



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

SIRP α / HEK293 Recombinant Cell Line

Catalog # 60689

Background

SIRP α (CD172A, PTPNS1) is a member of the signal-regulatory-protein (SIRP) family, which are transmembrane immunoglobulin receptors that negatively regulate receptor tyrosine kinase-coupled signaling processes. SIRP α is expressed predominantly on macrophages and dendritic cells. Interaction with its ligand CD47 mediates signals that blocks phagocytosis (also known as the "don't eat me" signal).

Description

Recombinant HEK293 cell constitutively expressing full length human SIRP α (#NM_080792). Surface expression is confirmed by flow cytometry.

Host Cell

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

Format

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

Storage

Store in liquid nitrogen immediately upon receipt.

Application

This cell line is useful for SIRP α :CD47 protein binding analyses and screening for antibodies or inhibitors of CD47.

Mycoplasma Testing

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

Application References

1. Lee WY *et.al.* (2010) The Role of *cis* Dimerization of Signal Regulatory Protein α (SIRP α) in Binding to CD47. *J. Biol. Chem.* **285**: 37953-37963

Culture Medium

Thaw Medium 1 (BPS Bioscience #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Thermo Fisher, Cat. #26140079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

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

Growth Medium 1L (BPS Bioscience #79555): Thaw Medium 1 (BPS Bioscience, Cat. #60187) plus 2 µg/ml Puromycin Dihydrochloride (Thermo Fisher, Cat. #A1113803).

Recommended Culture Condition

Frozen Cells: Prepare a 50 ml conical tube with 10 ml of pre-warmed Thaw Medium 1 (**no puromycin**). 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 of the vial into Thaw Medium 1 (no puromycin). Avoid pipetting up and down, and gently rock the conical tube.

Spin the cells down at 150 x g for 5 min. Discard the medium and re-suspend the cell pellet in fresh Thaw Medium 1 (no puromycin). Transfer the entire content to a T25 flask to distribute the cells. Incubate the cells in a humidified 37°C incubator with 5% CO₂. After 48-72 hours of incubation, change to fresh Thaw Medium 1 (**no puromycin**), without disturbing the 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 Growth Medium 1L (**contains Puromycin Dihydrochloride**) 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. 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% CO₂. 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.

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

Quality Assurance

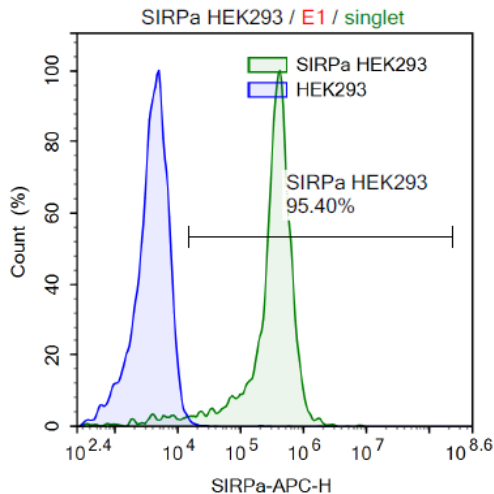


Figure 1. Human SIRP α expression in HEK293 cells

Flow cytometry showed APC-conjugated anti-human CD172a (SIRP α) antibody (Clone REA144; Miltenyi #130-099-785) detects SIRP α on SIRP α / HEK293 cells (green), using HEK293 cells as a negative control (blue).

Vector and Sequence

Human SIRP α (#NM_080792) was cloned into the MCS of pLVX-EF1a-IRES-puro vector (Clontech, #631988).

AA Sequence

```
MEPAGPAPGRLGPLLCLLLAASCAWSGVAGEEELQVIQPDKSVLVAAGETATLRCTATSLIPVG
PIQWFRGAGPGRELIYNQKEGHFPRVTTVSDLTKRNNMDFSIRIGNITPADAGTYCVKFRKGS
PDDVEFKSGAGTELSVRAKPSAPVVS GPAARATPQHTVSFTCESHGFSPRDITLKWFKNGNEL
SDFQTNVDPVGESVSYSIHSTAKVVLTRDVDHSQVICEVAHVTLQGDPLRGTANLSETIRVPPT
LEVTQQPVRAENQVNVTCQVRKFYPQRLQLTWLENGNVSR TETASTVTENKDGTYNWMSWL
LVNVAHRDDVKLTCQVEHDGQPAVSKSHDLKVAHPKEQGSNTAENTGSNERNIYIVGVV
CTLLVALLMAALYLVRIRQKKAQGSTSSTRLHEPEKNAREITQDTNDITYADLNLPKGKKPAPQA
AEPNNHTEYASIQTSPQPA SEDTLTYADLDMVHLNRTPKQPAPKPEPSFSEYASVQVPRK
```

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

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 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 Bioscience products to your laboratory. BPS Bioscience 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.

Related Product

<u>Related Product</u>	<u>Cat. #</u>	<u>Size</u>
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml
SIRP-α (CD172a), His-tag (Human)	71145	100 µg
SIRP-α (SIRP alpha), His-tag, Biotin-labeled (Human)	71138	50 µg
SIRP-γ (CD172g), Fc fusion, Biotin-labeled (Human)	71236	50 µg
CD47, His-tag (Human)	71127	100 µg
CD47, Fc-Fusion, Strep-Tag (Human) HiP™	71292	100 µg
CD47, Fc fusion (Human) HiP™	71177	100 µg
CD47, Fc fusion, Biotin-labeled (Human) HiP™	71169	50 µg
CD47 - HEK293 Cell Line	71249	2 vials
CD47/TCR-Activator CHO-K1 Cell Line	60602	2 vials
CD47:SIRP-α[Biotinylated] Inhibitor Screening Assay Kit	72044	96 rxns.
CD47:SIRP-γ[Biotinylated] Inhibitor Screening Assay Kit	72059	96 rxns.

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