

TREVIGEN[®] Instructions

For Research Use Only. Not For Use In Diagnostic Procedures

Cultrex[®] HA-R-Spondin1-Fc 293T Cells

Cat# 3710-001-01

Stably transfected 293T cells that express murine Rspo1 with an N-terminal HA epitope tag and fused to a C-terminal murine IgG2a Fc fragment.

HA-R-Spondin1-Fc 293T Cells

Stably transfected 293T cells that express murine Rspo1 with an N-terminal HA epitope tag and fused to a C-terminal murine IgG2a Fc fragment.

Cat# 3710-001-01

Table of Contents

Page

I.	Introduction.....	1
II.	Precautions and Limitations	1
III.	Materials Supplied	1
IV.	Materials/Equipment Required But Not Supplied	1
V.	Reagent Preparation.....	3
VI.	Protocol.....	3
VII.	References	6
VIII.	Troubleshooting.....	7
IX.	Related Products Available From Trevigen	8

Legal Notice:

The purchase of this material conveys to the recipient the non-transferable right to use the purchased amount of the product and all replicates and derivatives only for research purposes conducted by the members of the receiving laboratory. The recipient cannot sell or otherwise transfer (a) this product (including replicates and derivatives), (b) its components, or (c) materials made, discovered or identified using this product or its components, to a third party or otherwise use this product (including replicates and derivatives) or its components or materials made using this product or its components for Commercial Purposes. The buyer agrees that any activity undertaken with the product and replicates or derivatives will be conducted in compliance with all applicable guidelines, laws and regulations. Commercial Purposes means any activity by a party for consideration of any kind and includes but is not limited to: (1) use of the product (including replicates and derivatives) or its components in manufacturing; (2) use of the product (including replicates and derivatives) or its components to provide a service, information or data for a third party; (3) use of the product (including replicates and derivatives) or its components for therapeutic, diagnostic or prophylactic purposes; or (4) resale of the product (including replicates and derivatives) or its components, whether or not such product or its components are resold for use in research. "Replicates" means any biological or chemical material that represents a substantially unmodified copy of the material such as, but not limited to, material produced by growth of cells. "Derivatives" means material created from research materials (including replicates, derivatives, conditioned media and purified growth factors) that is substantially modified to have new properties.

This material is understood to be experimental in nature and may have hazardous properties. IT IS BEING SUPPLIED TO RECIPIENT WITH NO WARRANTIES, EXPRESS OR IMPLIED, INCLUDING ANY WARRANTY OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. RECIPIENT ASSUMES ALL RISK AND RESPONSIBILITY IN CONNECTION WITH THE RECEIPT, HANDLING, STORAGE, DISPOSAL, TRANSFER AND USE OF THIS MATERIAL INCLUDING WITHOUT LIMITATION TAKING ALL APPROPRIATE SAFETY AND HANDLING PRECAUTIONS TO MINIMIZE HEALTH OR ENVIRONMENTAL RISK. PURCHASER AGREES THAT ANY ACTIVITY UNDERTAKEN WITH THE MATERIAL AND REPLICATES OR DERIVATIVES WILL BE CONDUCTED IN COMPLIANCE WITH ALL APPLICABLE GUIDELINES, LAWS AND REGULATIONS. TREVIGEN makes no representations that the use of this material will not infringe any patent or proprietary rights of third parties.

I. Introduction

Roof plate-specific Spondin-1 (R-Spondin 1 or RSPO1), also known as CRISTIN3, is a 27 kDa secreted activator protein that belongs to the R-Spondin family. R-Spondins positively regulate Wnt/beta-catenin signaling, most likely by acting as a ligand for LGR4-6 receptors and an inhibitor for ZNRF3. R-Spondin-1 induces proliferation of intestinal crypt epithelial cells, increases intestinal epithelial healing, and supports intestinal epithelial stem cell renewal [1-5]. The 293T cell line is stably transfected to express murine Rspo1 with an N-terminal HA epitope tag and fused to a C-terminal murine IgG2a Fc fragment.

II. Precautions and Limitations

1. Successful and consistent results are dependent upon the quality and degree of characterization of the cells under investigation. Highly passaged cells may undergo both genotypic and phenotypic changes that render them an inadequate *in vitro* model for specific investigations. We recommend for all studies that highly qualified low passage number cells are used to ensure reliable and reproducible results.
2. For Research Use Only. Not for use in diagnostic procedures.
3. This cell line is not known to harbor any agent known to cause disease in healthy adult humans. Handle as a potentially biohazardous material under at least a Biosafety Level 1 containment. This cell line has NOT been screened for Hepatitis B, human immunodeficiency viruses or other adventitious agents. Trevigen recommends that appropriate safety procedures be used when handling all cell lines, especially those derived from human or other primate material. Trevigen assumes no liability for damage resulting from handling or contact with these products.

