

**MANUAL**



# Alpha SureFire® Ultra™ HV Multiplex

## Phospho (Eu) + Total (Tb) Target

<b>Part Number</b>	<b>MPSU-XXXX-X-HV</b>
<b>Assay Points</b>	<b>100 (96 well format)</b>

This is a generic manual for the Alpha SureFire® Ultra™ Multiplex HV Phospho + Total kits.

For assay-specific information, relating to Kit Specificity, Control Lysates and Representative Data, please refer to the Technical Data Sheet of the kit, also available from [www.revvity.com](http://www.revvity.com).

**Note: For kit handling and disposal information see pages 4-6 of this manual**

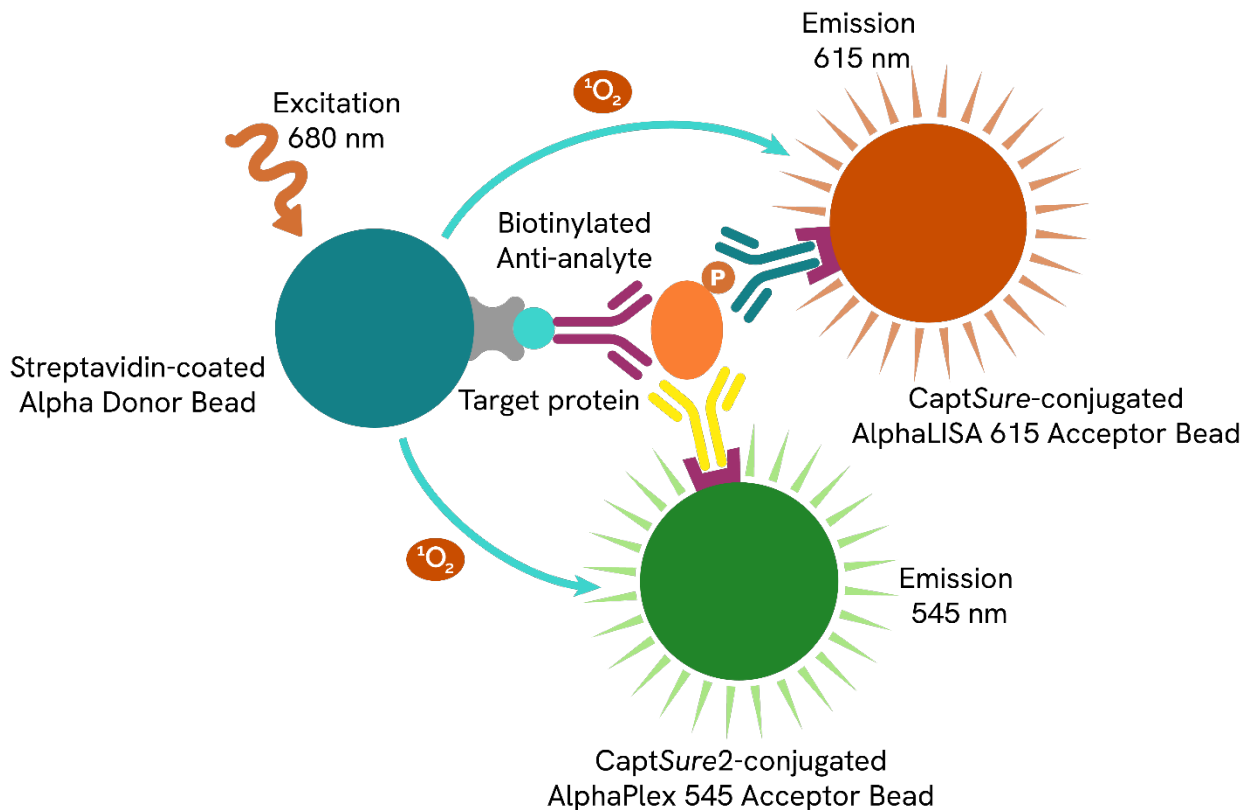
## ASSAY PRINCIPLE

The Alpha SureFire® Ultra™ Multiplex HV Phospho + Total assay kits allow for the rapid and sensitive detection of phosphorylated and total cellular proteins from the same protein. This Alpha Multiplex measurement is carried out in the same assay plate well from a single sample of cell lysate and is achieved by the use of two types of Alpha Acceptor beads that emit at distinct wavelengths (545nm and 615nm).

