

Echelon Biosciences Inc.

Sphingosine 1-Phosphate Assay Kit (S1P – ELISA)

K-1900 (96 tests)

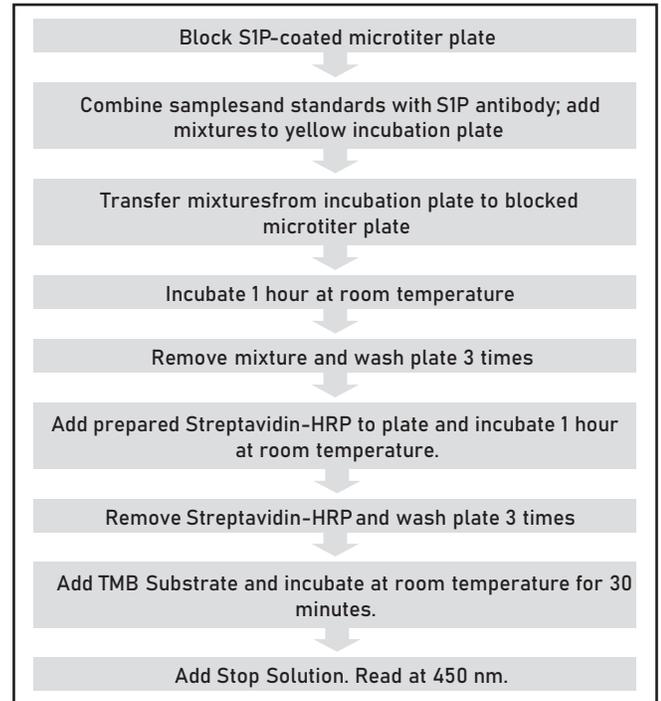
Support: echelon@echelon-inc.com

Description: The S1P ELISA is for the quantification of S1P in biological samples.

Materials Provided

Catalog #	Description	Quantity
K-1901	Anti-S1P Antibody (Clone #LT1002)	Lyophilized powder
K-1903	Block Solution	30 mL
K-1904	Delipidized Serum (DHS)	15 mL
K-1905	Sphingosine 1-Phosphate (S1P) Standard	Lyophilized powder
K-1907	S1P Coated Microtiter Plate (12 X 8 strip-well plate)	1 Plate
K-PTAB	PBS Tablets	4 Tablets
K-PBST	Diluent Buffer	5 mL
K-SEC3h	Streptavidin HRP	80 µL
K-TMB1	TMB Substrate	12 mL
K-STOPt	1N Sulfuric Acid	8 mL
---	Yellow U-bottom Mixing Plate	1 Plate
---	Plate Seals Clear Acetate Sheet, 1 side adhesive	2 Seals

Quick Protocol



Additional Materials Provided by User

- Pipettes (capable of delivering between 5 and 1,000 µL with appropriate tips)
- Multichannel pipettes
- Absorbance microplate reader capable of reading at 450 nm
- Plate Shaker (optional)

Storage Upon receipt, store the S1P-ELISA kit at -20°C. Under proper storage conditions, the kit components are stable for at least 6 months from date of receipt. Opened and reconstituted solutions are less stable. All components and solutions should be protected from excessive light and heat.

Notice to Purchaser: Echelon Biosciences products are sold for research and development purposes only and are not to be incorporated into products for resale without written permission from Echelon Biosciences. This kit is covered by a U.S. Patent and includes a limited, non-transferable license under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. For inquiries email echelon@echelon-inc.com

Echelon Biosciences Inc.

Background

Sphingosine 1-Phosphate (S1P) is a crucial bioactive lipid mediator and key component of the sphingolipid signaling cascade. S1P, produced by the enzymes, Sphingosine Kinase 1 and 2 (SphK1, SphK2), initiates a proliferative¹, pro-angiogenic², and anti-apoptotic³ sequence of events that contributes to the progression of cancer, diabetes, and osteoporosis. SphK1 has been shown to be up-regulated in a variety of cancers and recent research suggests that S1P itself, is a potent tumorigenic growth factor that may be a novel biomarker for early stage cancer detection^{4,5}. S1P levels are tightly regulated by enzymes S1P Lyase (SPL), involved in its degradation, and two S1P-specific phosphatases and three lipid phosphatases that are involved in its synthesis. Additionally, S1P levels and its actions are controlled through a family of cell surface G protein receptors (S1PR1-5). These receptors have been implicated in a variety of developmental and disease related processes; such as immune⁶ and vascular regulation⁷. S1P is also known for its direct role as a second messenger during inflammation⁷ and SPHK/S1P/S1P-receptor signaling has been implicated in a variety of pathophysiological conditions and diseases such as atherosclerosis, cancer, diabetes, multiple sclerosis, sepsis, and so forth. Interference of these routes has the potential for the treatment of cancer, chronic inflammatory disorders and autoimmune diseases.