III. Materials Supplied

<u>Component</u>	<u>Quantity</u>	<u>Storage</u>	<u>Catalog #</u>
HA-R-Spondin1-Fc 293T Cells	1 Vial (10 ⁶ Cells)	Liquid Nitrogen***	3710-001-01

***Shipped on Dry Ice, immediately thaw for use, or for long term storage place in vapor phase of liquid nitrogen.

IV. Materials/Equipment Required But Not Supplied

Equipment

1. 1 - 20 µl, 20 - 200 µl, and 200 - 1000 µl pipettors
2. Class 2 Biosafety Hood
3. 37°C CO₂ incubator
4. 37°C Water Bath
5. Hemocytometer or other means to count cells

6. Inverted standard or phase microscope
7. Pipette aid
8. Liquid Nitrogen Storage
9. Low speed swinging bucket centrifuge and tubes for cell harvesting
10. Cell freezing container that allows for slow freezing of cells
(e.g. Fisher Scientific cat#15-350-50)

Reagents

1. Cell Culture Medium: DMEM High Glucose (Thermo Fisher Scientific cat# 11965-092 or equivalent)
2. Conditioning Medium: Advanced DMEM/F12 (Thermo Fisher Scientific cat# 12634-010)
3. Cell Harvesting Reagent (Trypsin, Dispase, or equivalent)
4. Fetal Bovine Serum
5. 100X Penicillin-Streptomycin supplement for Media
(Thermo Fisher Scientific cat# 15140-122 or equivalent)
6. 100 mg/ml Zeocin (Thermo Fisher Scientific cat# 250-01)
7. GlutaMAX (Thermo Fisher Scientific cat# 35050061)
8. CD293 Medium (Thermo Fisher Scientific cat# 11913-019)
9. PBS (Mg²⁺, Ca²⁺ free) or HBSS, tissue culture grade
10. Trypan blue or equivalent viability stain
11. DMSO, tissue culture grade
12. 70% Ethanol
13. Sterile ddH₂O
14. Protein A Agarose Purification Kit (KPL cat# 553-50-00 or equivalent)

Disposables

1. Cell culture flasks, 25 cm², 75 cm², or 185 cm²
2. 525 cm² 3-layer flasks (Corning cat# 353143)
3. 15 ml tubes
4. 0.22 µm Filter Unit
5. 1 - 200 µl and 200 - 1000 µl pipette tips
6. 2, 5 and 10 ml serological pipettes
7. gloves

V. Reagent Preparation

These procedures should be performed in a biological hood utilizing aseptic technique to prevent contamination.

1. Basal Growth Medium

For 500 ml of Medium:

DMEM Medium	440 ml
Fetal Bovine Serum	50 ml
100X Penicillin-Streptomycin	5 ml
GlutaMAX	5 ml

Filter sterilize medium and store at 4°C for one month. Ensure medium is at room temperature or 37°C prior to use.

2. Selection Growth Medium (containing 300 µg/ml Zeocin)

For 50 ml of Medium:

Basal Growth Medium	50 ml
100 mg/ml Zeocin	150 µl

Prepare immediately before use. Scale as needed.

3. Freeze Medium

For 10 ml:

Basal Growth Medium	5 ml
FBS	3 ml
DMSO	2 ml

4. Expression Medium

For 1000 ml of Medium:

CD293 Medium	1,000 ml
GlutaMAX	10 ml

5. Conditioning Medium

For 500 ml of Medium:

Advanced DMEM/F12 Medium	500 ml
GlutaMAX	5 ml

VI. Protocol

These procedures should be performed in a biological hood utilizing aseptic technique to prevent contamination. Vessels should be sprayed down with 70% ETOH before placing in Tissue Culture Hood.

A. Thawing HA-R-Spondin1-Fc 293T Cells:

1. Pre-warm Basal Growth Medium (Section V.1) to 37°C by placing in H₂O bath or in Tissue Culture Incubator.
2. Immediately before use, remove the vial of cryopreserved cells from liquid nitrogen freezer and thaw quickly (3-4 minutes) in a 37°C H₂O bath. Ensure

cells are completely thawed before proceeding and do not leave cells at 37°C past thawing.

