

MANUAL



AlphaLISA™ SureFire® Ultra™

Part number:	ALSU-XXXX-X500	ALSU-XXXX-X10K	ALSU-XXXX-X50K
Assay points:	500	10,000	50,000

This is a generic manual for all the AlphaLISA™ SureFire® Ultra™ kits.

For assay-specific information, relating to Kit Specificity, Control Lysates and Representative Data, please refer to the Technical Data Sheet of the kit, available from www.revvity.com.

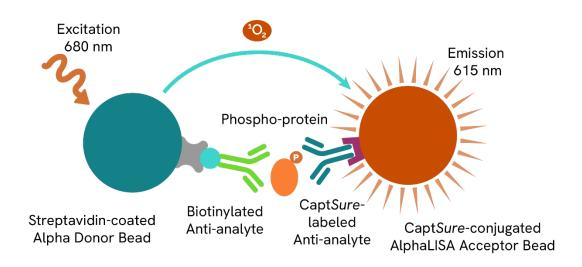
Note: For kit handling and disposal information see page 3 & 4 of this manual

ASSAY PRINCIPLE

The AlphaLISATM SureFire® UltraTM assay kits enable the rapid and sensitive detection of phosphorylated or total cellular proteins. The kit utilizes bead-based Alpha Technology. The Acceptor Bead is coated with a proprietary CaptSureTM agent to specifically immobilize the assay specific antibody which is labeled with a CaptSureTM tag. The Donor Bead is coated with streptavidin to capture the biotinylated antibody.

In the presence of a target protein, the two antibodies bring the Donor and Acceptor Beads in close proximity. This enables the generation of an Alpha signal upon illumination of Donor Beads by the Alpha-enabled plate reader, such as the EnVision™ Multilabel Plate Reader or Enspire™ and EnSight™ Multimode Plate Readers. The amount of light emission is directly proportional to the amount of target protein present in the sample. This assay system performs well in the presence of extraneous antibodies, such as antibody biotherapeutics, and can be used to screen such reagents.

This assay eliminates the need for laborious techniques, such as Western blotting or conventional ELISA. It is a homogeneous assay with no washing steps, minimal handling, short incubation times, enhanced signal-to-noise windows, better well-to-well reproducibility (lower CV%), and robotic operation if desired.



1

MAIN FEATURES

The AlphaLISA™ SureFire® Ultra™ assays are used to measure a phosphorylated or total protein in cellular extracts. The assays are an ideal system for the screening of modulators of receptor activation (e.g agonists and antagonists) as well as agents acting intracellularly, such as small molecule inhibitors of upstream events and also for targeted protein degradation approaches such as PROTAC. The assays measure changes in levels of cellular proteins in endogenous or engineered systems and can be applied to all type of cellular extracts, including cultured primary/non-primary cell lines and tissue lysates.

The kits are very sensitive and formulated to provide improved signal:noise assay windows and to perform without interference in the presence of extraneous antibodies.

KIT-SPECIFICITY / CONTROL LYSATE / REPRESENTATIVE DATA INFORMATION

The assay specific Technical Data Sheet and Certificate of Analysis (COA) are available on the website. Search for Lot Specific COA's from www.revvity.com

KIT CONTENTS

	Kit size			
	500 points	10,000 points	50,000 points	
Lysis Buffer (5X) *	1 x 12 mL	4 x 60 mL	3 x 400 mL	
Supplement** (only in kits with Lysis Buffer B or C (5X))	1 x 1.5 mL	1 x 30 mL	1 x 150 mL	
Activation Buffer *	1 x 0.8 mL	1 x 10 mL	1 x 50 mL	
Reaction Buffer 1 - ALSU	1 x 1.5 mL	1 x 28 mL	1 x 140 mL	
Reaction Buffer 2 - ALSU	1 x 1.5 mL	1 x 28 mL	1 x 140 mL	
Dilution Buffer	1 x 3 mL	1 x 60 mL	1 x 300 mL	
AlphaLISA™ CaptSure™ Acceptor Beads (2 mg/mL in PBS plus 0.03% Proclin-300)	1 x 0.06 mL	1 x 1.1 mL	1 x 5.5 mL	
Alpha Streptavidin Donor Beads (2 mg/mL in PBS plus 0.03% Proclin-300)	1 x 0.06 mL	1 x 1.1 mL	1 x 5.5 mL	
Positive Control Lysate	1 X Lyophilized tube			

The above volumes supplied are in excess to the actual volume required to perform assay.

^{*} Some kits contain assay-specific Lysis Buffer B or C (5X) and Activation Buffer B or C.