Assay Design

The S1P-ELISA is an enzyme-linked immunoassay designed for the in vitro measurement of S1P levels in human or animal biological fluids (blood, serum, synovial fluid, cell lysate, and tissue lysate).

The S1P-ELISA is a competitive ELISA in which the colorimetric signal is inversely proportional to the amount of S1P present in the sample. The samples are pre-mixed with the Anti-S1P Antibody while the Microtiter Plate is blocked. Once the block step is complete, the sample/Anti-S1P Antibody mixtures are added to the Blocked Microtiter Plate for competitive binding. Streptavidin-HRP and colorimetric detection is used to detect the Anti-S1P bound to the Microtiter Plate. The concentration of S1P in the sample is determined using a standard curve of known amounts of S1P. This assay should be read at 450 nm and requires 3.5 hours to run.

Assay Range: 0.0625 μ M to 2 μ M

Sensitivity: The lower detection limit of the assay is 0.06 μ M^{13,14}

Sample Volume: For duplicate data points 25 μ L of each sample is required.

Sample Type: This kit detects S1P from human or animal in serum, plasma, tissue homogenate and cell lysate. This assay is not species specific and has been tested with S1P from human, Bovine, equine, caprine, and mouse.

Further Validation Studies: Linearity studies were run by serially diluting samples at dilutions 1:15, 1:20, and 1:30 and comparing observed values with expected values. The observed recoveries ranged from 101%-109%¹³ Recovery of spiked S1P standards was tested by adding 3 different S1P concentrations to eight different human plasma samples with various levels of endogenous S1P. The spiked recovery ranged from 90% to 106%.¹³

Assay Performance

For best results, please follow the protocols provided. Not following the instructions may result in suboptimal performance of the kit and

failure to produce accurate data. All sample types must be diluted in the Delipidized Serum (K-1904) provided in the kit.

Health Hazard Data

The kit contains Delipidized Serum (catalog # K-1904). This reagent should be handled as a potentially bio-hazardous material. The Delipidized Serum was derived from Human blood donors who were individually tested and shown by FDA approved methods to be negative for antibodies to Human Immunodeficiency Virus (HIV) 1/2 and Hepatitis C Virus (HCV), non-reactive to HIV-1, Hepatitis B surface antigen (HBsAg) and Sexually Transmitted Diseases (RPR). Since no test method can offer complete assurance that infectious agents are absent, the Delipidized Serum should be handled following all universal precautions.

S1P-ELISA Assay Notes

1. Ensure samples are free from debris before adding to the plate.
2. All samples and standards MUST be diluted in Delipidized Serum (K-1904).
3. Be cautious of edge effects. The results from wells at the edge of the Microtiter Plate may vary from the interior wells of the plate. If concerned about the edge effect or have observed increased variation with the exterior wells, then do not use the exterior wells of the Microtiter Plate. To reduce edge effects ensure all reagents are at room temperature before use.
4. We recommend heparin or sodium citrate as the anticoagulant used in plasma samples. EDTA interferes with the S1P-ELISA.
5. Lipid cross reactivity was assessed by testing related lipids (DH-S1P, SPH, DH-SPH, CER, CIP, LPA, PAF, SM, PE, PS, DSPA, LPC, PC) at physiologically relevant levels (10 μ M). No cross reactivity was observed except with dihydrosphingosine 1-phosphate (DHS1P) and sphingosylphosphorylcholine (SPC). Until all factors have been tested, the possibility of interference cannot be excluded.
6. **When analyzing biologic samples we advise running a known normal (low) S1P sample and a disease (high) S1P sample in conjunction with your unknown samples. These will serve as positive and negative controls to aid in distinguishing between normal healthy samples and disease samples.**
7. This assay is optimized for detection of S1P in serum and plasma. Sample optimization is highly recommended for other sample types.
8. The S1P-ELISA has been tested with human, bovine, equine and caprine serum sources. The S1P antibody is not species specific. All sample types must be diluted in the Delipidized Serum (K-1904) provided in the kit.
9. Lysis and homogenization buffers may affect the assay. Always test internal lysis and homogenization buffers in the assay for interference before running samples. All samples and standards must be diluted in Delipidized Serum (K-1904) and lysis/homogenate buffers to reduce potential buffer effects.

