

Data Sheet

CD70, His-Avi-Tag

Human, Recombinant, N-terminal His-Avi-tag

Catalog#: 100240

Lot#: 190220

Conc.: 0.57 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°C . Avoid
 freeze/thaw cycles.

References:

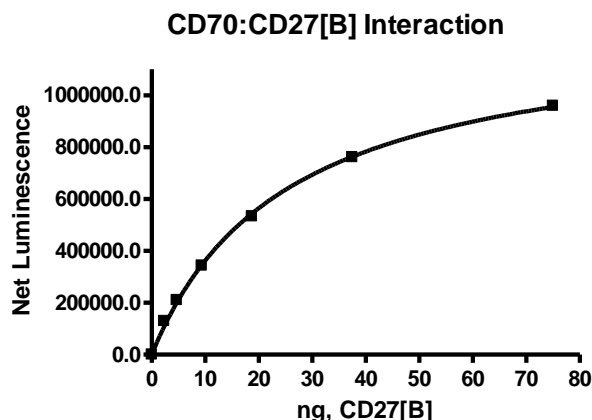
1. Arens, R. *et al.*, *Immunity*. 2001; **15**(5):
 801-812.
2. Law, C.L. *et al.*, *Canc. Res.* 2006; **66**:
 2328-2337.

Description: Human CD70, also known as CD70 antigen, Tumor Necrosis Factor Ligand Superfamily Member 7, TNFSF7, CD27 Ligand, Ki-24 antigen, CD27L, and CD27LG, GenBank Accession No. NM_001252, a.a. 39-193 (end) with N-terminal His-Avi-tag, expressed in a HEK293 cell expression system. MW = 20 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Conditions: Assay was performed according to BPS catalog #79695 with CD27 [B] enzyme titrated from 600 ng/rxn – 0 ng/rxn.

Applications: Useful for studying protein binding and for screening small molecules and antibodies.

Quality Assurance



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

Lane 1:

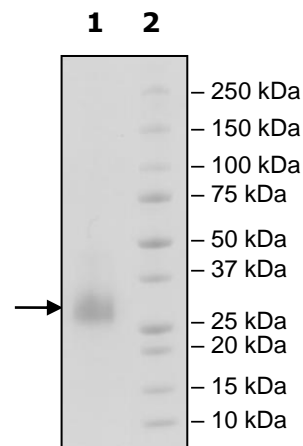
4 μg CD70

Lane 2:

Protein Marker

MW:
 20 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 **1.858.202.1401** 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: 1.858.202.1401
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 **1.858.202.1401** 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)