

Alpha SureFire® Ultra™ Multiplex

Terbium SureFire® Ultra™ Assay Kit

For multiplexing with AlphaLISA® SureFire® Ultra™ (ALSU) assay kits

Assay Kits Manual

| Assay Points | Catalog # |
|--------------|----------------|
| 500 | TBSU-XXXX-X500 |
| 10 000 | TBSU-XXXX-X10K |
| 50 000 | TBSU-XXXX-X50K |

This Manual is a generic manual for all the 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.perkinelmer.com

For Research Use Only. Not for use in Diagnostic Procedures.

Note: For kit handling and disposal information see page 5 of this manual

Alpha SureFire® Ultra™ Multiplex

Assay Principle

The Alpha SureFire Ultra Multiplex kits allow the rapid, sensitive, and quantitative detection of two phosphoprotein targets in each well of an assay plate. This Multiplex measurement is achieved by the use of two types of Alpha Acceptor beads that emit at distinct wavelengths (Terbium, 545nm and Europium, 615nm).

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

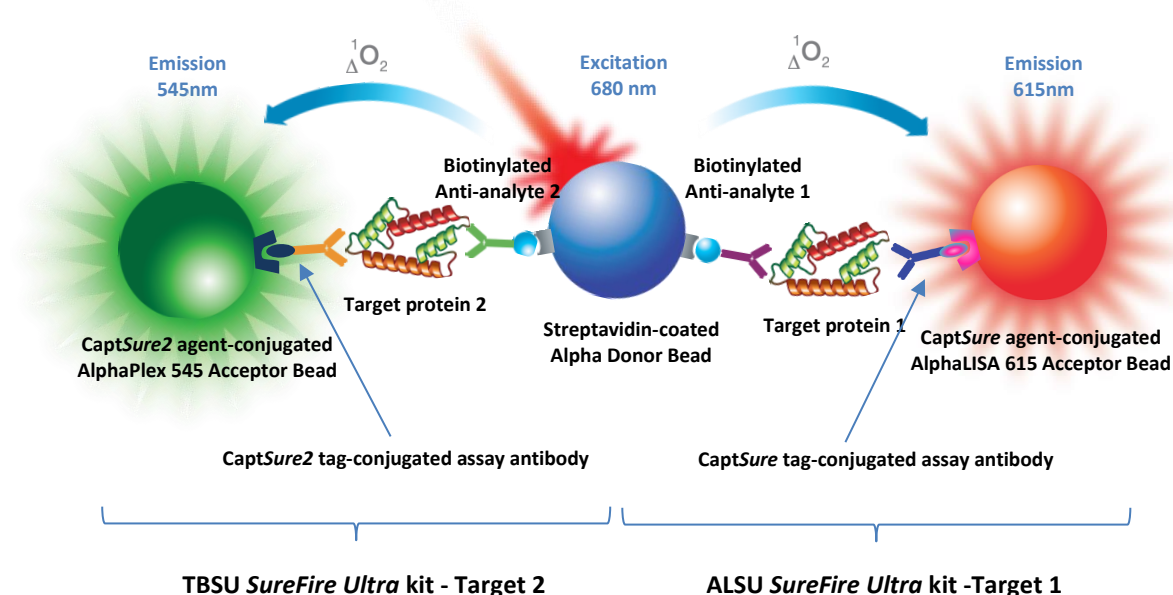
Terbium SureFire Ultra Assay kit for multiplexing with an AlphaLISA SureFire Ultra assay kit:

The Terbium SureFire Ultra (TBSU) assay kits contain reagents to be combined with those of an AlphaLISA SureFire Ultra (ALSU) assay kit, to allow multiplexed measurement of two different phosphoprotein targets. Antibodies in each TBSU or ALSU kit measure a single target. Combining two such kits allows the highly flexible selection of two phosphoprotein targets to be measured.