Protocol for the detection of S1P

Please read this entire section and assay notes before beginning. This protocol has been developed for duplicate reaction points. If singlet or triplicate points are required, the protocol will need to be adjusted accordingly. To begin, place the Streptavidin HRP (K-SEC3h) and Anti-S1P Antibody (K-1901) on ice. Allow the remaining kit components to warm to room temperature before use.

Reagent Preparation

1. **PBS Buffer:** Add 1 PBS Tablet (K-PTAB) to 200 mL DI water. If using plate washer, mix all 4 PBS tablets (K-PTAB) with 800

Echelon Biosciences Inc.

mL DI water. Mix until tablet(s) are completely dissolved. Store prepared PBS Buffer at room temperature.

- S1P Standard Curve** Prepare 100 μM S1P Standard by adding 50 μL DI water to the lyophilized S1P Standard (K-1905). Vortex to mix and place on ice. The S1P standard curve has been designed for serum and plasma samples. If your samples are expected to run outside of this curve, an S1P Standard curve with one standard point higher (4 μM) and/or lower (0.0313 μM) in concentration can be used. Cell lysate and tissue homogenate samples tend to have low S1P; therefore the standard curve (step 2.b) has been adjusted accordingly. The hydrated S1P Standard can handle up to 3 freeze-thaw cycles and should be stored at -20°C when not in use.

a. S1P Standard Curve for the detection of S1P in Serum or Plasma samples Prepare the S1P Standard Curve by adding 12 μL of the 100 μM S1P Standard (K-1905) to 588 μL Delipidized Serum (K-1904) for 2 μM S1P. Then, from the 2 μM S1P make 1:2 serial dilutions using Delipidized Serum to obtain S1P Standards of 1, 0.5, 0.25, 0.125, 0.0625, and 0 μM . These prepared standards can be stored at room temperature for 1-2 hours.

b. S1P Standard Curve for the detection of S1P in Tissue Homogenate or Cell Lysis samples When using cell lysate or tissue homogenate samples a Standard Curve Diluent needs to be prepared. The diluent must contain the same amount of lysis or homogenate buffer as your prepared samples (step 3b; See Assay Notes 2 and 9). The Standard Curve Diluent is used to dilute the S1P Standards and controls. If a smaller or larger dilution is required for your samples the diluent will need to be adjusted as well.

i. Standard Curve Diluent (1 in 10 sample dilution) Prepare by adding 600 μL of Lysis or Homogenate Buffer to 5,400 μL of Delipidized Serum (K-1904).

ii. S1P Standard Curve Prepare by adding 6 μL of the 100 μM S1P Standard (K-1905) to 594 μL Standard Curve Diluent (step 2.b.i) for 1 μM S1P. From the 1 μM S1P, make 1:2 serial dilutions, using the Standard Curve Diluent, to obtain S1P Standards of 0.5, 0.25, 0.125, 0.0625, 0.0313 and 0 μM . These prepared standards can be set at room temperature for 1-2 hours.

3. S1P Sample Preparation

a. Serum and Plasma Samples We recommend a 1:10 dilution of serum and plasma samples in the provided Delipidized Serum (K-1904). For duplicate data points, add 20 μL of sample to 180 μL of Delipidized Serum. This is your working S1P sample.

Important Note: Plasma samples collected in EDTA interfere with the S1P ELISA. We recommend heparin or sodium citrate as the anticoagulant.

b. Cell Lysate or Tissue Homogenate For tissue homogenate and cell culture samples we recommend $> 30 \mu\text{g}$ of total protein/well. Cell lysate and tissue homogenate samples tend to have low amounts of S1P. Therefore, it is highly suggested to determine the amount of total

protein/well needed for your sample before you run your experiments. Concentration of the sample may be required.

i. Tissue Homogenate and Cell Culture Lysate Samples (30 $\mu\text{g}/\text{well}$):

- Dilute the tissue homogenate or cell lysis samples to 4 $\mu\text{g}/\mu\text{L}$ total protein with lysis or homogenate buffer.

