

G-LISA[®] Activation Assays

Technical Guide



Important Notice

This technical guide is an accompaniment to the G-LISA kit Protocol and is designed to provide additional information that may be of use when designing or troubleshooting an activation assay.

Content

Section I:	Introduction	4
Section II:	Preparation of Lysates for G-LISA Assays	
	Recommended lysate amount per assay	5
	Lysis buffer compatibility between assays	6
	Growth and treatment of tissue culture cell lines	7-13
	Preparation of lysates from 2D tissue culture cell lines	14
	Preparation of lysates from 3D tissue culture cell lines	15
	Preparation of lysates from tissue samples	15
	Equalization of cell lysate concentrations	16-17
Section III:	Assay technical tips	
	Use of a multichannel pipettor	18
	Plate shaker recommendations	18
	Vortex of samples after binding buffer addition	18
	Spectrophotometer / Luminometer settings	18-19
Section IV:	Data Analysis	
	Excel format	20-21
	Experimental record Sheet	22
	Plate Record Template	23
Section V:	References	24-26

Section I: Introduction

Technical Guides are an accompaniment to the G-LISA kit Protocols and are designed to provide additional information to the end user that may be of use when designing or troubleshooting an activation assay.

Go online to www.cytoskeleton.com for the latest Technical Guide updates.

For assistance, contact Technical Support at 303-322-2254 or e-mail tservice@cytoskeleton.com.

Section II: Preparation of Lysates for G-LISA

Recommended lysate amounts per assay

The G-LISA range of activation assays generally requires 6.25-50 µg of lysate per assay. Table 1 summarizes the recommended concentrations of lysate for each small G-protein (GTPase) target. The values in Table 1 are recommended starting points for optimizing a given G-LISA assay. Lysates that are too far above the recommended range may contain a high level of GAPs (GTPase activation proteins) that can inactivate the target GTPase, even in lysis buffer. Final lysate concentrations will vary depending upon experimental factors such as;

- A) Total amount of a given GTPase in your cell line or tissue: In general cell lines or tissue that have a high endogenous level of the target GTPase will give a more robust response to an activator.
- B) Degree of activation achieved under your experimental conditions: As a general guideline roughly 2-10% of total cellular small G-protein is activated in response to a given stimuli.

Table 1: Recommended lysate amounts for G-LISA assays

Small G-protein target	Recommended lysate amount per G-LISA assay (µg)	Recommended lysate concentration range (mg/ml)	Volume of lysate per assay (µl)
Arf1	12.50 - 25.00	0.50 - 1.00	25
Arf6	12.50 - 25.00	0.50 - 1.00	25
Cdc42	7.50 - 50.00	0.15 - 1.00	50
RalA	6.25 - 25.00	0.25 - 1.00	25
Rac1	12.50 - 50.00	0.25 - 1.00	50
Rac1,2,3	6.25 - 25.00	0.25 - 1.00	25
Ras (pan)	6.25 - 25.00	0.25 - 1.00	25
RhoA	10.00 - 50.00	0.40 - 2.00	25

Section II: Preparation of Lysates for G-LISA

Lysis buffer compatibility between assays

Lysis buffer composition is critical to optimal performance of a G-LISA assay. In most assays a standard lysis buffer composition (GL36) has been used (see Table 2). This allows the end user to test various GTPase activities in the same lysate.

For Arf6 analysis one can dilute a GL36 lysate 1:1 with water to achieve the correct assay lysis buffer.

It should be noted that the Cdc42 lysis buffer (GL35) is not compatible with GL36 lysates.

Table 2: List of G-LISA Lysis buffers

Small G-protein target	Lysis Buffer in Kit	Notes
Arf1	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.
Arf6	GL70	Buffer composition is 0.5X GL36
Cdc42	GL35	Identical components to GL36, composition varies as follows; Tris (1X GL36), MgCl ₂ (8X GL36), NaCl (2X GL36), IGEPAL (1X GL36), SDS (5X GL36)
RalA	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.
Rac1	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.
Rac1,2,3	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.
Ras	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.
RhoA	GL36	Standard G-LISA buffer, compatible with most G-LISA assays. Composed of a proprietary formulation of Tris pH 7.5, MgCl ₂ , NaCl, IGEPAL and SDS.

Section II: Preparation of Lysates for G-LISA

Growth and treatment of tissue culture cells

The health and responsiveness of your cell line is the single most important parameter for the success and reproducibility of GTPase activation assays. Growth conditions vary widely between cell lines, therefore the reader is referred to Tables 3A-H for helpful references regarding GTPase activation in a variety of cell lines. Optimal GTPase activation/inhibition conditions in a given cell line will need to be determined empirically.