The terbium “Alpha 545 Acceptor Beads” in the TBSU kit are conjugated with a “CaptSure2” agent, which binds a specific CaptSure2 tag on one of the assay antibodies. In combination with the biotinylated antibody provided, the 545 nm signal generated will report levels of the phosphoprotein of interest (Target 2). The TBSU kit is to be used in conjunction with the reagents in an ALSU kit, such that the ALSU kit will report that target of interest by Eu emission at 615nm (Target 1). This is shown diagrammatically below.

The protocol details how to combine these kits to allow multiplexed measurement of the desired targets.

Alpha SureFire® Ultra™ Multiplex Technology



General Information on the Terbium *SureFire*® *Ultra*™ Assay kit

The Terbium *SureFire Ultra* Assay (TBSU) kit is used to measure the phosphorylation of a target of interest plus another target of choice from the AlphaLISA *SureFire Ultra* (ALSU) list of kits. The TBSU kit must be used in conjunction with an ALSU kit, and **can not** be used alone to measure a single target.

Multiplexing is achieved through the use of two Alpha Acceptor beads. The first (545nm Tb) in this TBSU kit will report on the levels of the designated target of interest. The second (615nm Eu) Acceptor bead from another AlphaLISA *SureFire Ultra* kit reports on the protein designed to be measured using that kit.

The Donor beads in this kit are identical to those in a standard AlphaLISA *SureFire Ultra* kit.

Control lysates are provided in each kit to allow testing of the signal generated in the two channels (545 and 615 nm).

This kit has been formulated to provide optimal signal:background (i.e. S:B) assay windows, and to perform without interference in the presence of extraneous antibodies.

Unless otherwise indicated, the antibodies in each TBSU kit are identical to the antibodies to the same target in the corresponding ALSU kit, except for the CaptSure or CaptSure2 tag.

The assay utilizes the bead-based Alpha Technology, and requires an Alpha Technology-compatible plate reader capable of reading dual emission wavelengths. See www.perkinelmer.com/AlphaPlex 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.

Kit-Specificity Information / Control Lysate Information / Representative Data

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

Kit Contents

| | Kit Size | | |
|--|--|---------------|---------------|
| | 500 points | 10,000 points | 50,000 points |
| Reaction Buffer 3 – TBSU <i>(Biotinylated anti-target antibody)</i> | 1 x 0.15 mL | 1 x 2.8 mL | 1 x 14 mL |
| Reaction Buffer 4 – TBSU <i>(CaptSure2-tagged anti-target antibody)</i> | 1 x 0.15 mL | 1 x 2.8 mL | 1 x 14 mL |
| Alpha 545 CaptSure2 Acceptor Beads (2mg/mL in PBS plus 0.05% Proclin-300) | 1 x 0.06 mL | 1 x 1.1 mL | 1 x 5.5 mL |
| Alpha Streptavidin Donor Beads (2mg/mL in PBS plus 0.05% Proclin-300) | 1 x 0.06 mL | 1 x 1.1 mL | 1 x 5.5 mL |
| Positive Control Lysate | 1 lyophilized tube to be re-dissolved with 250 µL H ₂ O | | |

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

IMPORTANT: The components of this kit will be used in conjunction with the components of another standard AlphaLISA SureFire Ultra (ALSU) kit to provide multiplexing. On page 8 we detail how to set up the assay solutions from both kits.

Kit Content Handling Conditions

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 <http://www.perkinelmer.com>

Lysis Buffer (5X) - *Ultra* (from AlphaLISA *SureFire Ultra* kit)

Lysis Buffer (5X) - *Ultra* 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.

Lysis Buffer B (5X) and Lysis Buffer C (5X) are assay specific and should not be inter-changed.

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.

Activation Buffer (from AlphaLISA *SureFire Ultra* kit)

Activation Buffer precipitates at 4°C. To re-dissolve, warm to 37°C and mix before each use.

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 should 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 Certificate of Analysis for the recommended dilution in the linear range of the assay.

