species | Isotype:**

Human | IgG1

**Formulated in:** 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Endotoxin Level:** <1 EU/μg

**Stability:** Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

**References:**

1. Doody, R., *et al.*, *New Eng. J. Med.* 2014; **370**: 311-321.
2. Farlow. M., *et al.*, *Science Direct.* 2012; **8(4)**: 261-271.

**Description:**

Purified, recombinant, humanized Anti-β-Amyloid, Plaque Neutralizing (IgG1) antibody was expressed in HEK293 cells. Heavy chain MW = 51 kDa; light chain MW = 26 kDa. This antibody reacts with human β-amyloid and has not been tested for cross-reactivity with other species.

**Purification:**

Protein A affinity chromatography from HEK293 supernatants.

**Applications:**

Useful as a reference antibody for screening β-Amyloid neutralizing antibodies.

### Quality Assurance

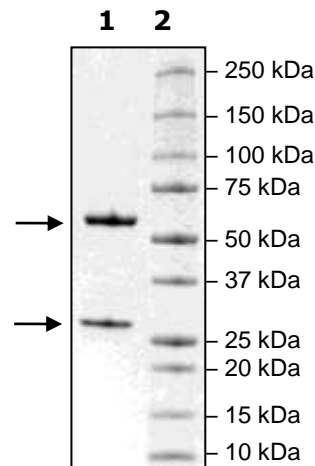
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
5 μg Anti-β-Amyloid

**Lane 2:**  
Protein Marker

**MW:**  
Heavy chain 51 kDa +  
glycans  
Light chain 26 kDa

**Purity:** >90%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)