Some general guidelines are given below;

- 1) When possible, untreated samples should have low basal cellular levels of GTPase activity (controlled state). For example, serum starvation often results in low basal levels of GTPase activity and leads to a much greater response to a given activator (see Tables 3A-H).
- 2) GTPase activation is often very transient, lasting only seconds to minutes. A time courses of activation/inhibition and titration of activating or inhibiting factors are highly recommended when establishing assay conditions.
- 3) When possible cells should be checked for their responsiveness (responsive state) to a known stimulus (see Tables 3A-H).
- 4) Poor culturing technique such as sub-culturing of cells that have previously been allowed to become overgrown can result in essentially non-responsive cells. For example, Swiss 3T3 cells grown to >70% confluency should not be used for Rho GTPase activation studies.
- 5) Cell confluency can affect the cellular response of GTPases to activators/inhibitors. For example RhoA activation generally requires sub confluent cells for a robust response to activators. Confluency restrictions vary widely between different cell lines and different GTPases (see Tables 3A-H).
- 6) Once an inducible GTPase response has been achieved, it is particularly important to maintain consistent experimental conditions. An Experimental Record Sheet is included in this Guide (Data Analysis section) and is useful for tracking key experimental parameters.

Section II: Preparation of Lysates for G-LISA

Table 3A: Transfection-based GTP-Arf1 modulation

Gene transfected	Treatment	Cell type used	Response	Assay Type	Ref.
Arf1-HA + GBF1	5-25 μ g/ml BFA	Cos7 cells	GBF1 increased Arf1-GTP levels 2-3 fold and this activation was not inhibited by BFA.	GGA3 pull-down	1
HCV NS5A	N/A	Huh7 cells	Hepatitis C virus NS5A protein expression reduced Arf1-GTP levels	GGA3 pull-down	2
Ephrin-B1	2 μ g/ml EphB2-Fc	Capan-1 cells Panc-1 cells	Treatment for 1.5 hr with EphB2-Fc resulted in elevated Arf1-GTP levels	GGA3 pull-down	3
Enterovirus 3A protein	N/A	Cos7 cells	Expression of enterovirus 3A protein decreased total Arf1-GTP levels by 60%	GGA3 pull-down	4

Table 3B: Non-transfection-based GTP-Arf1 modulation

Arf1 Modulator	Treatment	Cell type used	Response	Assay Type	Ref.
Formyl methionyl-lysyl-phenylalanine (fMLF)	5 ng/ml	Polymorphonuclear neutrophils (PMNs), PLB-985 cells	Treatment for 2 minutes with fMLF resulted in ~5-fold activation of Arf1 in PMNs.	GGA3 pull-down	5
QS11 (ARFGAP1 inhibitor)	1-2.5 μ M	NIH 3T3 cells	Treatment for 36 hr resulting in dose-dependent activation of Arf1 and Arf6	GGA3 pull-down	6
Poliovirus	10 PFU/cell	HeLa cells	Time-dependent increase in Arf1-GTP with a maximal activation of ~3.75-fold	GGA3 pull-down	7
Ephrin-B1	2 μ g/ml EphB2-Fc	SUIT-4 cells	Treatment for 1.5 hr with EphB2-Fc resulted in elevated Arf1-GTP levels	GGA3 pull-down	3

Section II: Preparation of Lysates for G-LISA

Table 3C: Modulators of GTP-Arf6 levels in cells

Arf6 Modulator	Treatment	Cell type used	Response	Assay Type	Ref.
Fibronectin (Fn)	5 ng/ml Fn coated plates	Mouse embryonic fibroblasts (MEFs)	Control cells harvested and kept in suspension for 90 min. followed by replating to Fn-coated plates for 15-30 min. resulted in a 4-fold activation of Arf6	GGA3 pull-down	8
Fibronectin (Fn)	5 ng/ml Fn coated plates	NIH 3T3 cells	Control cells harvested and kept in suspension for 60 min. followed by replating to Fn-coated plates for 5 min. resulted in a modest activation of Arf6	GGA3 pull-down	9
Formyl methionyl-lysyl-phenylalanine (fMLF)	5 ng/ml	Polymorphonuclear neutrophils (PMNs), PLB-985 cells	Treatment for 2 min. with fMLF resulted in ~10-fold activation of Arf6 in PMNs, and a ~ 3.5-fold activation in PLB-985 cells	GGA3 pull-down	5
QS11 (ARFGAP1 inhibitor)	1-2.5 μ M	NIH 3T3 cells	Treatment for 36 hr resulting in dose-dependent activation of Arf6	GGA3 pull-down	6
Collagen	10 μ g/ml	Platelets	Time-dependent decrease in Arf6-GTP levels with collagen treatment	GGA3 pull-down	10
Insulin	100 nM	HepG2 cells	Treatment of cells for 20 min. with insulin resulted in ~3-fold increase in Arf6-GTP	GGA3 pull-down	11
HGF	10 ng/ml	MDCK cells	Time-dependent increase in Arf6-GTP up to 16 hr. with ~30-fold maximal increase in Arf6-GTP	GGA3 pull-down	12
Semaphorin 3E	100 ng/ml	SVEC cells	Time-dependent ~2-fold increase in Arf6-GTP	GGA3 pull-down	13

