

Data Sheet

CD47, His-tag

Human, Recombinant, C-Terminal His-tag

Catalog #: 71127

Lot #: 151125-G3 **Conc.:** 3.0 mg/ml

Formulated in: 10 mM Phosphate, pH 7.4,
 138 mM NaCl, 2.7 mM KCl

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

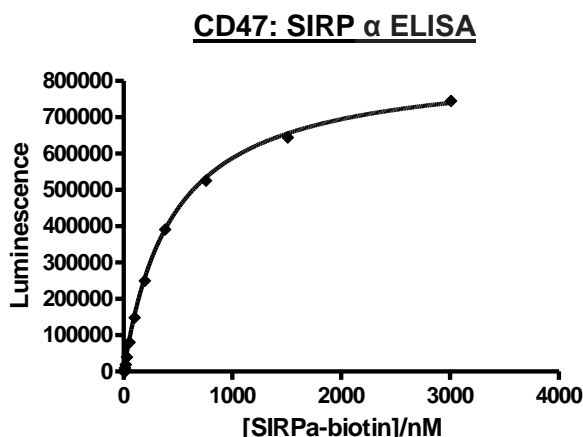
1. Lavender, K.G., et al., *Blood*. 2013 Dec; **122(25)**:4013-20.
2. Zhao, X., et al., *Mol Ther*. 2013 Oct **21(10)**:1811.

Description: Human Recombinant CD47 protein, also known as MER6, with C-terminal His-Tag, GenBank Accession No. NM_001777, a.a. 19-139 expressed in a HEK293 cell expression system. MW = 15 kDa (monomer). This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Assay Description: Coated wells in assay plate with either PBS (No Coat) or 250 ng/well of CD47-His in 50 μl PBS (overnight at 4 deg). Mixed various concentrations of SIRP α -biotin (0-3000nM) in assay buffer and added to assay plate wells. Incubated for 2h at RT. Detected binding with Strep-HRP generated luminescence. Assay Buffer is PBS + 0.1% BSA.

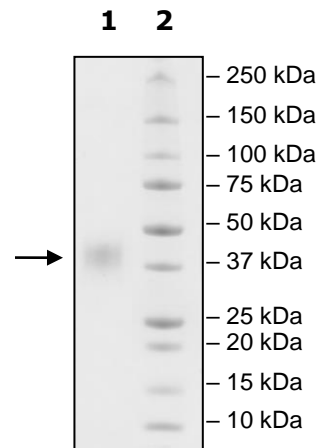
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
 2 μg CD47
Lane 2:
 Protein Marker
MW: 15 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

Please visit our website at: 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

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)