The two distinct Alpha Acceptor beads report their binding to a distinct antigen through their association with specific assay antibodies, as indicated below.

### Single target - Phospho + Total Assay kits

The Alpha 615 Acceptor Bead is coated with the proprietary CaptSure™ agent, which binds the CaptSure-tagged anti-phospho target antibody. The Alpha 545 Acceptor Bead is coated with the proprietary CaptSure2 agent, which binds the CaptSure2 tagged anti-total target protein antibody. The Donor Bead binds the biotinylated anti-total target protein antibody.



## MAIN FEATURES

The Alpha *SureFire*<sup>®</sup> *Ultra*<sup>™</sup> *HV* Multiplex assay kits are used to measure both the phosphorylation and total levels of endogenous signaling proteins in cellular extracts. The assay is 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 signal transduction. 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 615nm (Eu) signal corresponds to the phosphorylated protein analysis, and the 545nm (Tb) signal corresponds to the total protein analysis.

This kit has been formulated to provide improved signal:noise assay windows, and to perform without interference in the presence of extraneous antibodies.

The AlphaLISA<sup>™</sup> *SureFire*<sup>®</sup> *Ultra*<sup>™</sup> *HV* Multiplex assays are carried out using a higher volume of reagents than the standard AlphaLISA<sup>™</sup> *SureFire*<sup>®</sup> *Ultra*<sup>™</sup> Multiplex assays, and in 96 well plate format rather than 384 well format. These kits are useful for those laboratories where handling of the larger volumes is more appropriate for the laboratory equipment.

The assay utilizes the bead-based Alpha Technology, and requires an Alpha Technology-compatible plate reader capable of reading dual emission wavelengths. See [www.revivity.com](http://www.revivity.com) for more information about the AlphaPlex technology and download the "AlphaPlex Quick Start Guide" and the "AlphaPlex Assay Development Guide" to find guidance about filters and mirrors selection, instrument protocol and channels crosstalk correction. It is to be noted that, as the analytes recognized by both assays (i.e. the phosphorylated protein and the total protein) cannot be dissociated, it is not possible to omit one or the other analyte for the establishment of the channels crosstalk correction, but one or the other type of acceptor beads needs to be omitted instead. i.e. all the assay components but the Alpha 615 beads must be assembled to establish the crosstalk of the Alpha 545 beads into the 615 nm channel, and all the assay components but the Alpha 545 beads must be assembled to establish the crosstalk of the Alpha 615 beads into the 545 nm channel.

## 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.revivity.com](http://www.revivity.com).

## KIT CONTENTS

	100 points
Lysis Buffer (5X)*	1 x 12 mL
Supplement** (only in kits with Lysis Buffer B or C (5X))	1 x 1.5 mL
Activation Buffer *	1 x 0.3 mL
Reaction Buffer 1 - MPSU (Biotinylated anti-Total antibody)	1 x 0.6 mL
Reaction Buffer 2 - MPSU (CaptSure™ tagged anti-phospho antibody)	1 x 0.6 mL
Reaction Buffer 3 - MPSU (CaptSure2™ tagged anti-Total antibody)	1 x 0.6 mL
Dilution Buffer	1 x 1.8 mL
Alpha 615 CaptSure™ Acceptor Beads (2mg/mL in PBS plus 0.03% Proclin-300)	1 x 0.045 mL
Alpha 545 CaptSure2™ Acceptor Beads (2mg/mL in PBS plus 0.03% Proclin-300)	1 x 0.045 mL
Alpha Streptavidin Donor Beads (2mg/mL in PBS plus 0.03% Proclin-300)	1 x 0.045 mL
Positive Control Lysate	1 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.