- Further dilute the 4 $\mu\text{g}/\mu\text{L}$ lysis sample 1:10 in Delipidized Serum (K-1904) for a 0.4 $\mu\text{g}/\mu\text{L}$ total protein concentration. For duplicate data points, add 20 μL of sample in 180 μL of Delipidized Serum. This is your working S1P sample.

- Anti-S1P Antibody Preparation** Keep the Anti-S1P Antibody (K-1901) on ice. Add 300 μL DI water to rehydrate the lyophilized Anti-S1P Antibody (K-1901). Briefly vortex and place on ice until fully dissolve. Briefly vortex and verify that the pellet is fully dissolved. This stock solution is stable for at least one freeze-thaw cycle and a month's time at -20°C . To prepare Working Anti-S1P Antibody, add 260 μL of the Anti-S1P Antibody stock to 3.5 mL Diluent Buffer (K-PBST). Mix and place on ice until use. **This reagent is not stable at the working concentration and should be prepared immediately before use.**

Working Streptavidin HRP Dilute the Streptavidin HRP

- (K-SEC3h) by adding 20 μL Streptavidin HRP (K-SEC3h) to 12 mL Block Solution (K-1903). Mix well. **Prepare immediately before use.** The Streptavidin HRP (K-SEC3h) can withstand 3 freeze-thaw cycles and should be stored at -20°C . **Save the remaining Block Solution for step 2 in the Assay Procedure.**

Assay Procedure

- Remove S1P-coated Microtiter Plate (K-1907) from plastic bag. Block Microtiter Plate by adding 150 μL of Block Solution (K-1903) to each well. Save the remaining Block Solution (K-1903) for dilution of the Streptavidin HRP (K-SEC3h). Place plate seal on Microtiter Plate and incubate at room temperature for 1 hour. Immediately proceed to the next step.
- Combine your Working S1P Samples or S1P Standards with the Working Anti-S1P Antibody in the Yellow Mixing Plate. This step is written for duplicate data points. For single data points, decrease the volumes listed by half. Use the Yellow Mixing Plate Layout as a guide.
 - Add 60 μL of Working Anti-S1P Antibody to each well of the Yellow Mixing Plate except the Blank control.
 - Add 180 μL of each S1P Standard or Working S1P Sample to the Yellow Mixing Plate according to your plate layout.
 - Carefully tap the plate to mix or place on plate shaker at a moderate speed. Once prepared, the Yellow Mixing Plate is stable for up to 60 minutes at room temperature. Cover with acetate plate sealer if needed.

Only half of the wells in the Yellow Mixing Plate should be utilized. Each well, once mixed, will be transferred in duplicate to the Microtiter Plate. For this step, nothing should be added to the Blank control wells.

- After the 1 hour block step (step 1), remove the Block Solution from the Microtiter Plate and wash 3 times with 200 $\mu\text{L}/\text{well}$



Echelon Biosciences Inc.

Yellow Mixing Plate Layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	2 μ M SIP Standard	-Empty-	Sample # 1	-Empty-	Sample # 9	-Empty-	Sample # 17	-Empty-	Sample # 25	-Empty-	Sample # 33	-Empty-
B	1 μ M SIP Standard	-Empty-	Sample # 2	-Empty-	Sample # 10	-Empty-	Sample # 18	-Empty-	Sample # 26	-Empty-	Sample # 34	-Empty-
C	0.5 μ M SIP Standard	-Empty-	Sample # 3	-Empty-	Sample # 11	-Empty-	Sample # 19	-Empty-	Sample # 27	-Empty-	Sample # 35	-Empty-
D	0.25 μ M SIP Standard	-Empty-	Sample # 4	-Empty-	Sample # 12	-Empty-	Sample # 20	-Empty-	Sample # 28	-Empty-	Sample # 36	-Empty-
E	0.13 μ M SIP Standard	-Empty-	Sample # 5	-Empty-	Sample # 13	-Empty-	Sample # 21	-Empty-	Sample # 29	-Empty-	Sample # 37	-Empty-
F	0.06 μ M SIP Standard	-Empty-	Sample # 6	-Empty-	Sample # 14	-Empty-	Sample # 22	-Empty-	Sample # 30	-Empty-	Sample # 38	-Empty-
G	0 μ M SIP Standard	-Empty-	Sample # 7	-Empty-	Sample # 15	-Empty-	Sample # 23	-Empty-	Sample # 31	-Empty-	Sample # 39	-Empty-
H	Blank	-Empty-	Sample # 8	-Empty-	Sample # 16	-Empty-	Sample # 24	-Empty-	Sample # 32	-Empty-	Sample # 40	-Empty-

