

## Data Sheet

### CD19, Fc-Fusion (IgG1), Avi-Tag

Human, Recombinant, C-terminal Fc-Avi-Tag

**Catalog #:** 79472

**Lot#:** 180221

**Conc.:** 0.5 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

1. Brentjens, R.J., *et al.*, *Science Translational Medicine*. 2013 Mar 2013;**5(177)**: 177.
2. Carter, R.H., *et al.*, *Science*. 1992 Apr 1992;**256(5053)**: 105-107.

**Description:**

Human CD19, also known as B-lymphocyte antigen CD19, CVID3, and Cluster of Differentiation 19, GenBank Accession No. NM\_001770, a.a 21-291, with C-terminal Avi-Tag<sup>TM</sup> fused to the Fc portion of Human IgG1 and expressed in a HEK293 cell expression system. MW = 59 kDa. This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

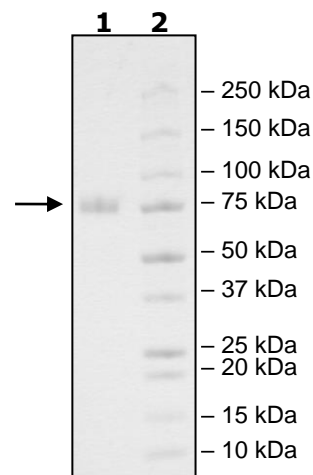
**Application:**

Useful as a positive control for western blot studies.

## **Quality Assurance**

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

**Lane 1:**  
4 $\mu\text{g}$  CD19  
**Lane 2:**  
Protein Marker  
**MW:**  
59 kDa + glycans  
**Purity:**  $\geq 90\%$



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

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