<b>Unopened kit</b>		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.
<b>Opened kit</b>	Lysis Buffer (5X) (including B and C) Reaction Buffer 1 - MPSU Reaction Buffer 2 - MPSU Reaction Buffer 3 - MPSU Dilution Buffer	Store at 4°C
	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
Half AreaPlate-96, white Opaque assay plate <sup>(1)</sup>	Revvity Inc.	6002290	2x25
Half Area AlphaPlate-96, Light Gray Opaque assay plate <sup>(2)</sup>	Revvity Inc.	6002350	2x25
Half Area AlphaPlate-96, Light Gray Opaque, Sterile, TC-Treated assay plate <sup>(3)</sup>	Revvity Inc.	Custom product	1x50
Half Area ViewPlate-96, White with clear bottom, Sterile, TC-Treated assay plate <sup>(4)</sup>	Revvity Inc.	6005760	1x40
White adhesive seal for the bottom of microplates <sup>(5)</sup> .	Revvity Inc.	6005199	1X55
Spectraplate-96, Clear, sterile TC-treated plate <sup>(6)</sup>	Revvity Inc.	6005650	50/box
TopSeal-A 384, clear adhesive sealing film	Revvity Inc.	6050185	100/box
Envision™ or Ensign™ Alpha-reader with adequate AlphaPlex filters (see table below)	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](http://www.revvity.com).

**Table:** AlphaPlex Optics for EnVision Multilabel Reader – for complete information about how to set an AlphaPlex reading, please refer to the AlphaPlex Guides available at [www.revvity.com](http://www.revvity.com).

	Description	Catalog #	Barcode	Recommendations
Mirrors	AlphaScreen	2101-4010	444	For Tb and Eu single and sequential reading ; not for Sm
	AlphaPlex Single Tb-Eu-Sm	2102-5910	605	Preferred mirror for all sequential AlphaPlex applications
	AlphaPlex Dual Tb-Eu	2102-5900	653	For <b>simultaneous</b> duplexing of Tb with Eu
Filters	AlphaScreen	2100-5710	244	Suitable for AlphaPlex single plexing, not for multiplexing
	Resorufine/Amplex Red	2100-5570	124	Suitable for Tb single plexing and Tb/Eu duplexing.
	Europium	2100-5090	203	Preferred filter for all Eu applications and multiplexing
	AlphaPlex Tb	2100-5930	701	Preferred filter for all Tb applications and multiplexing

TGRSU237.8

SureFire is a registered trademark of TGR BioSciences Pty Ltd, Australia. SureFire Ultra and CaptSure are trademarks of TGR BioSciences Pty Ltd, Australia. Revvity, OptiPlate and AlphaLISA are registered trademarks of Revvity, Inc.

## 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.revivity.com](http://www.revivity.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 must not 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 must NOT 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 assay positive control only 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