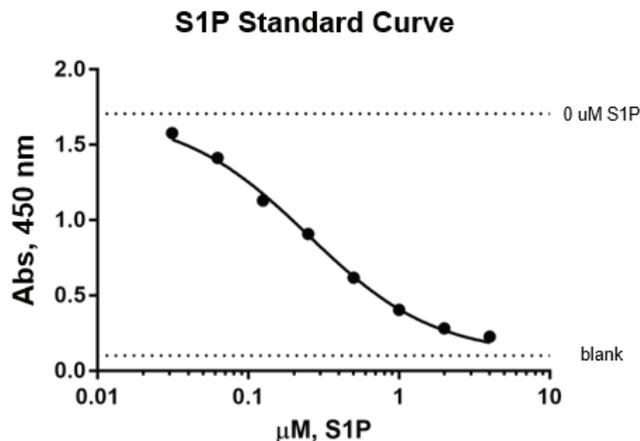
- PBS Buffer. At the end of each wash step, ensure all PBS Buffer is removed from the plate by inverting the plate and blotting it on absorbent paper.
- Transfer 100 μ L of the sample/standard-antibody mixture from the Yellow Mixing Plate, in duplicate, to each well of the Microtiter Plate. Use the Microtiter Plate Layout as a guide. This transfer is easily accomplished with a multi-channel pipette. For best results, prime the pipette 3-6 times before transferring from the Yellow Mixing Plate to the Microtiter Plate. Once the transfers are complete place a plate seal on the Microtiter Plate and incubate at room temperature for 1 hour.

Microtiter Plate Layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	2 μ M SIP Standard	2 μ M SIP Standard	Sample # 1	Sample # 1	Sample # 9	Sample # 9	Sample # 17	Sample # 17	Sample # 25	Sample # 25	Sample # 33	Sample # 33
B	1 μ M SIP Standard	1 μ M SIP Standard	Sample # 2	Sample # 2	Sample # 10	Sample # 10	Sample # 18	Sample # 18	Sample # 26	Sample # 26	Sample # 34	Sample # 34
C	0.5 μ M SIP Standard	0.5 μ M SIP Standard	Sample # 3	Sample # 3	Sample # 11	Sample # 11	Sample # 19	Sample # 19	Sample # 27	Sample # 27	Sample # 35	Sample # 35
D	0.25 μ M SIP Standard	0.25 μ M SIP Standard	Sample # 4	Sample # 4	Sample # 12	Sample # 12	Sample # 20	Sample # 20	Sample # 28	Sample # 28	Sample # 36	Sample # 36
E	0.13 μ M SIP Standard	0.13 μ M SIP Standard	Sample # 5	Sample # 5	Sample # 13	Sample # 13	Sample # 21	Sample # 21	Sample # 29	Sample # 29	Sample # 37	Sample # 37
F	0.06 μ M SIP Standard	0.06 μ M SIP Standard	Sample # 6	Sample # 6	Sample # 14	Sample # 14	Sample # 22	Sample # 22	Sample # 30	Sample # 30	Sample # 38	Sample # 38
G	0 μ M SIP Standard	0 μ M SIP Standard	Sample # 7	Sample # 7	Sample # 15	Sample # 15	Sample # 23	Sample # 23	Sample # 31	Sample # 31	Sample # 39	Sample # 39
H	Blank	Blank	Sample # 8	Sample # 8	Sample # 16	Sample # 16	Sample # 24	Sample # 24	Sample # 32	Sample # 32	Sample # 40	Sample # 40

- Wash the Microtiter Plate 3 times with PBS Buffer (see step 3).
- Add 100 μ L of Working Streptavidin HRP to each well (including the Blank control wells) of the Microtiter Plate. Place a plate seal on the Microtiter Plate and incubate at room temperature for 1 hour.
- Wash the Microtiter Plate 3 times with PBS Buffer (see step 3).
- Add 100 μ L/well of the TMB Substrate (K-TMB1) and incubate for 30 minutes in a dark location. Then add 50 μ L 1N Sulfuric Acid (K-STOPt) to each well of the Microtiter Plate to stop the reaction. Read Microtiter Plate absorbance at 450 nm.