^{**} An additional Supplement B or C is provided in kits with Lysis Buffer B or C (5X)only. This supplement must be added to the respective Lysis Buffer B or C (5X) before use. See Storage and Handling Conditions and Buffer Preparation sections for details.

STORAGE AND HANDLING CONDITIONS

Expiry date indicated on kit box

Unopened k	it	Store at 4°C. DO NOT freeze the kit. The Reaction Buffer contains antibodies and freeze/thaw cycles can lead to a loss of activity.
	Lysis Buffer (5X) (including B and C) Reaction Buffer 1 - ALSU Reaction Buffer 2 - ALSU Dilution Buffer	Store at 4°C.
Opened kit	Supplement B or C (only in kits with Lysis Buffer B or C (5X))	Precipitates at 4°C. To re-dissolve, warm to 37°C and mix before each use. Alternatively, once re-dissolved can be stored at room temperature with no loss in activity.
	Activation Buffer (including B and C)	Precipitates at 4°C. To re-dissolve, warm to 37°C and mix before each use. Alternatively, once re-dissolved can be stored at room temperature with no loss in activity.
	Acceptor/Donor Beads	Store at 4°C, in the dark zip lock bag or box provided. Before use, mix well by vortexing vigorously for 5-10 seconds, then give a quick spin (e.g 500rpm for 1 second) to ensure bead suspension is at the bottom of the tube/bottle.
	Positive Control Lysate	Store at -20°C.

MATERIALS REQUIRED BUT NOT PROVIDED

Item	Suggested source	Part #	Size
Optiplate-384, White Opaque assay plate (1)	Revvity Inc.	6007290	50/box
AlphaPlate-384, Light Gray Opaque assay plate (2)	Revvity Inc.	6005350	50/box
CulturPlate-384, White Opaque, Sterile, TC-Treated (3)	Revvity Inc.	6007680	50/box
ViewPlate-384, White with clear bottom, Sterile, TC-Treated (4)	Revvity Inc.	6007480	40/box
White adhesive seal for the bottom of microplates ⁽⁵⁾ .	Revvity Inc.	6005199	1X55
Spectraplate-96, Clear, sterile TC-treated plate (6)	Revvity Inc.	6005650	50/box
TopSeal-A 384, clear adhesive sealing film	Revvity Inc.	6050185	100/box
Envision™ or Ensight™ Alpha-reader	Revvity Inc.	-	-

(1) Plates used for the immunoassay or for the one-plate protocol (from cell seeding to immunoassay) using suspension cells; (2) Same as (1) but optimal if cross-talk needs to be reduced; (3) Plates for assays run in a 1-plate protocol (from cell seeding to immunoassay) using adherent cells; (4) Same as (3) but with the possibility to check cells by microscopy, in this case a white adhesive seal should be stuck to the bottom of the plate before plate reading; (5) This seal can be used to turn the clear bottom of microplates opaque; (6) Plates used to seed and stimulate cells before Lysis and transfer of lysate in an immunoassay plate. For more assay plates options, please go to www.revvity.com.

KIT CONTENT INFORMATION

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are generic for all kits and available from **www.revvity.com**

Lysis Buffer (5X)

Lysis Buffer (5X) is a proprietary mixture of buffers, detergents and generic phosphatase inhibitors (Orthovanadate, Pyrophosphate and sodium fluoride), optimized for lysis of a broad range of cells without the excessive release of nuclear DNA. It does not contain protease inhibitors. Additives can be supplemented to the lysis buffer as required for particular cell types and may include excipients such as protease inhibitors or extra detergents. These will need to be tested on a case-by-case basis.

All Lysis Buffers contain Triton X-100, otherwise known as p-tert-octylphenol ethoxylate, which must be disposed of as Controlled Waste in accordance with Local Regulations.

Lysis Buffers (5X) B and C are assay-specific and <u>must not</u> be inter-changed.

Lysis Buffer color may vary from clear-yellow-green. The visual appearance has no impact on performance.

Supplement B or C

An additional Supplement B or C is provided in kits with Lysis Buffer B (5X) or Lysis Buffer C (5X). This supplement must be added to respective Lysis Buffer B or C (5X) before use.

Activation Buffer

Activation Buffers B and C are assay-specific and must not be interchanged.

Alpha Streptavidin Donor Beads

Alpha Streptavidin Donor Beads are light-sensitive. All Alpha assays using the Donor Beads should be performed under subdued laboratory lighting (< 100 lux). Green filters (LEE 090 filters (preferred) or Roscolux filters #389 from Rosco, or the equivalent) can be applied to light fixtures. The Donor Beads <u>must NOT</u> be used under red/orange light as can be found in photographic work darkrooms because red light (680 nm) excites the beads. All other assay reagents can be used under normal light conditions.