<p><b>1X Lysis Buffer</b></p>	<p>Dilute 5X Lysis Buffer <b>5-fold</b> in deionized water.</p> <p>For example: For 10 mL of 1X Lysis Buffer: Add 2 mL of 5X Lysis Buffer to 8 mL deionized water.</p>
<p><b>5X Lysis Buffer B or C</b></p>	<p><b>Supplement must be added before use.</b> Dilute Supplement B or C <b>10-fold</b> in respective 5X Lysis Buffer B or C.</p> <p>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.</p>
<p><b>1X Lysis Buffer B or C</b></p>	<p>Dilute 5X Lysis Buffer B or C (with supplement) <b>5-fold</b> in deionized water.</p> <p>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.</p>
<p><b>Acceptor Mix</b></p> <p>Reaction Buffer 1 + Reaction Buffer 2 + Reaction Buffer 3 + Activation Buffer + 615 Acceptor Beads + 545 Acceptor Beads</p>	<p>Combine Reaction Buffers 1, 2 and 3 in equal parts. Dilute Activation Buffer <b>25-fold</b> in final volume. Dilute each Acceptor bead <b>50-fold</b> in final volume.</p> <p>For example: For 300 µL of Acceptor Mix: Combine 92µL Reaction Buffer 1, 92 µL of Reaction Buffer 2 and 92 µL of Reaction Buffer 3. Then add 12µL Activation Buffer. Then add 6µL 615 Acceptor Bead and 6µL 545 Acceptor Bead.</p> <p>The Acceptor mix should be made up and used within 30min for best results.</p>
<p><b>Donor Mix</b></p> <p>Dilution buffer + Donor Beads</p>	<p>Dilute Donor beads <b>50-fold</b> in Dilution buffer</p> <p>For example: For 300 µL of Donor Mix: Add 6 µL Donor Beads to 294 µL of Dilution Buffer</p> <p>The Donor Mix should be made up and used within 30 minutes for best results. <b>Prepare and use under low light conditions.</b></p>
<p><b>Positive Control Lysate</b></p>	<p>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.</p>



# Alpha SureFire® Ultra™ HV Multiplex PHOSPHO + TOTAL ASSAY PROTOCOL

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

### Cell Seeding

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

### Cell Treatment

2. Remove culture media and stimulate the cells with 50 µL agonists prepared in serum-free media (25 µL for 384-well plates). (*If testing antagonists, prior to stimulation remove culture medium and replace with 50 µL serum-free media containing antagonists (25 µL for 384-well plates)*). Return cells to 37°C incubator for desired time. 1 hour is often sufficient for signal transduction inhibitors, and 5-20 minutes for receptor agonists.

**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% BSA)

### Lysate Preparation

3. To lyse cells, remove medium from wells, and add freshly prepared 1X Lysis Buffer (50- 100 µL for a 96 well plate, 20-25 µL for a 384 well plate). Agitate on a plate shaker (~350 rpm) for 10 minutes at room temperature.
4. Take 30 µL of the lysate and transfer to a 96-well 1/2AreaPlate™ for assay. Add 30 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.*