3. Aseptically, transfer the thawed cells to an empty 15 ml conical tube. Wash ampoule with 1 ml of warm Basal Growth Medium and add to thawed cells. Add 1 ml of warm Basal Growth Medium to 15 ml conical tube containing cells, gently swirling to mix between drops. Total Volume should be 3 ml.
4. Centrifuge cells at 200 x g for 3 minutes.
5. Remove supernatant gently to avoid disturbing cell pellet and resuspend cell pellet in 6 ml of fresh Basal Growth Medium.
6. Transfer cell suspension to a sterile T25 (25 cm²) Tissue Culture Flask.
7. Place Tissue Culture Flask/Dish in 5% CO₂ Tissue Culture Incubator at 37°C.
8. Change medium in flasks using freshly prepared and pre-warmed Selection Growth Medium (Section V.2) the next day.

B. Passaging HA-R-Spondin1-Fc 293T Cells:

Note: HA-R-Spondin1-Fc 293T Cells may be sequentially passaged under selection into larger culture vessels to expand cell numbers. We recommend that cultures be maintained at densities between 40 – 90% confluent for optimal growth and survival.

1. Selection Growth Medium should be changed every 2-3 days. Cells should be passaged when 80-90% confluent for optimal growth rate/efficiency. We recommend splitting cells at a density of 1:4 to 1:6. Cells have a doubling time of approximately 24 hrs.
2. Prepare Selection Growth Medium on day of use. Warm Basal Growth Medium to 37°C by placing in 37°C H₂O bath or in Tissue Culture Incubator. In a sterile container, add the required volume (10-12 ml/flask) of Basal Growth Medium and Zeocin (final concentration of 300 µg/ml).
3. Remove medium from T25 flask containing the HA-R-Spondin1-Fc 293T Cells.
4. Gently wash flask with 5 ml of sterile 1X PBS (Ca²⁺ and Mg²⁺ free) and remove PBS.
5. Add 1 ml of warmed Trypsin to the flask and place at 37°C for 3-5 minutes until cells are no longer attached to plate. Add 3 ml of Basal Growth Medium to flask to inactivate Trypsin and transfer to 15 ml conical tube.
6. Centrifuge cells at 200 x g for 3 minutes.
7. Remove supernatant gently to avoid disturbing cell pellet and resuspend cell pellet in 2 ml of fresh Selection Growth Medium.
8. Add 1 ml of cell suspension to 11 ml of Selection Growth Medium and transfer to one T75 flask.

C. Freezing Cells

1. In general, one confluent T75 flask will provide cells for 10-12 vials of 1 x 10⁶ cells per vial.
2. Prepare 2X Freeze Medium, (see Section V.3) according to volume required. Typically, 0.5 ml of 2X Freeze Medium is mixed with 0.5 ml of 2x10⁶ cells.
3. To prepare cells remove medium from T-75 flask containing HA-R-Spondin1-Fc 293T Cells.
4. Gently wash flask with 5-10 ml of sterile 1X PBS (Ca²⁺ and Mg²⁺ free) and remove PBS.

5. Add 2 ml of pre-warmed Trypsin to each flask, and place at 37°C for 2-3 minutes until cells are no longer attached to the plate. Add 4 ml of Basal Growth Medium to flask to inactivate Trypsin and transfer to 15 ml conical tube.
6. Centrifuge cells at 200 x g for 3 minutes.
7. Remove supernatant gently to avoid disturbing cell pellet and resuspend cell pellet in 2 ml of Basal Growth Medium.
8. Count cells on hemocytometer (per standard protocol) and dilute cells to 2 x 10⁶ cells/ml in Basal Growth Medium.
9. Add equal volume of 2X Freeze Medium to the cells, mix gently and aliquot 1 ml of cells into labeled cryovials.
10. For initial storage, first place cryovials on ice for 15-30 minutes. Transfer cells to specialized cell freezing container and place in -80°C freezer overnight.
11. For long term storage, transfer cells (next day) to liquid nitrogen freezer to ensure long-term viability.

D. HA-R-Spondin1-Fc Expression and Purification

1. After at least five days in Selection Growth Medium, HA-R-Spondin1-Fc cells may be expanded in Basal Growth Medium until they reach 80% confluence in the desired culture vessel. We recommend using a 3-layer flask (Corning cat# 353143) to express and purify HA-R-Spondin1-Fc. A 3-layer flask has a surface area of 525 cm², accommodates 150-200 ml of medium, and can produce up to 300 µg of HA-R-Spondin-1-Fc.
2. Once the cells were passaged to the 3-layer flask and reach 80% confluence, change the medium to CD293 medium containing GlutaMAX, and continue to culture for 7-10 days. The cells will detach from the cell culture vessel and grow in suspension after several days.
3. Collect the supernatant, and centrifuge at 3,000 x g for 15 minutes at 4 °C to remove cells and debris.
4. Filter the supernatant through 0.22 µm filter at 4 °C.
5. Purify HA-R-Spondin1-Fc using the Fc tag via Protein A Agarose Purification.