Section II: Preparation of Lysates for G-LISA

Table 3D: Modulators of GTP-Cdc42 levels in cells

Cdc42 Modulator	Treatment	Cell type used	Response	Assay Type	Ref.
ML-7	20 μ M	Primary rat microglia	Inhibition of Pyk2 mediated Cdc42 activation	Cdc42 G-LISA	14
Epidermal Growth Factor	50 ng/ml	Mouse mesenchymal stem cells	1.5 fold activation of Cdc42	Cdc42 G-LISA	15
Ethanol	30 mM	Rat hippocampal neurons	35% reduction in Cdc42 activation	Cdc42 G-LISA	16

Table 3E: Modulators of GTP-Rac levels in cells

Rac Activator	Treatment	Cell Line Used	Response	Assay Type	Ref.
Epidermal Growth Factor	50 ng/ml	US7MG human glioblastoma	1.5 fold activation after 5 minutes with 2D cultures. 1.3 fold activation in 3D cultures	Rac G-LISA [®]	17
MCP-1	10 ng/ml	Murine alveolar macrophages	Maximal activation at 4h	Rac G-LISA [®]	18
Heregulin beta1	0-30 ng/ml	Breast cancer cell lines	Dose-dependent activation	PAK-PBD pulldown assay	19
Interleukin-3	5 μ g/ml for 5 min	Primary bone marrow derived mast cells	2 fold increase over control	PAK-PBD pulldown assay	20

Section II: Preparation of Lysates for G-LISA

Table 3F: Modulators of GTP-RalA levels in cells

RalA Modulator	Treatment	Cell type used	Response	Assay Type	Ref.
Epidermal growth factor	50 ng/ml	Cos7	FRET activity and lamellipodia induction	FRET	21

Table 3G: Modulators of GTP-Ras levels in cell

Ras Activator	Treatment	Cell line used	Response	Assay Type	Ref.
Epidermal Growth Factor	100 ng/ml 5 minutes	HeLa	Dose-dependent activation	Raf1-RBD pulldown assay	22
Hepatocyte Growth Factor	100 ng/ml 5 minutes	HeLa	Dose-dependent activation	Raf1-RBD pulldown assay	22
IL-3	50 ng/ml 5 minutes	BaF3	14 fold activation after 5 minutes	Raf1-RBD pulldown assay	24
aCD3 + aCD28	5 mg/ml each 10 minutes	Jurkat	Dose-dependent activation	Raf1-RBD pulldown assay	24
PMA ionomycin	100 ng/ml 500 ng/ml 10 minutes	Jurkat	Dose-dependent activation	Raf1-RBD pulldown assay	24

Section II: Preparation of Lysates for G-LISA

Table 3H: Modulators of GTP-Rho levels in cell

Rho Activator	Treatment	Cell line used	Response	Assay Type	Ref.
Lysophosphatidylcholine (<i>lysophospholipid</i>)	100 nM	Human melanocyte	After 1 min, 1.5 fold increase over control. Activation peaked at 5 min (5.3 fold increase) and declined to baseline by 60 min.	RhoA G-LISA (Cat.# BK124)	25
Calpeptin (Cat.# CN01)	0.5 µM/ml	Uterine myocyte	After 15 min treatment, 2.4 fold increase over vehicle control	RhoA G-LISA (Cat.# BK124)	26
Bombesin	10 nM	Swiss 3T3 cells	Maximal activation after 1 min which is sustained for at least 30 min	Actin morphology	27
Calpeptin	100 µg/ml	REF-52 fibroblast & Swiss 3T3 cells	Maximal activation after 5 to 10 min with extended activation time up to 30 min, decreasing thereafter to basal levels after 60 min.	Actin morphology	28
Colchicine (<i>microtubule destabilizer</i>)	10 µg/ml	Swiss 3T3 cells, adherent or suspension	Maximal activation of 2-4 fold activation after 30 min	Rhotekin-RBD pulldown	29
Cytochalasin D (<i>actin filament destabilizer</i>)	0.5 µg/ml	Swiss 3T3 cells, adherent or suspension	Maximal activation of 2-3 fold after 60 min	Rhotekin-RBD pulldown	29
Fibronectin (<i>extracellular matrix protein</i>)	Culture plate is coated with fibronectin	Swiss 3T3 cells	Biphasic regulation after plating cells on fibronectin-coated plates. Initial period of low RhoA activity (10-20 min) followed by a 2-7 fold activation peaking at 60-90 min and then dropping to basal levels after 6 h.	Rhotekin-RBD pulldown	29