Positive Control lysate

The Positive Control lysates are prepared from various cell types, which have been cultured and prepared to optimize the activation of the intracellular pathway of interest. The lysate is intended for use as an <u>assay positive control only</u> and should not be used for the absolute quantification of a particular protein or phosphorylated target. The lysate can be further diluted with Lysis Buffer (1X) and used to give an indication of the expected signal range for a given assay.

See the COA for the recommended dilution in the linear range of the assay.

BUFFER PREPARATION

Prepare volumes as needed and discard excess

1X Lysis Buffer	Dilute 5X Lysis Buffer 5-fold in deionized water. For example: For 10 mL of 1X Lysis Buffer: Add 2 mL of 5X Lysis Buffer to 8 mL deionized water.	
5X Lysis Buffer B or C	Supplement must be added before use. Dilute Supplement B or C 10-fold in respective 5X Lysis Buffer B or C. For example: For 2 mL of 5X Lysis Buffer B or C: Add 0.2 mL of Supplement B or C to 1.8 mL of respective 5X Lysis Buffer B or C.	
1X Lysis Buffer B or C	Dilute 5X Lysis Buffer B or C (with supplement) 5-fold in deionized water. For example: For 10 mL of 1X Lysis Buffer B or C: Add 2 mL of 5X Lysis Buffer B or C (with supplement) to 8 mL deionized water.	
Acceptor Mix Reaction Buffer 1 + Reaction Buffer 2 + Activation Buffer + Acceptor Beads	Combine Reaction Buffer 1 and Reaction Buffer 2 in equal parts . Dilute Activation Buffer 25-fold in final volume. Dilute Acceptor Beads 50-fold in final volume. For example: For 300 µL of Acceptor Mix: Combine 141 µL of Reaction Buffer 1 and 141 µL of Reaction Buffer 2. Then add 12 µL of Activation Buffer. Then add 6 µL of Acceptor Beads. The Acceptor Mix should be made up and used within 30 minutes for best results.	
Donor Mix Dilution Buffer + Donor Beads	Dilute Donor Beads 50-fold in Dilution Buffer. For example: For 300 μL of Donor Mix: Add 6 μL Donor Beads to 294 μL of Dilution Buffer. The Donor Mix should be made up and used within 30 minutes for best results. Prepare and use under low light conditions.	
Positive Control Lysate	Reconstitute with 250µL deionized water. Reconstituted lysate can generally be stored effectively at -20°C in single use aliquots but should be tested on a case by case basis. Dilute as required with 1X Lysis Buffer.	

AlphaLISA™ SureFire® Ultra™ ASSAY PROTOCOL

A. 2-Plate Assay - assay protocol for adherent cells

Cell Seeding

1. Seed cells (200 μ L of cells for 96 well plates, 50 μ L for 384 well plates) in tissue culture plates. Incubate at 37°C overnight in serum-containing media.

Cell Treatment

2. Remove culture media and replace with 45 μ L of antagonists/inhibitors prepared 1X in <u>serum-free</u> media (22.5 μ L for 384-well plates).

Note: Peptidic agonists and antagonists can often stick to plastic surfaces. To minimize this effect, dilute in serum-free media containing a suitable carrier protein (e.g. 0.1% protease-free BSA).

3. Return cells to incubator at 37°C for 5 min to 2 hours.

Note: 1 hour is often sufficient for signal transduction inhibitors and 5 - 20 minutes for receptor antagonists. For short incubation times, the plate can stay at room temperature.

4. Stimulate the cells by the addition of 5 μ L of 10X agonist prepared in <u>serum-free</u> media (2.5 μ L for 384-well plates). If not testing antagonists, directly add 50 μ L of 1X agonists prepared in <u>serum-free</u> media (25 μ L for 384-well plates).

Note: Optimal agonist stimulation time is often between 5 and 20 minutes.

Lysate Preparation

- 5. To lyse cells, remove medium from wells and add freshly prepared 1X Lysis Buffer (50-100 μ L for a 96 well plate, 20-25 μ L for a 384 well plate). Agitate on a plate shaker (~350 rpm) for 10 minutes at room temperature.
- 6. Take 10 µL of the lysate and transfer to a 384-well Optiplate[™] for assay. Add 10 µL of Control lysates/1X Lysis Buffer to separate wells. We recommend testing a serial dilution of Control lysate in 1X Lysis Buffer. See the COA for recommended dilution in the linear range of assay.