E. HA-R-Spondin1-Fc Conditioned Medium

1. After at least five days in Selection Growth Medium, HA-R-Spondin1-Fc cells may be expanded in Basal Growth Medium until they reach 80% confluence in the desired culture vessel. We recommend using a 3-layer flask (Corning cat# 353143) to produce HA-R-Spondin1-Fc conditioned medium. A 3-layer flask has a surface area of 525 cm², and can produce 150-200 ml of conditioned medium.
2. Once the cells were passaged to the 3-layer flask and reach 80% confluence, change the medium to Advanced DMEM/F12 containing GlutaMAX, and continue to culture for 7-10 days. The cells will detach from the cell culture vessel and grow in suspension after several days.
3. Collect the supernatant, and centrifuge at 3,000 x g for 15 minutes at 4 °C to remove cells and debris.
4. Filter the supernatant through 0.22 µm filter at 4 °C.

5. Verify the presence and titrate HA-R-Spondin1-Fc in the medium by western blot using an anti-R-Spondin1 and/or an anti-HA tag antibody, the expected size should be approximately 70-75 kDa.

Note: If desired, the activity of the conditioned medium can be determined by performing a TopFlash assay (protocol available at www.trevigen.com)

VII. References

1. Ootani, A., et al., *Sustained in vitro intestinal epithelial culture within a Wnt-dependent stem cell niche*. Nat Med, 2009. **15**(6): p. 701-706.
2. Barker, N., et al., *Lgr5+ve Stem Cells Drive Self-Renewal in the Stomach and Build Long-Lived Gastric Units In Vitro*. Cell Stem Cell, 2010. **6**(1): p. 25-36.
3. Sato, T., et al., *Single Lgr5 stem cells build crypt-villus structures in vitro without a mesenchymal niche*. Nature, 2009. **459**(7244): p. 262-265.
4. Sato, T. and H. Clevers, *Growing Self-Organizing Mini-Guts from a Single Intestinal Stem Cell: Mechanism and Applications*. Science, 2013. **340**(6137): p. 1190-1194.
5. Jung, P., et al., *Isolation and in vitro expansion of human colonic stem cells*. Nat Med, 2011. **17**(10): p. 1225-7.

VIII. Troubleshooting

PROBLEM	CAUSE	ACTION
Poor viability from initial freeze	Improper thawing of cells	<p>Ensure medium is added slowly to re-equilibrate the cells from freeze medium</p> <p>Ensure cells were removed from freeze medium immediately after vial has been thawed</p> <p>Ensure vial of cells was thawed at 37°C</p> <p>Fresh medium was warmed to 37°C</p>
Poor proliferation	<p>Fetal Bovine Serum not optimal for cell growth</p> <p>Media not optimal for cell growth</p> <p>Frequency of medium change</p> <p>CO₂ incubator not humidified</p> <p>No gas exchange is allowed by flask</p>	<p>Try alternative lot/source of FBS</p> <p>Ensure medium is of the proper formulation</p> <p>Ensure medium is changed every 2-3 days</p> <p>Ensure pH of medium fresh medium has not changed</p> <p>Add sterile water to CO₂ incubator per manufactures instructions</p> <p>Ensure cap is loosened to allow air gas or use vented flask</p>
Contamination of Cells	<p>Contaminated Medium</p> <p>Improper aseptic technique</p> <p>Hood is working improperly</p> <p>Contaminated CO₂ Incubator</p>	<p>To prevent contamination, filter medium through a 0.22 µm filter before use</p> <p><i>Never use contaminated medium once cloudy or after microorganisms are visible under the microscope</i></p> <p>Spray down hands, reagents and hood with 70% ethanol before opening any flasks</p> <p>Check to make sure blower is on and functioning</p> <p>Ensure hood is currently certified</p> <p>Wipe down hood with 70% ethanol</p> <p>Ensure CO₂ incubator is free of microbial growth</p>
Loss of HA-R-Spondin-1-Fc expression	<p>Not growing cells in Zeocin</p> <p>Frequency of Medium Change</p>	<p>Ensure Zeocin was added to basal medium just before addition to cells</p> <p>Ensure medium is being changed every 2-3 days.</p>

IX. Related Products Available From Trevigen

Related Products:

Catalog#	Description	Size
3750-001-01	Cultrex [®] Organoid Progenitor Cells; Mouse Small Intestine	500 µl
3700-100-01	Cultrex [®] Organoid Harvesting Solution	100 ml
3433-005-R1	Cultrex [®] Reduced Growth Factor BME-R1, PathClear [®]	5 ml
3432-005-01	Cultrex [®] Basement Membrane Extract, PathClear [®]	5 ml
3532-005-02	Cultrex [®] Basement Membrane Extract, Type 2, PathClear [®]	5 ml
3533-005-02	Cultrex [®] Reduced Growth Factor BME, Type 2, PathClear [®]	5 ml
3632-005-02	Cultrex [®] Basement Membrane Extract, Type 3, PathClear [®]	5 ml
3445-005-01	Cultrex [®] 3-D Culture Matrix [™] BME, PathClear [®]	5 ml
3446-005-01	Cultrex [®] 3-D Culture Matrix [™] Laminin I	5 ml
3447-020-01	Cultrex [®] 3-D Culture Matrix [™] Collagen I	100 mg
3434-005-02	Cultrex [®] Stem Cell Qualified RGF BME, PathClear [®]	5 ml
3415-001-03	Cultrex [®] Stem Cell Qualified Human BME, PathClear [®]	1 mg
3400-010-03	Cultrex [®] Stem Cell Qualified Laminin I, PathClear [®]	1 mg
3420-001-03	Cultrex [®] Stem Cell Qualified Human Fibronectin, PathClear [®]	1 mg
3420-001-03	Cultrex [®] Stem Cell Qualified Human Vitronectin, PathClear [®]	200 µg
3400-010-01	Cultrex [®] Mouse Laminin I	1 mg
3400-010-02	Cultrex [®] Mouse Laminin I, PathClear [®]	1 mg
3410-010-01	Cultrex [®] Mouse Collagen IV	1 mg
3440-100-01	Cultrex [®] Rat Collagen I	100 mg
3442-050-01	Cultrex [®] Bovine Collagen I	50 mg
3420-001-01	Cultrex [®] Human Fibronectin, PathClear [®]	1 mg
3416-001-01	Cultrex [®] Bovine Fibronectin, NZHD*	1 mg
3421-001-01	Cultrex [®] Human Vitronectin, PathClear [®]	50 µg
3417-001-01	Cultrex [®] Bovine Vitronectin, NZHD*	50 µg

Related Assays and Kits:

Catalog#	Description	Size
3500-096-K	Cultrex® 3D Spheroid Cell Invasion Assay	96 samples
3510-096-K	Cultrex® 3D Spheroid Fluorometric Proliferation/Viability Assay	96 samples
3511-096-K	Cultrex® 3D Spheroid Colorimetric Proliferation/Viability Assay	96 samples
3470-096-K	Cultrex® In Vitro Angiogenesis Assay, Tube Formation Kit	96 samples
3471-096-K	Cultrex® In Vitro Angiogenesis Assay, Endothelial Cell Invasion Kit	96 samples
3450-048-SK	Cultrex® Directed In Vivo Angiogenesis Assay (DIVAA™) Starter Kit	48 samples
3450-048-K	Cultrex® DIVAA™ Kit	48 samples
3450-048-IK	Cultrex® DIVAA™ Inhibition Kit	48 samples
3465-024-K	Cultrex® 24 well Migration Cell Assay	24 inserts
3455-024-K	Cultrex® 24 well BME Cell Invasion Assay	24 inserts
3456-024-K	Cultrex® 24 well Laminin I Cell Invasion Assay	24 inserts
3457-024-K	Cultrex® 24 well Collagen I Cell Invasion Assay	24 inserts
3458-024-K	Cultrex® 24 well Collagen IV Cell Invasion Assay	24 inserts
3465-096-K	Cultrex® 96 well Migration Cell Assay	96 samples
3455-096-K	Cultrex® 96 well BME Cell Invasion Assay	96 samples
3456-096-K	Cultrex® 96 well Laminin I Cell Invasion Assay	96 samples
3457-096-K	Cultrex® 96 well Collagen I Cell Invasion Assay	96 samples
3458-096-K	Cultrex® 96 well Collagen IV Cell Invasion Assay	96 samples
3448-020-K	Cultrex® 3-D Culture Cell Harvesting Kit	20 samples

*New Zealand Herd Derived

The product accompanying this document is intended for research use only and is not intended for diagnostic purposes or for use in humans.

Trevigen, Inc.

8405 Helgerman Ct. Gaithersburg, MD 20877

Tel: 1-800-873-8443 • 301-216-2800

Fax: 301-560-4973

e-mail: info@trevigen.com

www.trevigen.com



Please
Recycle

P10/20/16v1