Storage Conditions Upon Receipt

Expiry date indicated on kit box.

| | | |
|---------------------|----------------------------------|--|
| Unopened kit | | 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. |
| Opened kit | Reaction Buffer 3 - <i>Ultra</i> | Store at 4°C. |
| | Reaction Buffer 4 - <i>Ultra</i> | |
| | Acceptor/Donor Beads | Store at 4°C, in the dark zip lock bag or box provided. |
| | Positive Control Lysate | Store at 4°C or for long term storage at -20°C. |

Materials Required But Not Provided

| Item | Suggested source | Catalog # | Size |
|--|------------------|-----------|---------|
| Optiplate™-384, White Opaque assay plate ⁽¹⁾ | PerkinElmer Inc. | 6007290 | 50/box |
| AlphaPlate-384, Light gray assay plate ⁽²⁾ | PerkinElmer Inc. | 6005350 | 50/box |
| CulturPlate-384, white, sterile, with lid, 384-well ⁽³⁾ | PerkinElmer Inc. | 6007680 | 50/box |
| SpectraPlate-96, TC, clear, 96-well, tissue culture treated, sterile, with lids ⁽⁴⁾ | PerkinElmer Inc. | 6005650 | 50/box |
| TopSeal-A 384, clear adhesive sealing film | PerkinElmer Inc. | 6050185 | 100/box |
| Envision® Alpha-reader with adequate AlphaPlex filters (see table below) | PerkinElmer Inc. | - | - |

(1) Plates used for the immunoassay in the 2-plates assay; (2) Same as (1) but optimal if cross-talk needs to be reduced; (3) plates for assays run in a 1-plate assay (from cell seeding to immunoassay); (4) plates used to seed and stimulate cells before Lysis and transfer of lysate in an immunoassay plate in the 2-plates assay.

For more assay plates options, please go to www.perkinelmer.com/microplates

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.perkinelmer.com/AlphaPlex

| | 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 simultaneous 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 |

Buffer Preparation and Subsequent Storage Conditions

1. Prepare sufficient Acceptor Mix from the ALSU AlphaLISA *SureFire Ultra* kit as you would for a standard single target assay, according to that kit manual.
2. The Terbium *SureFire Ultra* Assay kit reagents will be added to the pre-prepared ALSU Acceptor mix from step 1 to complete the Multiplex Acceptor Mix.
3. We recommend that you combine the Donor beads from both ALSU and TBSU kits into a single tube prior to use. Donor Mix will then be made up as indicated below.

| Components | Preparation |
|--|--|
| 1X Lysis Buffer Note: Lysis Buffer from ALSU kit | Dilute 5X Lysis buffer in deionised water to a final concentration of 1X For example: for 10 mL of 1X Lysis Buffer, add: 2 mL of 5X Lysis Buffer – <i>Ultra</i> to 8 mL deionised water. Discard unused 1X buffer. |
| <u>Multiplex Acceptor Mix</u> Prepare Acceptor Mix as normal from ALSU <i>SureFire Ultra</i> kit. <u>To this:</u> Add TBSU Reaction Buffer 3 (1:20) Add TBSU Reaction Buffer 4 (1:20) Vortex Add CaptSure2 Alpha 545 Acceptor Beads (1:50) Vortex See flowchart table | Important: Combine the reagents in the order indicated below <u>For example: for 60 samples (require minimum 300 µL of Duoplex Acceptor Mix):</u> Make up 300 µL of ALSU kit Acceptor Mix, according to that kit's protocol. To this ALSU Acceptor Mix add 15 µL of Reaction Buffer 3 and 15 µL of Reaction Buffer 4. Mix combined reagents by vortex. Then add 6µL CaptSure2 Alpha 545 Acceptor Beads. Mix again by vortex. The Multiplex Acceptor mix should be made up and used within 30min for best results. Excess mix should be discarded. |
| <u>Donor Mix*</u> Alpha Donor beads (1:25) (4 parts or 4%) Note: Dilution buffer from ALSU kit See flowchart for table * Prepare and use under low-light conditions. | Dilute Donor beads 25-fold in Dilution buffer <u>For example: for 60 samples (require minimum 300 µL of Donor Mix):</u> Add 12 µL Donor Beads to 288 µL of Dilution Buffer. Note: Dilution Buffer from ALSU kit. The Donor mix should be made up and used within 30min for best results. Excess mix should be discarded. |
| Positive control lysate | Reconstitute with deionised water. Store at -20°C in single use aliquots and use within 3 months. Dilute as required. |