Quantification of Samples:



S1P Standards

μ M SIP	pmol/well SIP
4.0	300
2.0	150
1.0	75
0.5	37.5
0.25	18.75
0.125	9.75
0.0625	4.5
0.03125	2.25

Generate a best fit curve for the S1P standards in order to interpolate relative sample values. The S1P standard curve (below) shows an 8 point S1P curve that was generated using non-linear regression analysis with GraphPad Software. A semi log [Sigmoidal dose-response (variable slope)] analysis was utilized. For best results, constrain the standard curve top & bottom using the 0 μ M S1P & Blank controls. The S1P concentration, for all samples,

should include the dilution factor utilized in step 3 of reagent preparation.

If you are testing tissue homogenate or cell lysate samples the S1P standard curve should be graphed in pmol/well S1P (see table above). The interpolated sample values should then be normalized with grams of total protein or mgs of tissue.

Echelon Biosciences Inc.

Preparation of Plasma

Materials and Equipment

- Human blood sample.
- Vacutainer tubes containing anticoagulant
- Serological pipettes of appropriate volumes (sterile), centrifuge tubes, cryovials
- Benchtop centrifuge (NOT refrigerated) with swing-out rotor and appropriate carriers

Procedure

1. Draw blood into vacutainer tube(s) containing anticoagulant following local standard operating procedures (this may vary depending on manufacturer). Be sure to draw the full volume to ensure the correct blood-to-anticoagulant ratio.
2. Invert vacutainer tubes carefully 10 times to mix blood and anticoagulant and store at room temperature until centrifugation.
3. Samples should undergo centrifugation immediately. This should be carried out for a minimum of 10 minutes at 1000-2000 RCF (generally 1300 RCF) at room temperature (refer to speeds and times recommended by manufacturer). Do not use brake to stop centrifuge.
4. This will give three layers: (from top to bottom) plasma, leucocytes (buffy coat), erythrocytes. Carefully aspirate the supernatant (plasma) at room temperature and transfer to a centrifuge tube. Do not to disrupt the cell layer or transfer any cells.
5. Inspect plasma for turbidity. Turbid samples should be centrifuged and aspirated again to remove remaining insoluble material.
6. Aliquot plasma into cryovials and store at -80°C . Ensure that the cryovials are adequately labeled with the relevant information, including details of additives present in the blood.

Important Note: Plasma samples collected in EDTA interfere with the S1P ELISA. We recommend heparin or sodium citrate as the anticoagulant.

Preparation of Serum

Materials and Equipment

- Human blood sample
- Vacutainer tubes (containing either no additive or a clot activator)
- Serological pipettes of appropriate volumes (sterile), centrifuge tubes, cryovials
- Benchtop centrifuge (NOT refrigerated) with swing-out rotor and appropriate carriers

Procedure

1. Draw whole blood into vacutainer tube(s) containing no anticoagulant following local standard operating procedures. Draw approximately 2½ times the volume needed for use e.g. 10 mL blood for 4 mL serum.
2. Incubate in an upright position at room temperature for 30-45 min (no longer than 60 min) to allow clotting. If using a clot-activator tube, invert carefully 5-6 times to mix clot activator and blood before incubation.
3. Centrifuge for 15 minutes at manufacturer's recommended speed (usually 1,000-2,000 RCF). Do not use brake to stop centrifuge.
4. Carefully aspirate the supernatant (serum) at room temperature and transfer into a centrifuge tube, taking care not to disturb the cell layer or transfer any cells. Use a clean pipette for each tube. Inspect the serum for turbidity. Turbid samples should be centrifuged and aspirated again to remove remaining insoluble matter.
5. Aliquot into cryovials and store at -80°C .