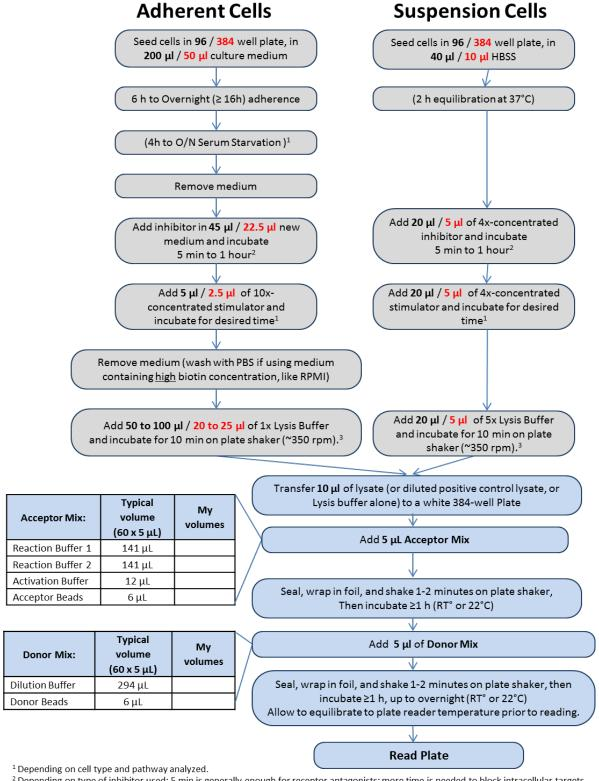
SureFire Ultra Assay

- 7. Add 5 μ L of Acceptor Mix to wells. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature.
- 8. Add $5 \,\mu\text{L}$ of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature in the dark.

Note: Longer incubation may give greater sensitivity. Plates can be incubated overnight if required.

9. Read plate on an Alpha Technology-compatible plate reader, using standard AlphaLISA settings.

AlphaLISA® SureFire® Ultra™: 2-plates / 2-incubation assay flowchart



² Depending on type of inhibitor used: 5 min is generally enough for receptor antagonists; more time is needed to block intracellular targets.

³ May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of lysate before pipetting.

AlphaLISA™ SureFire® Ultra™ ASSAY PROTOCOL

B. <u>1 Plate Assay</u> - assay protocol for suspension cells, and for high-throughput applications.

Cell Seeding

1. Harvest cells by centrifugation and re-suspend cells in HBSS at a suitable cell density. We recommend 10^7 cells/mL as a starting point. Seed 4 μ L of cells/well into a 384-well white opaque culture plate (e.g. Revvity Cat # 6007680).

Note: As engaging less cells per well can result in increased signal to background ratios, it is important to optimize this factor.

Cell Treatment

2. If using test agents/inhibitors, add 2 μ L/well of 3X inhibitors prepared in HBSS.

Note: Peptidic agonists and antagonists can often stick to plastic surfaces. To minimize this effect, dilute in HBSS containing a suitable carrier protein (e.g. 0.1% protease-free BSA).

3. Return cells to incubator at 37°C for 5 min to 2 hours.

Note: 1 hour is often sufficient for signal transduction inhibitors and 5-20 minutes for receptor antagonists. For short incubation times, the plate can stay at room temperature.

4. Stimulate cells with agonists by addition of 2 μ L/well of 4X agonist stock in HBSS containing 0.1% BSA. The final volume in the wells should be 8 μ L. (If no antagonists were used in step 2, stimulate the cells with 4 μ L/well of 2X agonist, to give a final volume in the wells of 8 μ L).

Note: Optimal agonist stimulation time is often between 5 - 20 minutes.

Lysate Preparation

5. To lyse the cells add 2 μL/well of 5X Lysis Buffer. Add 10 μL of Control lysate/ 1X lysis Buffer to separate wells. We recommend testing a serial dilution of Control lysate in 1X Lysis Buffer. See the COA for recommended dilution in the linear range of assay.