Dual target measurement using combined Terbium *SureFire® Ultra™* Assay kit + AlphaLISA® *SureFire® Ultra™* target assay kit

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 - *Ultra* (50-100 µL for a 96 well plate, 25 µL for a 384 well plate). Agitate on a plate shaker (~350 rpm) for 10 minutes at room temperature.
4. Take 10 µL of the lysate and transfer to a 384-well Optiplate™ for assay. Add 10 µL 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.*

Alpha SureFire Ultra Multiplex Assay

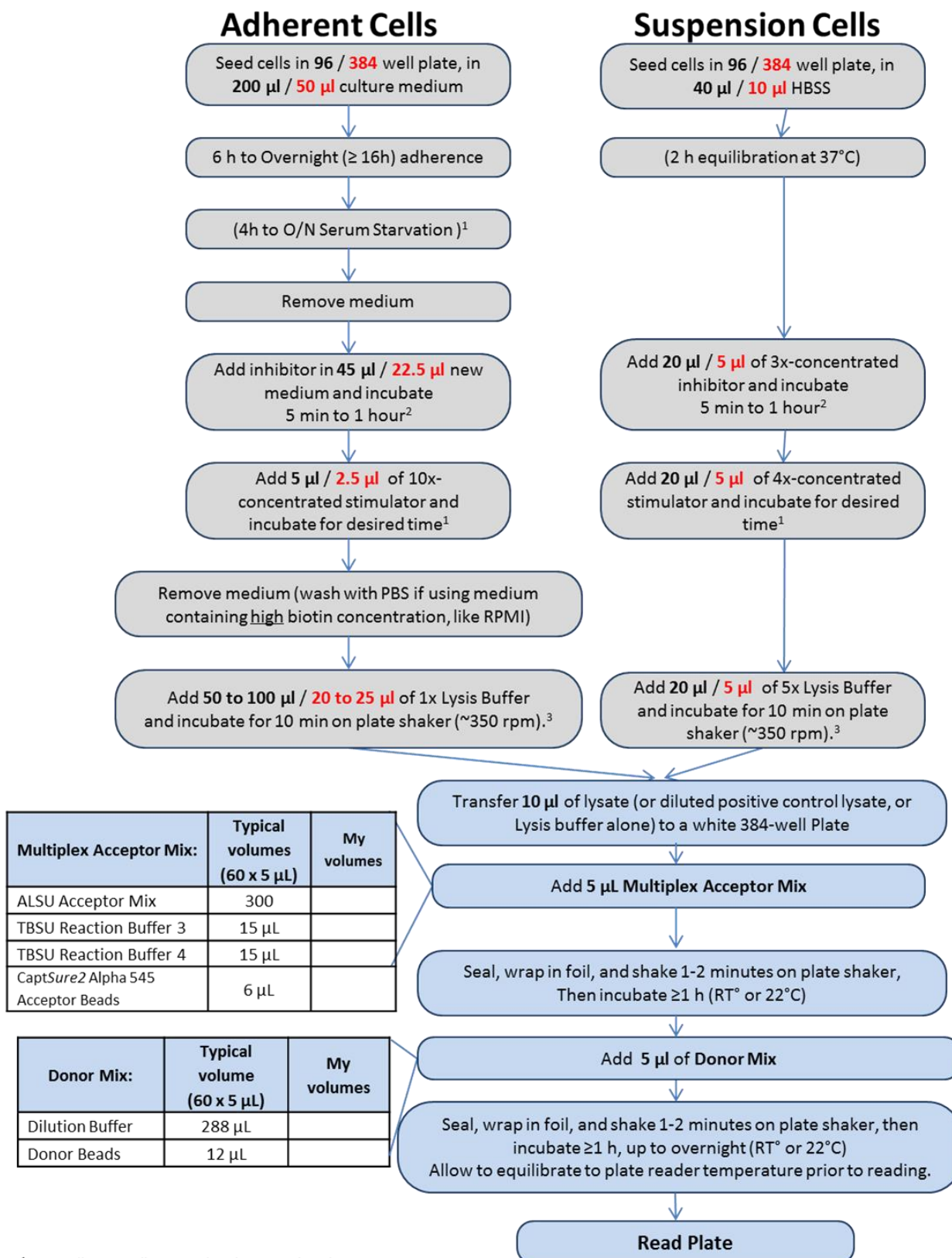
5. Add 5 µL of Multiplex Acceptor Mix to wells. Seal plate with Topseal-A adhesive film. Incubate for 1 hour at room temperature.
6. Add 5 µL of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature in the dark.