Preparation of Tissue Homogenate

The following protocol was utilized for the determination of S1P concentrations in embryonic and adult kidneys. This protocol was provided by an outside collaborator and has not been validated internally

Procedure

1. Kidneys were homogenized and sonicated in the following homogenization buffer: 20 mM Tris-HCl, pH 7.4; 20% glycerol; 1 mM β -mercaptoethanol; 1 mM EDTA; 1 mM Naorthovanadate; 15 mM NaF; 1 mM PMSF; protease inhibitor cocktail (Sigma); 0.5 mM deoxyypyridoxine; 40 mM β -glycerophosphate.
2. Total protein concentration was measured.
3. The homogenates were then frozen at -80°C until analysis.

Preparation of Cultured Cell Lysate

The S1P ELISA was tested with cell lysates from PC-3, DU-145, and LNCaP cell lines from ATCC. The cell lysate protocol was provided by Dr. Shahriar Koochekpour, School of Medicine, LSU Health Sciences Center, New Orleans, LA.

Analysis Notes

- No cross reactivity with base cell culture media (DME high Glucose, RPMI 1640, SF900 II SFM, BacVector Insect Media, Dulbecco's PBS, HBSS)
- Lipid extraction and sonication do not work – no detectable S1P.
- Media contains FBS. Use of serum free media is recommended

Procedure

1. All reagents should be freshly prepared and all steps should be performed on ice.
2. 75 cm flasks were cultured up to 80-85% confluence, either serum starved for 24 hour or wash with PBS X 2.
3. Cells were lysed in 400 μL of Lysis Buffer with 20 mM PIPES, 150 mM NaCl, 1 mM EGTA, 1% V/V Triton X-100, 1.5 mM MgCl_2 , 0.1% SDS, 1 mM NaOrthovanadate, 1X Protease inhibitor cocktail (without EDTA), pH 7. See Lysis buffer preparation in table below.
4. The clarified lysate was frozen immediately at -80°C . Protein concentration was measured by the BCA method.
5. Diluted cell lysate samples (1:10 in Delipidized Serum, K-1904) are then analyzed with the S1P-ELISA Assay.

Lysis Buffer: Prepare and store at 4°C

Reagent	Stock Concentration	For 100 mL	For 500 mL
PIPES	20 mM	604.8 mg	3.024 g
NaCl	150 mM	876.6 mg	4.383 g
EGTA	1 mM	38.04 mg	190.2 mg
Triton X-100	1% V/V	1 mL	5 mL
MgCl_2	1.5 mM	30.5 mg	152.5 mg
Add immediately before use			
Reagent	Stock Concentration	For 5 mL	
SDS	10%	50 μL	
Protease inhibitor Cocktail	25X	200 μL	
Na-Orthovanadate	1 M	5 μL	



Echelon Biosciences Inc.

Lysis buffer preparation

1. Mix solution with stir bar for 20 minutes and adjust pH to 7.4 using NaOH or HCl until a clear solution is obtained.
2. Filter the solution using a 0.45µm filter. This is stable for 6-12 months and must be stored on ice or 4°C.
3. To make the final "lysis buffer" add 50 µL of 10% SDS, 200 µL of 25X protease inhibitor cocktail and 5 µL 1 M NaOrthovanadate for each 5 mL of working lysis buffer. Vortex mixture and place on ice.

References

General References

1. Dolivera A, Spiegel S. Sphingosine-1-phosphate as a second messenger in cell proliferation induced by PDGF and FCS mitogens. *Nature*. 1993;365:557-560.
2. Anelli V, et al. Role of sphingosine kinase-1 in paracrine/trans-cellular angiogenesis and lymphangiogenesis in vitro. *FASEB J*. 2010;24:2727-2738.
3. Cuvillier O, et al. Suppression of ceramide-mediated programmed cell death by sphingosine-1-phosphate. *Nature*. 1996;381:800-803.
4. Pyne NJ, Pyne S. Sphingosine 1-phosphate and cancer. *Nat Rev Cancer*. 2010;10:489-503.
5. Guan H, et al. Sphingosine Kinase 1 Regulates the Akt/FOXO3a/Bim Pathway and Contributes to Apoptosis Resistance in Glioma Cells. *PLoS One*. 2011;6:e19946.
6. Bao, M., Chen, Z., Xu, Y., Zhao, Y., Zha, R., Huang, S., Liu, L., Chen, T., Li, J., Tu, H., and He, X. (2012) Sphingosine kinase 1 promotes tumour cell migration and invasion via the SIP/EDG1 axis in hepatocellular carcinoma. *Liver international : official journal of the International Association for the Study of the Liver* 32, 331-338
7. Schwalm, S., Pfeilschifter, J., and Huwiler, A. (2013) Sphingosine-1-phosphate: a Janus-faced mediator of fibrotic diseases. *Biochimica et biophysica acta* 1831, 239-250