Section II: Preparation of Lysates for G-LISA

Table 3H cont.: Modulators of GTP-Rho levels in cell

RhoA Modulator	Treatment	Cell type used	Response	Assay Type	Ref.
Lysophosphatidic acid (LPA) <i>(serum lipid & G-protein coupled receptor agonist)</i>	1 µg/ml	Swiss 3T3 cells, adherent & suspension	Maximal activation of 2-6 fold after 1 min then dropping to basal after 30 min	Rhotekin-RBD pulldown	29
Lysophosphatidic acid (LPA)	1 µM	N1E-115 neuronal cells	Maximal activation of 3-5 fold after 3 min	Rho-kinase pull down assay	30
Nocodazole <i>(microtubule destabilizer)</i>	10 µM	MG63 human osteosarcoma cell & HeLa cells	Maximal activation of 2-3 fold activation after 30 min	Actin morphology & Rhotekin-RBD pulldown	31
Serum	5 - 10%	Swiss 3T3 cells, adherent or suspension	Maximal activation of 2-6 fold (10%) & 2-3 fold (5%) after 1-5 min	Rhotekin-RBD pulldown	29
Sphingosine-1-phosphate <i>(serum lipid & G-protein coupled receptor)</i>	1 µg/ml	Swiss 3T3 cells, adherent or suspension	Maximal activation of 2-3 fold after 2 min for 3T3 cells and 20 min for HUVEC cells	Rhotekin-RBD pulldown	29, 32
Thrombin <i>(protease, G-protein coupled receptor)</i>	10 nM	HUVEC human endothelial primary cells	Maximal activation of 14 fold after 2 min, dropped to basal after 30 min	Rhotekin-RBD pulldown	32
Vinblastine <i>(microtubule destabilizer)</i>	50 µM	MG63 human osteosarcoma cells	Maximal activation after 30 min	Actin morphology	31

Section II: Preparation of Lysates for G-LISA

Preparation of lysates from 2D tissue culture cell lines

Activated GTPases are a labile entity and the bound GTP is susceptible to hydrolysis by GAPs during and after cell lysis, resulting in GTPase inactivation. Rapid processing at 4° C is essential for accurate and reproducible results. Detailed instructions for processing lysates from 2D tissue culture cells are given in the G-LISA assay Protocol manuals. The following guidelines are applicable to all G-LISA assays.

- 1) Wash cells in PBS: Normalization of total cell lysate concentration between samples is required for G-LISA assays. Washing cells in PBS allows removal of protein species from tissue culture media which could interfere with accurate lysate concentration determination.
- 2) Complete removal of PBS after washing: It is important to maintain the composition of the cell lysis buffer between samples, therefore complete removal of PBS, preferably with an aspirator, should be performed prior to cell lysis.
- 3) Use cold reagents for processing cells: The PBS used for washing cells and cell lysis buffer should be at 4°C. The cold buffers will help minimize hydrolysis of GTPases during processing. It should be noted, in some cell lines (e.g. 3T3 cells) Cdc42 activation appears to be affected by the temperature of the wash solution. In such cases room temperature PBS may be used. The need to use room temperature PBS should be determined empirically. In all cases, cold lysis buffer supplemented with protease inhibitors should be used.
- 4) It is recommended to process a "Test plate" prior to the experimental plates. This is used to determine the volume of lysis buffer required to achieve the lysate concentrations recommended in Table 1.
- 5) Work quickly to process lysates: GTPase proteins are labile entities and will hydrolyze over time. It is recommended to take no more than 15 minutes to process a sample and to process samples sequentially.
- 6) Snap freeze aliquots of lysate in liquid nitrogen: It is often impractical to process fresh lysates and use them immediately in an activation assay. In particular, if time-points are being taken or multiple samples are being processed, parallel processing is impractical and unnecessary. We highly recommend sequential preparation of samples followed by protein concentration determination, aliquoting and snap freezing aliquots in liquid nitrogen for later analysis. Lysates are stable for several months at -70°C.
- 7) Keep lysates at 4°C during processing: Lysates should be embedded in ice during processing. This helps to minimize changes in signal over time.
- 8) Thaw lysates quickly and DO NOT re-freeze: Thawing of cell lysates prior to use in the G-LISA[®] assay should be in a room temperature water bath, followed by rapid transfer to ice and immediate use in the assay. Transfer to ice while 1/3 of lysate remains frozen as remainder will thaw quickly on ice. Discard any unused lysate.