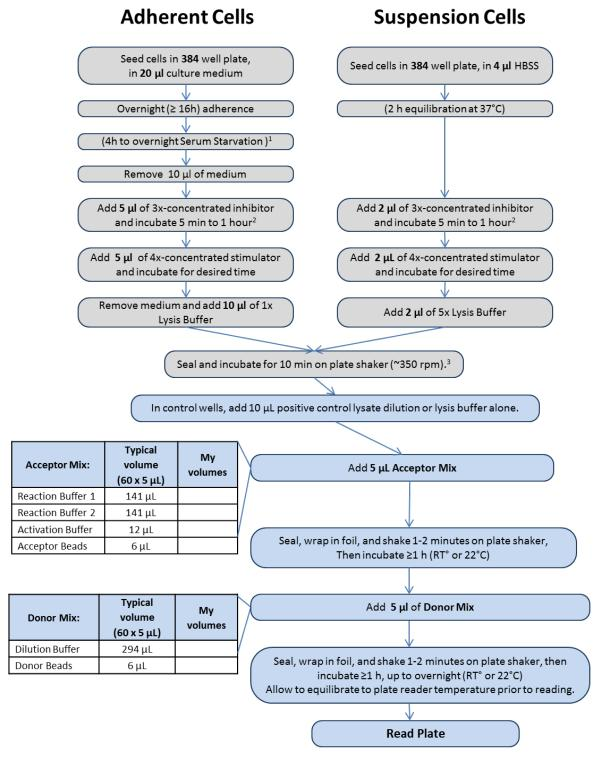
SureFire Ultra Assay

- 6. Add 5 μ L of Acceptor Mix to wells. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature.
- 7. Add 5 μ L of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature **in the dark**.

Note: Longer incubation may give greater sensitivity. Plates can be incubated overnight if required.

8. Read plate on an Alpha Technology-compatible plate reader, using standard AlphaLISA settings.

AlphaLISA® SureFire® Ultra™: 1-plate / 2-incubation assay flowchart



¹ Depending on cell type and pathway analyzed.

² Depending on type of inhibitor used: 5 min is generally enough for receptor antagonists; more time is needed to block intracellular targets.

³ May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of lysate before pipetting.

SUPPLEMENTARY BUFFERS AND BEADS

If using the standard protocol, sufficient amount of buffers and beads are provided in the kit. However, in case the standard protocol would be modified, more buffers or beads may be needed. In this case, you can order additional buffers and beads using the following part numbers:

ltem	Suggested source	Part numbers	Size
	Revvity Inc.	ALSU-LB-10mL	10 mL
Lysis Buffer (5X)	Revvity Inc.	ALSU-LB-100mL	100 mL
	Revvity Inc.	ALSU-LBB-10mL	10 mL
Lysis Buffer B (5X) *	Revvity Inc.	ALSU-LBB-100mL	100 mL
	Revvity Inc.	ALSU-LBC-10mL	10 mL
Lysis Buffer C (5X) *	Revvity Inc.	ALSU-LBC-100mL	100 mL
	Revvity Inc.	ALSU-AB-10mL	10 mL
Activation Buffer	Revvity Inc.	ALSU-AB-100mL	100 mL
	Revvity Inc.	ALSU-ABB-10mL	10 mL
Activation Buffer B	Revvity Inc.	ALSU-ABB-100mL	100 mL
	Revvity Inc.	ALSU-ABC-10mL	10 mL
Activation Buffer C	Revvity Inc. ALSU-ACAB-0.06mL	100 mL	
			10 mL
Dilution Buffer	Revvity Inc. ALSU-ABDB-100mL Revvity Inc. ALSU-DB-100mL Revvity Inc. ALSU-ASDB-0.06mL Revvity Inc. ALSU-ASDB-1.2mL Revvity Inc. ALSU-ACAB-0.06mL Revvity Inc. ALSU-ACAB-0.06mL Revvity Inc. ALSU-ACAB-1.2mL	100 mL	
	Revvity Inc.	ALSU-LB-10mL ALSU-LBB-100mL ALSU-LBB-100mL ALSU-LBB-100mL ALSU-LBC-10mL ALSU-LBC-100mL ALSU-AB-100mL ALSU-AB-100mL ALSU-ABB-100mL ALSU-ABB-100mL ALSU-ABB-100mL ALSU-ABC-10mL ALSU-ABC-10mL ALSU-ABC-10mL ALSU-ABC-10mL ALSU-ABC-10mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ABC-100mL ALSU-ASDB-0.06mL ALSU-ASDB-0.06mL ALSU-ACAB-0.06mL ALSU-ACAB-1.2mL	60 μL
Alpha Streptavidin Donor Beads -2 mg/mL	Revvity Inc.	ALSU-ASDB-1.2mL	1.2 mL
,,	Revvity Inc.	ALSU-ASDB-6mL	6 mL
Revvity Inc.	ALSU-ACAB-0.06mL	60 μL	
AlphaLISA™ CaptSure™ Acceptor Beads - 2 mg/mL	Revvity Inc.	ALSU-ACAB-1.2mL	1.2 mL
2 mg/ m2	Revvity Inc.	ALSU-ACAB-6mL	6 mL

^{*}Includes respective Supplement B or C.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

www.revvity.com



