

Data Sheet

Anti-CD20, Agonist Antibody

Recombinant, Human Monoclonal Antibody
Anti-CD20

Catalog #: 71209

Lot#: 190122

Conc.: 2.75 mg/ml

Host Species | Isotype:

Murine/Human | IgG1k

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

Endotoxin level: < 0.05 EU/ μ g

Stability: Stable for at least 6 months at -80°C.
Avoid freeze/thaw cycles.

References:

1. Ganjoo KN, *et al.*, *Leuk Lymphoma*. 2014 April 9..
2. Siegal T., *J Clin Neurosci*. 2014 April;**21(5)**:709-715.

Description: Murine/human chimeric Anti-CD20 monoclonal antibody which features the constant region of the human IgG1 isotype and the variable region of rituximab, a.a 142-184, expressed in a HEK293 cell expression system, MW = 51 kDa (heavy chain), MW = 25 kDa (light chain).

EC50: 0.24 μ g/ml

Assay Conditions: WIL2-S cells were treated with anti-CD20 + 10% rabbit complement. Cell viability was determined using AlamarBlue® Indicator.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for the detection of CD20 and binding to B cells or other cells expressing CD20 on their surface.

Quality Assurance

Specific Activity

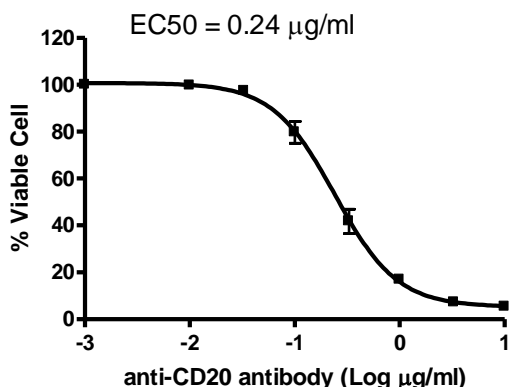


Figure 1. WIL2-S cells treated with various concentrations of CD20 Agonist Antibody (BPS Cat. #71209) + 10% rabbit complement. WIL2-S cell viability was measured using AlamarBlue® Indicator.

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