

## Data Sheet

### Anti-CD3 Agonist Antibody

Recombinant, Human Monoclonal

Catalog #: 71274-1

Lot #: 161122                      Conc.: 3.04 mg/ml

#### Host Species | Isotype:

Human | IgG1

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

**Stability:** Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles. Store diluted antibody with carrier protein (BSA 0.1 – 0.5%) at 4°C, stable for 2 weeks.

#### References:

1. Chatenoud, L., *et al.* PNAS. 1994; **91(1)**: 123-127.
2. Herold, K., *et al.*, NEJM. 2002; **346**: 1692-1698.

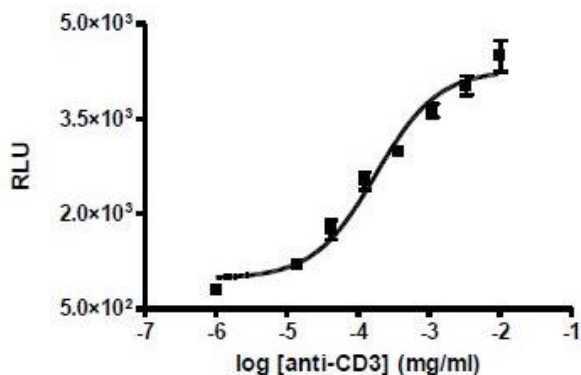
**Description:** Recombinant human antibody recognizing the CD3 binding region on immune cells.

**Purification:** Protein A affinity chromatography from HEK293 supernatants.

**Application:** Useful for the detection of CD3 and binding to immune cells expressing CD3 on their surface and useful as a functional grade antibody for T cell activation and expansion.

## Quality Assurance

### Specific Activity



**Figure 1.** IL-2 luciferase Jurkat reporter cells (BPS #60481) were incubated with anti-CD3 (BPS #71274) in SF-RPMI for 24 hours. BPS luciferase reagent was added and luminescence was quantified. RLU = Relative Luminescence Unit normalized to wells with RPMI only.

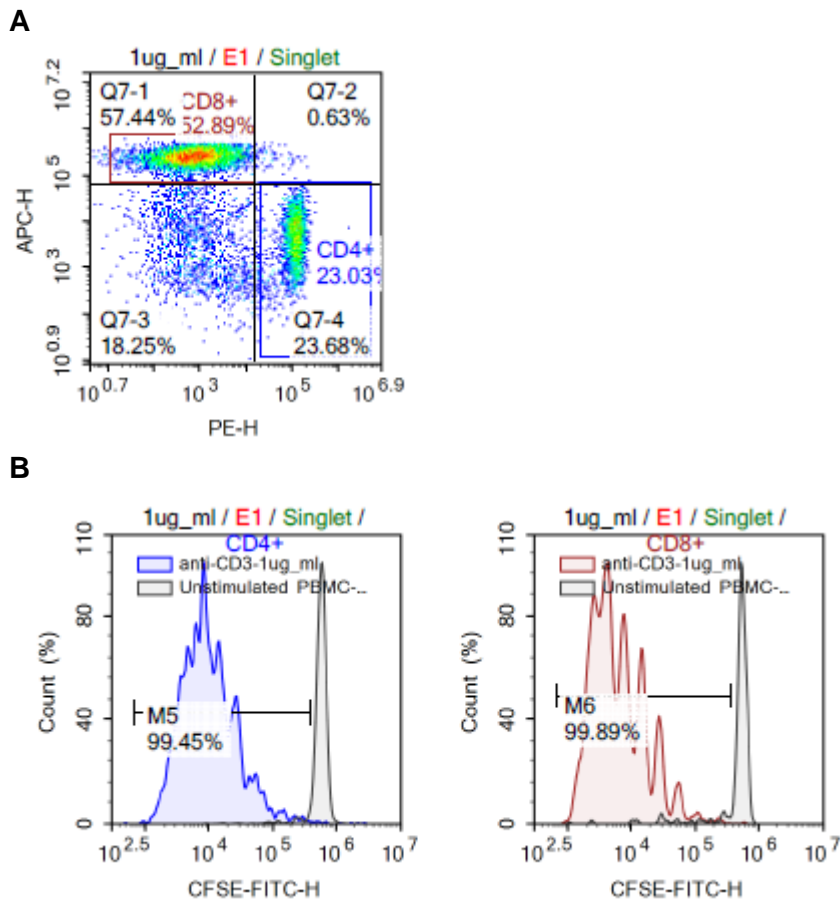
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

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

### Anti-CD3 antibody Functional Study Validation



**Figure 2. T cell division and expansion in response to anti-CD3 antibody**

PBMC (BPS Catalog # ) were treated with 1ug/ml of anti-CD3 antibody (BPS Catalog #71274-2) for 5 days in LGM-3™ medium (Lonza Catalog # CC-3211) supplemented with 10% FBS and hIL-2. CD4+ and CD8+ populations were identified (A) and level of cell proliferation is assessed using CellTrace™ CFSE Cell Proliferation kit (Thermo Fisher Catalog #C34554). Blue = proliferated CD4+ and red = proliferated CD8+ cells. Unstimulated controls are shown in gray.

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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)