Anti-SIP Antibody (LT1002) Reference

8. Nicole O'Brien¹, S. Tarran Jones, David G. Williams, H. Brad Cunningham, Kelli Moreno, Barbara Visentin, Angela Gentile, John Vekich, William Shestowsky, Masao Hiraiwa, Rosalia Matteo, Amy Cavalli, Douglas Grotjahn, Maria Grant, Geneviève Hansen, Mary-Ann Campbell, Roger Sabbadini. Production and Characterization of Monoclonal Anti-Sphingosine-1-Phosphate Antibodies *J Lipid Res* 2009, 50, 2245-57.

Selected SIP-ELISA (K-1900) References

Cell culture/cancer cells (human)

9. Olivera A, Spiegel S. Sphingosine-1-phosphate as a second messenger in cell proliferation induced by PDGF and FCS mitogens. *Nature*. 1993, 365:557-560.

SIP in Kidney Homogenates (Results validated by liquid chromatography-tandem mass spectrometry)

10. Kirby, R. J., Y. Jin, et al. (2009). "Dynamic regulation of sphingosine-1-phosphate homeostasis during development of mouse metanephric kidney." *Am J Physiol Renal Physiol* 296(3): F634-641. Lai, W. Q., A. W. Irwan, et al. (2008). "Anti-inflammatory effects of sphingosine kinase modulation in inflammatory arthritis." *J Immunol* 181(11): 8010-8017.

SIP in serum and synovial fluid (Mouse)

11. Cao, C., Y. Gao, et al. (2010). "The efficacy of activated protein C in murine endotoxemia is dependent on integrin CD11b." *J Clin Invest* 120(6): 1971-1980.

SIP in serum and plasma (Human)

12. Watson, L., K. Tullus, et al. (2012). "Increased Serum Concentration of Sphingosine-1-phosphate in Juvenile-onset Systemic Lupus Erythematosus." *J Clin Immunol*.
13. Ardawi M-SM, Rouzi AA, Al-Senani NS, Qari MH, Elsamanoudy AZ, Mousa SA. High Plasma Sphingosine 1-phosphate Levels Predict Osteoporotic Fractures in Postmenopausal Women: The Center of Excellence for Osteoporosis Research Study. *J Bone Metab*. 2018;25(2):87-98.

14. Bae SJ, Lee SH, Ahn SH, Kim HM, Kim BJ, Koh JM. The circulating sphingosine-1-phosphate level predicts incident fracture in postmenopausal women: a 3.5-year follow-up observation study. *Osteoporosis Int*. 2016;1-9. doi: 10.1007/s00198-016-3565-z.

SIP in cell lysate /Media (cell lines: A549,U937)

15. Schnitzer, S. E., A. Weigert, et al. (2009). "Hypoxia enhances sphingosine kinase 2 activity and provokes sphingosine-1-phosphate-mediated chemoresistance in A549 lung cancer cells." *Mol Cancer Res* 7(3): 393-401.
16. Wong, L., S. S. Tan, et al. (2009). "Synthesis and evaluation of sphingosine analogues as inhibitors of sphingosine kinases." *J Med Chem* 52(12): 3618-3626.

Related Products

Products	Catalog Number
Assays and Enzymes	
SIP Activity Assay	K-3500
SIP Inhibitor Screen	K-4400
Sphingosine Kinase 1 and 2	E-K068, E-K069
Antibodies	
Anti-SIP	Z-P300

Products	Catalog Number
Lipids and Beads	
SIP	S-2000
Biotin-SIP	S-200B
SIP coated Beads	S-6110
Agonists	
FYT720	B-0720
FYT720-Phosphate	B-0721

