

Data Sheet

CD70, His-tag

Human, recombinant, N-terminal His-tag

Catalog#: 71178

Lot#: 140716-D **Conc.:** 1.1 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol

Stability: >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 .

Endotoxin Level: <1 EU/ μg

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, CD27-L, and CD27LG, GenBank Accession No. NM_001252, a.a. 39-193 with N-terminal His-tag, expressed in a HEK293 cell expression system. MW = 19.7 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

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

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

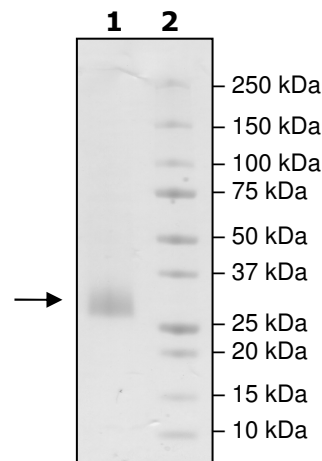
2 μg CD70

Lane 2:

Protein Marker

MW: 19.7 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



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

HHHHHHHHHHGGGSGGGSGGGSSIEGRQRFAQAQQQLPLESLGWDVAELQLNHTGP
QQDPRLYWQGGPALGRSFLHGPELDKGQLRIHRDGIYMVHIQVTLAICSSTTASRHP
TTLAVGICSPASRSISLLRLSFHQGCTIASQRLTPLARGDTLCTNLTGTLPSRNTDET
FGVQWVRP

Black: CD70 (a.a. 39-193(end))

Green: Linker

Red: 10xHis

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)