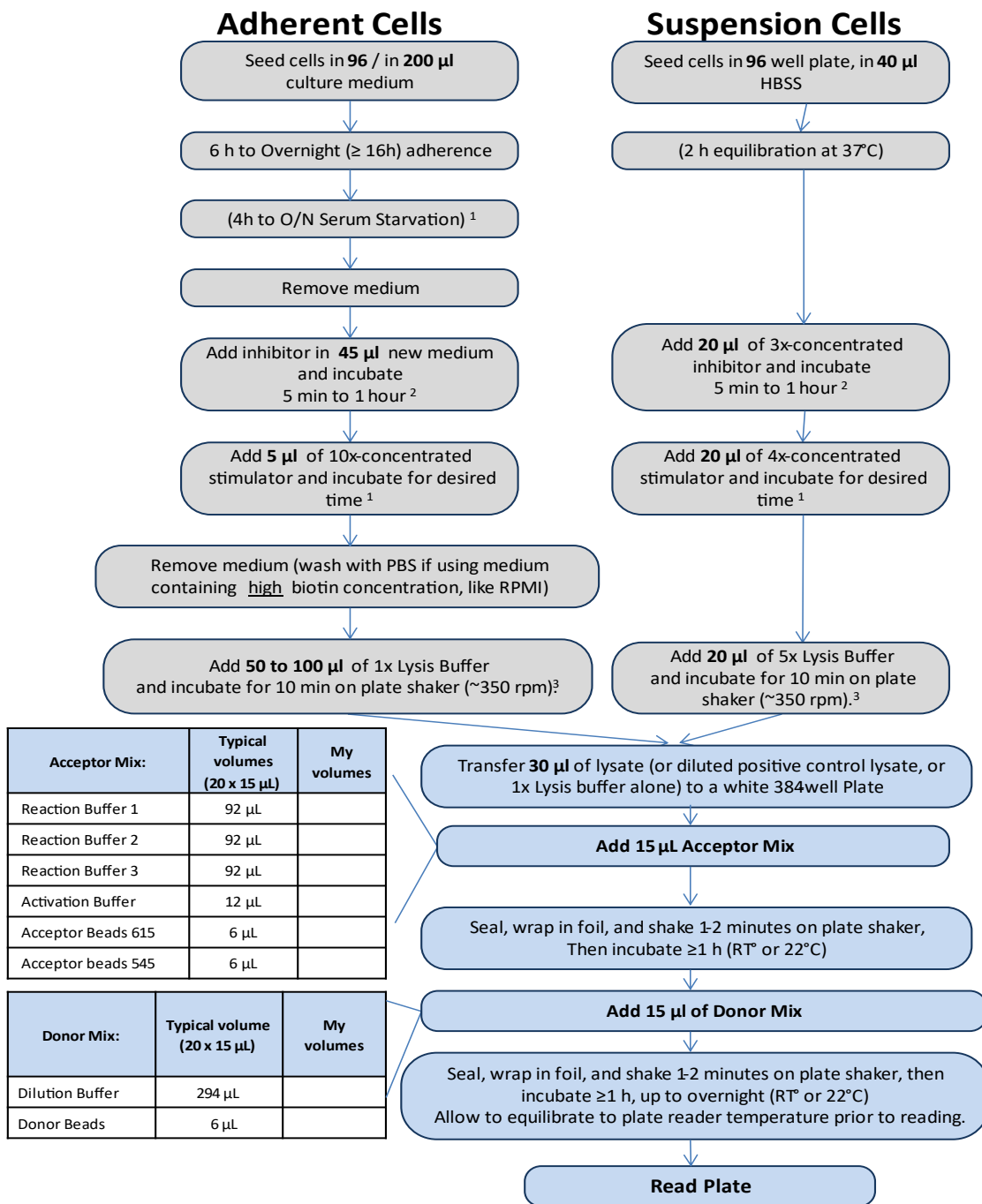
### Alpha SureFire Ultra Multiplex HV Assay

5. Add 15 µL of Acceptor Mix to wells. Seal plate with Topseal-A adhesive film. Incubate for 1 hour at room temperature.
6. Add 15 µL 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.

7. Read plate on an AlphaPlex Technology-compatible plate reader, using standard AlphaPlex settings.

**Alpha SureFire® Ultra™ HV Multiplex: 2-plates / 2-incubation assay flowchart**  
**Phospho / Total**



<sup>1</sup> Depending on cell type and pathway analyzed.

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

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

# Alpha SureFire® Ultra™ HV Multiplex PHOSPHO + TOTAL ASSAY PROTOCOL

## **B. 1 Plate Assay** – 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  $\mu$ L of cells/well into a 384-well white opaque culture plate.
2. If using test agents/inhibitors, add 2  $\mu$ L/well of 4X 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% BSA).

3. Return cells to incubator at 37°C for 1-2 hours.

### Cell Treatment

4. Treat cells with agonists/buffer by addition of 6  $\mu$ L/well of 4X agonist stock/buffer in HBSS containing 0.1% BSA. The final volume in the wells should be 24  $\mu$ L.

### Lysate Preparation

5. To lyse the cells, add 6  $\mu$ L/well of 5X Lysis Buffer. Add 30  $\mu$ L of Control lysates 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 15  $\mu$ 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 15  $\mu$ 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.