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 (see above).

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

Dual Targets



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.

Dual target measurement using combined Terbium *SureFire® Ultra™* Assay kit + AlphaLISA® *SureFire® Ultra™* target assay kit

B. 1 Plate Assay - assay protocol for non-adherent 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 (eg PerkinElmer Cat # 6007680).
2. If using test agents/inhibitors, add 2 μ 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. 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.)

Lysate Preparation

5. To lyse the cells, add 2 μ L/well of 5X Lysis Buffer - *Ultra*. Add 10 μ 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.

Alpha SureFire Ultra Multiplex Assay

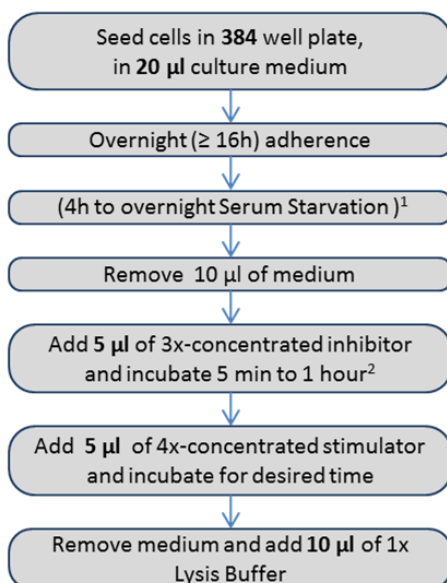
6. Add 5 μ L of Multiplex Acceptor Mix to wells. Seal plate with Topseal-A adhesive film. Incubate for 1 hour at room temperature.
7. Add 5 μ L of Donor Mix to wells under subdued light. Seal plate with Topseal-A adhesive film, and cover plate with foil. Incubate for 1 hour at room temperature in the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

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

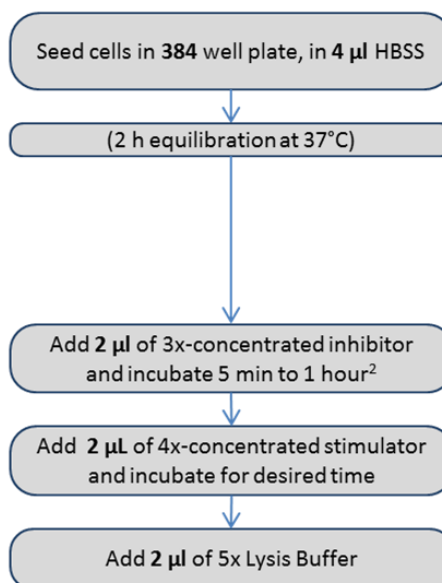
8. Read plate on an AlphaPlex Technology-compatible plate reader, using standard AlphaPlex settings (see above).

