



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

CD137L (4-1BBL) CHO-K1 Recombinant Cell Line

Catalog # 60523

Background

CD137L (4-1BBL; TNLG5A), a part of the tumor necrosis factor (ligand) superfamily (TNFSF9), is a type II glycoprotein that expresses on various antigen-presenting cells. It interacts with CD137 (4-1BB) on T cells following TCR stimulation to promote T cell expansion and cytokine production.

Description

Recombinant CHO-K1 cell constitutively expressing full-length human CD137L (NP_003802.1). Surface expression is confirmed by flow cytometry.

Application

This cell line is useful to stimulate human CD137 (4-1BB) on immune cells, and for use in binding assays for CD137.

Host Cell

Chinese Hamster Ovary Cells. Adherent epithelial cells.

Format

Each vial contains ~ 2 x 10⁶ cells in 1mL of 10% DMSO in FBS.

Storage

Store in liquid nitrogen immediately upon receipt.

Culture Medium

Thaw Medium 3 (BPS Bioscience, #60186): F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin

Growth Medium 3D (BPS Bioscience, #79539): F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 1 mg/ml G418

Recommended Culture Condition

Prepare a 50 ml conical tube and a T-25 culture flask with 5 ml of pre-warmed Thaw Medium 3. 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 contents to the conical tube with Thaw Medium 3 **without G418** and rock the tube the tube gently. Centrifuge the cells at 200 x g for 3 minutes. Re-suspend the cells in 5 ml of pre-warmed Thaw Medium 3 and transfer the entire contents to the T25 culture flask containing Thaw Medium 3 **without G418**. Avoid pipetting up and down, and gently rock the flask to distribute the cells. Incubate the cells in a

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

Please visit our website at: 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

humidified 37°C incubator with 5% CO₂. After 24-48 hours incubation, change to fresh medium without disturbing the attached cells. Continue to change the Thaw Medium 3 **without G418** every 2-3 days until reached desired confluency. If slow cell growth occurs during resuscitation, increase the FBS to 15% during the first week of culture.

Subculture: When cells reached 90% confluency, remove the media and wash twice with PBS (without Magnesium or Calcium). Treat cells with 2-3 ml of 0.25% trypsin/EDTA and incubate for 2-3 minutes at 37°C. After confirming cell detachment by a light microscope, add 10mL of pre-warmed Growth Medium 3D (**with G418**), and gently pipette up and down to dissociate cell clumps. Transfer cells to a 15 ml conical tube and centrifuge at 200xg for 5 minutes. Remove the media and resuspend cells in 10 ml of pre-warmed Growth Medium 3D (**with G418**). Dispense 2 mL of the cell suspension into a new T75 flask containing pre-warmed 18 ml Growth Medium 3D (**with G418**). Incubate cells in a humidified incubator as described above.

Freeze cells in freezing media (10% DMSO in FBS) when they reach 90% confluency. It is not recommended to use the cells after passage 20.

Mycoplasma Testing

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

Application References

1. Yoshimori, M., *et al.* (2014) CD137 Expression Is Induced by Epstein-Barr Virus Infection Through LMP1 in T or NK Cells and Mediates Survival Promoting Signals. *PLoS One* **9**: e112564.
2. Kim, H.J., *et al.* (2012) Reverse Signaling Through the Costimulatory Ligand CD137L in Epithelial Cells is Essential for Natural Killer Cell-Mediated Acute Tissue Inflammation. *Proc. Natl. Acad. Sci.* **109**: E13-E22
3. Salih, H.R., *et al.* (2000) Constitutive Expression of Functional 4-1bb (CD137) Ligand on Carcinoma Cells. *J Immunology* **165**: 2903-2910.

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

Please visit our website at: www.bpsbioscience.com

Quality Assurance

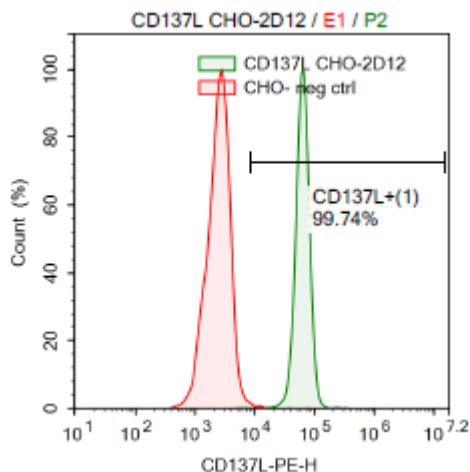


Figure 1. Human CD137L expression in CHO-K1 cells

Flow cytometry demonstrated PE-conjugated anti-human CD137L antibody (Clone 5F4; Biologend, Cat. # 311503) detects CD137L-positive cells (green), but not CHO-K1 wild-type cells used as a negative control (red).

Vector and Sequence

Human CD137L (NM_003811.3) was cloned into the MCS of pIRESneo3 vector (Clontech, Cat. #631621).

AA Sequence

MEYASDASLDPEAPWPPAPRARACRVLPWALVAGLLLLLLLLAAACAVFLACPWAVSGARASP
 GSAASPRLREGPELSPDDPAGLLDLRQGMFAQLVAQNVLIDGPLSWYSDPGLAGVSLTGGLS
 YKEDTKELVVAKAGVYYVFFQLELRVVAGEGSGSVSLALHLQPLRSAAGAAALALTVDLPPAS
 SEARNSAFGFQGRLLHLSAGQRLGVHLHTEARARHAWQLTQGATVLGLFRVTPPEIPAGLPSPR
 SE

License Disclosure:

Visit bpsbioscience.com/license for the label license and other key information about this product.

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

Please visit our website at: 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

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 |
| ONE-Step™ Luciferase Assay System | 60690-3 | 1 L |
| Thaw Media 3 | 60186 | 100 ml |
| Human CD137, Fc fusion (hIgG1) | 71170 | 100 µg |
| Human CD137, Fc fusion (hIgG1), biotinylated | 71171 | 50 µg |
| Mouse CD137, Fc fusion (hIgG1) | 71250 | 100 µg |
| Mouse CD137, Fc fusion (mIgG2a) | 71254 | 100 µg |
| Mouse CD137, Fc fusion (mIgG2a) , biotinylated | 71255 | 50 µg |

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

Please visit our website at: www.bpsbioscience.com