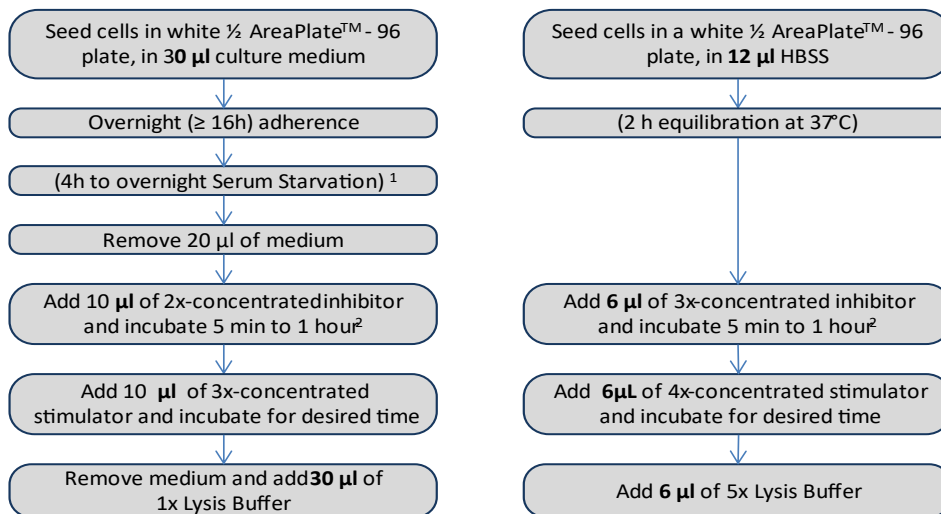
**Note:** Longer incubation may give greater sensitivity. Plates can be incubated overnight if required.

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

**Alpha SureFire® Ultra™ HV Multiplex: 1-plate / 2-incubation assay flowchart**  
**Phospho / Total**

**Adherent Cells**

**Suspension Cells**



Seal and incubate for 10 min on plate shaker (~350 rpm)<sup>3</sup>

In control wells, add 30 µL Control Lysate dilution or Lysis Buffer alone.

Acceptor Mix:	Typical volumes (20 x 15 µL)	My volumes
Reaction Buffer 1	92 µl	
Reaction Buffer 2	92 µl	
Reaction Buffer 3	92 µl	
Activation Buffer	12 µl	
Acceptor Beads 615	6 µl	
Acceptor beads 545	6 µl	

Add 15 µL Acceptor Mix

Seal, wrap in foil, and shake 1-2 minutes on plate shaker, Then incubate ≥1 h (RT<sup>o</sup> or 22°C)

Add 15 µl of Donor Mix

Seal, wrap in foil, and shake 1-2 minutes on plate shaker, then incubate ≥1 h, up to overnight (RT<sup>o</sup> or 22°C) Allow to equilibrate to plate reader temperature prior to reading.

Read Plate

Donor Mix:	Typical volume (20 x 15 µL)	My volumes
Dilution Buffer	294 µl	
Donor Beads	6 µL	

<sup>1</sup> Depending on cell type and pathway analyzed.