Alpha SureFire® Ultra™ Multiplex: 1-plate / 2-incubation assay flowchart Dual Targets

Adherent Cells



Suspension Cells



Seal and incubate for 10 min on plate shaker (~350 rpm).³

In control wells, add 10 µL positive control lysate dilution or lysis buffer alone.

| Multiplex Acceptor Mix: | Typical volumes (60 x 5 µL) | My volumes |
|------------------------------------|-----------------------------|------------|
| ALSU Acceptor Mix | 300 | |
| TBSU Reaction Buffer 3 | 15 µL | |
| TBSU Reaction Buffer 4 | 15 µL | |
| CaptSure2 Alpha 545 Acceptor Beads | 6 µL | |

Add 5 µL Multiplex Acceptor Mix

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

Add 5 µL of Donor Mix

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

Read Plate

| Donor Mix: | Typical volume (60 x 5 µL) | My volumes |
|-----------------|----------------------------|------------|
| Dilution Buffer | 288 µL | |
| Donor Beads | 12 µL | |

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 amounts of buffers and beads are provided in the kits. 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) - <i>Ultra</i> | PerkinElmer Inc. | ALSU-LB-10mL | 10mL |
| | PerkinElmer Inc. | ALSU-LB-100mL | 100mL |
| Activation Buffer - <i>Ultra</i> | PerkinElmer Inc. | ALSU-AB-10mL | 10mL |
| | PerkinElmer Inc. | ALSU-AB-100mL | 100mL |
| Dilution Buffer - <i>Ultra</i> | PerkinElmer Inc. | ALSU-DB-10mL | 10mL |
| | PerkinElmer Inc. | ALSU-DB-100mL | 100mL |
| Alpha Streptavidin Donor Beads -2mg/mL | PerkinElmer Inc. | ALSU-ASDB-0.06mL | 60µL |
| | PerkinElmer Inc. | ALSU-ASDB-1.2mL | 1.2mL |
| | PerkinElmer Inc. | ALSU-ASDB-6mL | 6mL |

Useful Links

For FAQ and troubleshooting, please go to:

www.perkinelmer.com/AlphaPlex

For a complete list of AlphaLISA *SureFire Ultra* and Alpha *SureFire Ultra* Multiplex kits, please go to:

www.perkinelmer.com/category/alpha-surefire-kits

For technical support please go to: www.perkinelmer.com/ASK

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer, Inc. warrants that, at the time of shipment, the products sold by it are free from defects in material and workmanship and conform to specifications which accompany the product. PerkinElmer Inc. makes no other warranty, express or implied with respect to the products, including any warranty of merchantability or fitness for any particular purpose. Notification of any breach of warranty must be made within 60 days of receipt unless provided in writing by PerkinElmer Inc. No claim shall be honored if the customer fails to notify PerkinElmer Inc. within the period specified. The sole and exclusive remedy of the customer for any liability of PerkinElmer Inc. of any kind including liability based upon warranty (express or implied whether contained herein or elsewhere), strict liability contract or otherwise is limited to the replacement of the goods or the refunds of the invoice price of goods. PerkinElmer Inc. shall not in any case be liable for special, incidental or consequential damages of any kind.

