

## Data Sheet

### **KIR2DL3 Fc-fusion, Biotin-labeled**

Human, recombinant, C-terminal Fc fusion  
(Human IgG1)

C-terminal Avi-Tag™

**Catalog #:** 71298

**Lot #:** 160407    **Conc.:** 1.64 mg/ml

**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. Moesta, A, *et al.*, *Immunology*. 2008 Mar; **180(6)**:3969-3979.
2. Maenaka, K., *et al.*, *Structure*. 1999 Jan; **7(4)** : 391-398.

**Description:** Human KIR2DL3, also known as Killer cell immunoglobulin-like receptor 2DL3, KIR2DS5 and KIRCL23. GenBank Accession No. NM\_015868. a.a. 22-245 fused at the C terminus to the Fc portion of human IgG1 followed by a C-termina Avi-tag™ and expressed in a HEK293 cell expression system. MW=53 kDa This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%. This protein runs at a higher MW due to glycosylation.

**Applications:** Useful as an antigen and in protein binding studies.

---

### **Quality Assurance**

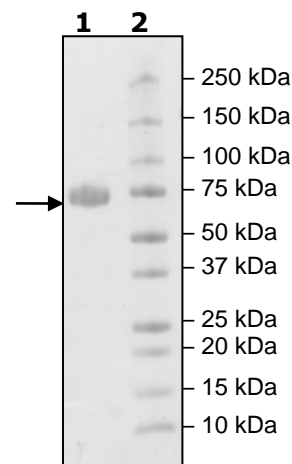
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
3 µg KIR2DL3

**Lane 2:**  
Protein Marker

**MW:** 53 kDa + glycans

**Purity:** ≥90%



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

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

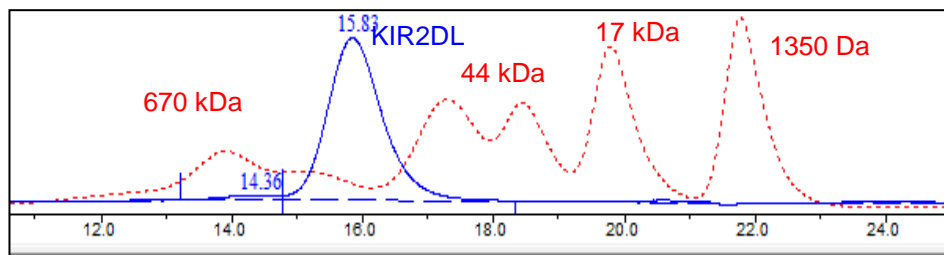
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** info@bpsbioscience.com

### Gel Filtration Curve



< 10% Aggregation

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)