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

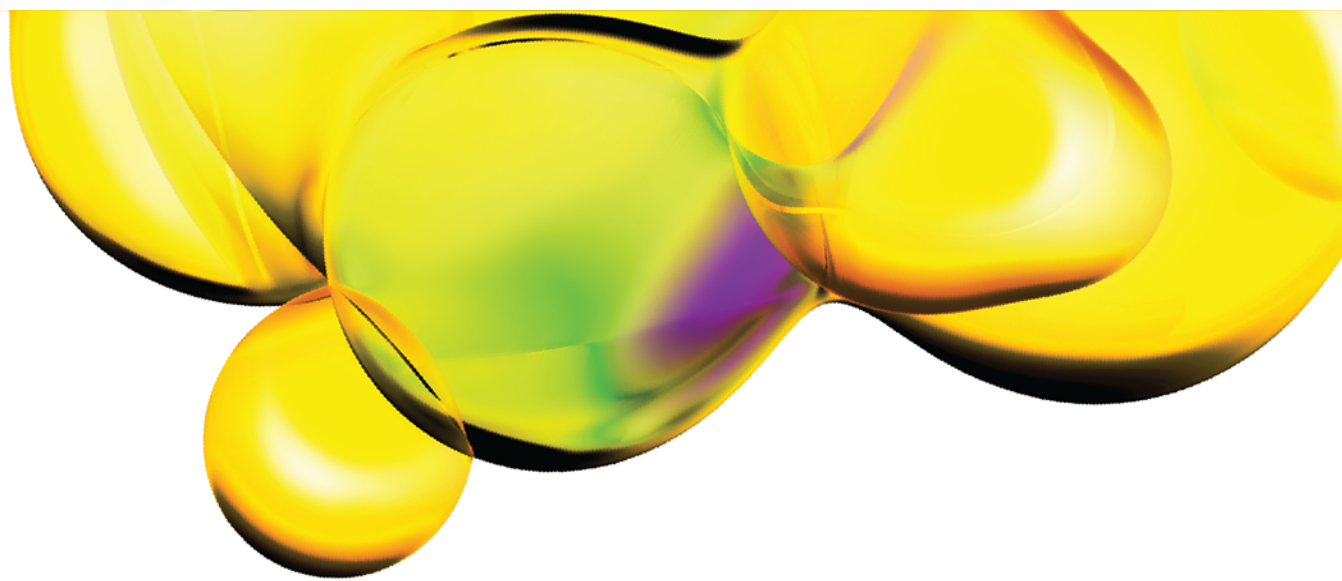
<sup>3</sup> May stop and freeze lysates at -20°C if desired. If doing this, re-shake after thawing to ensure homogeneity of lysate before pipetting.

## SUPPLEMENTARY BUFFERS AND BEADS

If using the standard protocol, sufficient amounts 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 catalog numbers:

Item	Suggested source	Catalog #	Size
Lysis Buffer (5X)	Revvity Inc.	ALSU-LB-10mL	10mL
	Revvity Inc.	ALSU-LB-100mL	100mL
Lysis Buffer B (5X)*	Revvity Inc.	ALSU-LBB-10mL	10 mL
	Revvity Inc.	ALSU-LBB-100mL	100 mL
Lysis Buffer C (5X)*	Revvity Inc.	ALSU-LBC-10mL	10 mL
	Revvity Inc.	ALSU-LBC-100mL	100 mL
Activation Buffer	Revvity Inc.	ALSU-AB-10mL	10mL
	Revvity Inc.	ALSU-AB-100mL	100mL
Activation Buffer B	Revvity Inc.	ALSU-ABB-10mL	10 mL
	Revvity Inc.	ALSU-ABB-100mL	100 mL
Activation Buffer C	Revvity Inc.	ALSU-ABC-10mL	10 mL
	Revvity Inc.	ALSU-ABC-100mL	100 mL
Dilution Buffer	Revvity Inc.	ALSU-DB-10mL	10mL
	Revvity Inc.	ALSU-DB-100mL	100mL
AlphaLISA™ CaptSure™ Acceptor Beads -2mg/ml	Revvity Inc.	ALSU-ACAB-0.06mL	60µL
	Revvity Inc.	ALSU-ACAB-1.2mL	1.2mL
	Revvity Inc.	ALSU-ACAB-6mL	6mL
Alpha Streptavidin Donor Beads -2mg/mL	Revvity Inc.	ALSU-ASDB-0.06mL	60µL
	Revvity Inc.	ALSU-ASDB-1.2mL	1.2mL
	Revvity Inc.	ALSU-ASDB-6mL	6mL
Alpha 545 (Tb) CaptSure2 Acceptor Beads -2mg/mL	Revvity Inc.	MPSU-CS2B-0.06mL	60µL
	Revvity Inc.	MPSU-CS2B -1.2mL	1.2mL
	Revvity Inc.	MPSU-CS2B -6mL	6mL

**\*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](http://www.revvity.com)

revvity



Revvity, Inc.  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.