List of Available TBSU kits

| Assay | Catalogue Numbers | | |
|---|-------------------|-----------------|-----------------|
| | 500 point | 10,000 point | 50,000 point |
| p-4E-BP1 (T37/46) | TBSU-P4EBP-A500 | TBSU-P4EBP-A10K | TBSU-P4EBP-A50K |
| p-Akt1/2/3 (T308) | TBSU-PAKT-A500 | TBSU-PAKT-A10K | TBSU-PAKT-A50K |
| p-Akt1/2/3 (S473) | TBSU-PAKT-B500 | TBSU-PAKT-B10K | TBSU-PAKT-B50K |
| Total Cofilin | TBSU-TCOF-A500 | TBSU-TCOF-A10K | TBSU-TCOF-A50K |
| p-CREB (S133) | TBSU-PCREB-A500 | TBSU-PCREB-A10K | TBSU-PCREB-A50K |
| p-EGF Receptor (Y1068) | TBSU-PEGFR-A500 | TBSU-PEGFR-A10K | TBSU-PEGFR-A50K |
| p-eIF2 α (S51) | TBSU-PEIF2-B500 | TBSU-PEIF2-B10K | TBSU-PEIF2-B50K |
| p-eIF4E (S209) | TBSU-PEIF4-A500 | TBSU-PEIF4-A10K | TBSU-PEIF4-A50K |
| p-ERK1/2 (T202/Y204) | TBSU-PERK-A500 | TBSU-PERK-A10K | TBSU-PERK-A50K |
| p-GSK3 β (S9) | TBSU-PGS3B-A500 | TBSU-PGS3B-A10K | TBSU-PGS3B-A50K |
| p-IGF-1 Receptor β (Y1135/1136) | TBSU-PIGFR-B500 | TBSU-PIGFR-B10K | TBSU-PIGFR-B50K |
| p-IKK α (S176/180) | TBSU-PIKKA-A500 | TBSU-PIKKA-A10K | TBSU-PIKKA-A50K |
| p-Insulin Receptor β (Y1150/1151) | TBSU-PINR-A500 | TBSU-PINR-A10K | TBSU-PINR-A50K |
| p-JNK1/2/3 (T183/Y185) | TBSU-PJNK-A500 | TBSU-PJNK-A10K | TBSU-PJNK-A50K |
| p-MEK1 (S218/222) | TBSU-PMEK1-A500 | TBSU-PMEK1-A10K | TBSU-PMEK1-A50K |
| p-mTOR (S2448) | TBSU-PMTOR-A500 | TBSU-PMTOR-A10K | TBSU-PMTOR-A50K |
| p-NF- κ B p65 (S536) | TBSU-PNFKB-A500 | TBSU-PNFKB-A10K | TBSU-PNFKB-A50K |
| p-p38 MAPK (T180/Y182) | TBSU-PP38-B500 | TBSU-PP38-B10K | TBSU-PP38-B50K |
| p-p70 S6K (T389) | TBSU-PP70-A500 | TBSU-PP70-A10K | TBSU-PP70-A50K |
| p-Ribosomal Protein S6 (S240/244) | TBSU-PS6R-A500 | TBSU-PS6R-A10K | TBSU-PS6R-A50K |
| p-SLP-76 (S376) | TBSU-PSLP-A500 | TBSU-PSLP-A10K | TBSU-PSLP-A50K |
| p-SMAD1 (S463/465) | TBSU-PSM1-A500 | TBSU-PSM1-A10K | TBSU-PSM1-A50K |
| p-SMAD3 (S423/425) | TBSU-PSM3-A500 | TBSU-PSM3-A10K | TBSU-PSM3-A50K |
| p-STAT1 (S727) | TBSU-PST1-B500 | TBSU-PST1-B10K | TBSU-PST1-B50K |
| p-STAT1 (Y701) | TBSU-PST1-A500 | TBSU-PST1-A10K | TBSU-PST1-A50K |
| p-STAT3 (Y705) | TBSU-PST3-A500 | TBSU-PST3-A10K | TBSU-PST3-A50K |
| p-STAT4 (Y693) | TBSU-PST4-A500 | TBSU-PST4-A10K | TBSU-PST4-A50K |
| p-STAT5 (Y694/699) | TBSU-PST5-B500 | TBSU-PST5-B10K | TBSU-PST5-B50K |
| p-STAT6 (Y641) | TBSU-PST6-A500 | TBSU-PST6-A10K | TBSU-PST6-A50K |
| p-SYK (Y525/526) | TBSU-PSYK-A500 | TBSU-PSYK-A10K | TBSU-PSYK-A50K |
| p-VEGF Receptor 2 (Y1175) | TBSU-PVGFR-A500 | TBSU-PVGFR-A10K | TBSU-PVGFR-A50K |