

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, also known as Anti-Solanezumab, 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-Solanezumab

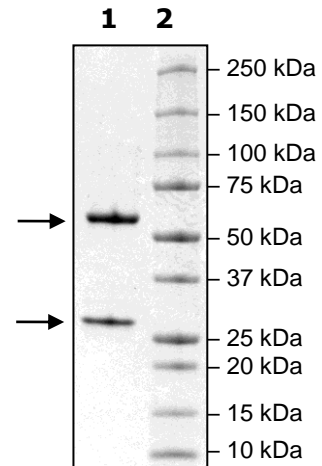
Lane 2:

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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com