



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## **Data Sheet**

### **CD137 (4-1BB) HEK293 Recombinant Cell Line**

### **Catalog # 60691**

#### **Background**

Human CD137 (4-1BB; TNFRS9) is an inducible costimulatory molecule that activates T cells. CD137-CD137L mediated signaling has been shown to be important for proliferation, effector functions and survival of T cells. CD137 is also expressed in NK and NKT cells. Antibodies targeting CD137 activation in immune cells have demonstrated potent anti-tumor effects in cancer patients.

#### **Description**

Recombinant HEK293 cell constitutively expressing full length human CD137 (Genbank #NM\_001561). Surface expression is confirmed by flow cytometry.

#### **Host Cell**

Human Embryonic Kidney cell line (HEK293). Adherent epithelial cells.

#### **Format**

Each vial contains ~ 2 x 10<sup>6</sup> cells in 1 ml of 10% DMSO in FBS.

#### **Storage**

Store in liquid nitrogen immediately upon receipt.

#### **Culture Medium**

**Thaw Medium 1** (BPS, #60187): MEM medium (Hyclone, #SH30024.01) supplemented with 10% FBS (Invitrogen, #26140-079), 1% non-essential amino acids (Hyclone, #SH30238.01), 1 mM Na pyruvate (Hyclone, #SH30239.01), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01)

**Growth Medium 1F** (BPS, #79540): Thaw Medium 1 (BPS, #60187) plus 100 µg/ml Hygromycin B (Thermo Fisher, #10687010).

#### **Recommended Culture Condition**

Frozen Cells: Prepare a 50 ml conical tube with 10 ml of pre-warmed Thaw Medium 1 (**no hygromycin**). Quickly thaw cells in a 37°C water bath with constant and slow agitation. Clean the outside of the vial with 70% ethanol and immediately transfer the entire content to Thaw Medium 1 (**no hygromycin**). Avoid pipetting up and down, and gently rock the conical tube.

Spin the cells down at 150 x g for 5 minutes. Discard the medium and re-suspend the cell pellet in fresh Thaw Medium 1 (**no hygromycin**). Transfer the entire content to a T25 flask to distribute the cells. Incubate the cells in a humidified 37°C incubator with 5% CO<sub>2</sub>. After 48-72 hours of incubation, change to fresh Thaw Medium 1 (**no hygromycin**), without disturbing the

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

**Tel:** 1.858.202.1401

**Fax:** 1.858.481.8694

**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

attached cells. Continue to change the medium every 2-3 days until the cells reach desired confluency. If slow cell growth occurs during resuscitation, increase FBS to 15% for the first week of culture. Begin adding Hygromycin B to Thaw Medium 1 (Growth Medium 1F) after the first passage.

*Subculture:* When cells reach 90% confluency, remove the medium and GENTLY wash once with PBS (without Magnesium or Calcium). These cells are loosely adherent and detach easily so do not re-suspend the PBS directly onto the cell surface. Treat cells with 2 ml of 0.25% trypsin/EDTA and incubate for 2-3 minutes at 37°C. After confirming cell detachment by light microscopy, add 10 ml pre-warmed medium and gently pipette up and down to dissociate cell clumps. Transfer cells to a 15 ml conical tube and centrifuge at 200 x g for 5 minutes. Remove the medium and re-suspend cells in 10 ml of pre-warmed Growth Medium 1F. Dispense 5 ml of the cell suspension into a new T75 flask containing 20 ml pre-warmed media. Incubate cells in a humidified 37°C incubator with 5% CO<sub>2</sub>. Freeze cells in freezing medium (10% DMSO in FBS) when cells reach 90% confluency. Cells have been demonstrated to be stable for at least 15 passages; BPS recommends preparing frozen stocks so cells are not used beyond passage 20.

### **Mycoplasma Testing**

This cell line has been screened using the MycoAlert™ Mycoplasma Detection Kit (Lonza, #LT07-118) to confirm the absence of Mycoplasma contamination. MycoAlert Assay Control Set (Lonza, #LT07-518) was used as a positive control.

### **Application**

This cell line is useful for CD137 binding assays or downstream CD137-dependent signaling events.

### **Application References**

1. McNamara II, J.O., *et.al.* (2008) Multivalent 4-1BB binding aptamers co-stimulate CD8+ T cells and inhibit tumor growth in mice. *J. Clin. Invest.* **118**: 376-386.
2. Ma, B.Y., *et.al.* (2005) The expression and regulatory role of OX40 and 4-1BB heterodimers in activated human T cells. *Blood.* **106**: 2002-2010.

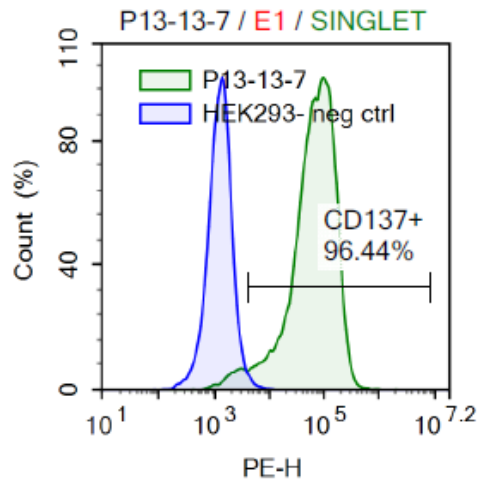
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

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

## Quality Assurance



**Figure 1. Expression of CD137 (4-1BB) protein validated by flow cytometry.** Flow cytometry showed PE-conjugated anti-human CD137 antibody (Clone 4B4-1, Biolegend, #309803) detects CD137 (4-1BB)-positive population (green), using wild-type HEK293 cells as a negative control (blue).

### Vector and Sequence

Human CD137 (NM\_001561) was cloned into pIRESHyg3.

```

MGNSCYNIVATLLLLVLNFERTRSLQDPCSNCPAGTFCDNNRNQICSPCPPNSFSSAGGQRTCD
ICRQCKGVFRTRKECSSTSNAECDCTPGFHCLGAGCSMCEQDCKQGQELTKKGCKDCCFGT
FNDQKRGICRPWTNCSLDGKSVLVNGTKERDVVCGPSPADLSPGASSVTPPAPAREPGHSPQ
IISFFLALTSTALLFLLFLLTLRFSVVKRGRKLLYIFKQPFMRPVQTTQEEDGCSCRFPEEEEEGG
CEL
  
```

**License Disclosure:** Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit the use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use, and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact sales@bpsbioscience.com for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

**Tel:** 1.858.202.1401

**Fax:** 1.858.481.8694

**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

### Related Products

	<u>Cat. #</u>	<u>Size</u>
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
CD137L (4-1BBL) CHO-K1 Cell Line	60523	2 vials
CD137L (4-1BBL) / TCR Activator- CHO-K1 Cell Line	60527	2 vials
CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit	72025	96 rxns.
CD137, Fc fusion (mIgG2a), Biotin-labeled (Mouse) HiP™	71255	50 µg
CD137, Fc fusion (mIgG2a), Avi-tag (Mouse) HiP™	71254	100 µg
CD137, Fc fusion (hIgG1) (Mouse)	71250	100 µg
CD137, Fc Fusion (Human) HiP™	71170	100 µg
CD137, Fc Fusion, Biotin-labeled (Human) HiP™	71171	50 µg
CD137L (4-1BB ligand), His-tag	71189	100 µg
Thaw Medium 1	60187	100 ml

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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