Section II: Preparation of Lysates for G-LISA

Preparation of lysates from 3D tissue culture cell lines

Cells grown in Matrigel or collagen gels can be assayed using the G-LISA format. The low amounts of cells in 3D cultures makes the conventional pull-down assay very difficult, variable and expensive.

Keely et al., 2007 (Methods in Enzymology, v426, p27) and Petroll et al., 2008 (J Cell Physiol, v217, p162) give in-depth descriptions using G-LISA technology and 3-D cultures. An additional modification we suggest is to increase the ratio of extract/binding buffer from the recommended 1:1 (v/v) to 1:3

Preparation of lysates from tissue samples

Tissue lysates can be used in G-LISA assays. Recommendations regarding tissue lysates are given below.

- 1) Ras superfamily GTPases are labile proteins that will hydrolyze bound GTP during sample handling. Tissues should therefore be processed quickly (2 min) and at 4°C. Tissues should be processed immediately using 4°C buffers or cut into small chunks (1 mm diameter), snap frozen in liquid nitrogen, and stored at -70°C for later processing.
- 2) Tissues can be homogenized using a micro-pestel on ice. Homogenates should be clarified by a 1 minute centrifugation at 4°C. Lysates can be used immediately in an activation assay or snap frozen in “experiment-sized” volumes. We strongly recommend snap-freezing lysates in liquid nitrogen.
- 3) When possible, tissues should be extracted in the respective G-LISA’s cell lysis buffer as this is the recommended buffer for the respective G-LISA assay (see Table 1).
- 4) It is recommended that lysis buffer be supplemented with protease inhibitors and phosphatase inhibitors. Recommended inhibitors include: Cytoskeleton protease inhibitor cocktail (Cat# PIC02), sodium fluoride (50 mM final), sodium pyrophosphate (20 mM final), p-Nitrophenyl phosphate (1 mM final), and microcystin LR (1 µM final).

Section II: Preparation of Lysates for G-LISA

Equalization of Cell Lysate concentrations

Equal protein concentration in all samples is a prerequisite for accurate comparison between samples in GTPase activation assays. Cell extracts should be equalized with ice-cold Lysis Buffer containing protease inhibitor cocktail to give identical protein concentrations. See Table 1 for recommended lysate protein concentrations for each G-LISA. Equalization is not necessary if the variation between samples is less than 10%.

The Precision Red Advanced Protein Assay Reagent, supplied in all G-LISA kits, is a simple one step procedure that results in a red to purple/blue color change characterized by an increase in absorbance at 600 nm.

The assay exhibits low variance in readings between different proteins of the same concentration and high reproducibility of the colorimetric response. The assay can also be performed in approximately 1-2 minutes. These properties are particularly valuable when applied to the labile lysates required for activation assays.

Quick Protein Concentration Method for 1 ml Cuvette (recommended)

- Add 20 μ l of each lysate or Lysis Buffer into disposable 1 ml cuvettes.
- Add 1 ml of Precision Red™ Advanced Protein Assay Reagent to each cuvette.
- Incubate for 1 min at room temperature.
- Blank spectrophotometer with 1 ml of Precision Red plus 20 μ l of Lysis Buffer at 600 nm.
- Read absorbance of lysate samples.
- Multiply the absorbance by 5 to obtain the protein concentration in mg/ml

Example Calculation

Assume a 20 μ l sample of cell lysate added to 1 ml of Precision Red reagent gives an absorbance reading of 0.1.

$$C = \frac{A}{\epsilon l} = \frac{0.1}{10 \times 1} \times 50 = 0.5 \text{ mg/ml}$$

Where c = protein concentration (mg/ml), A = absorbance reading, l = pathlength (cm), ϵ = extinction coefficient ($[\text{mg/ml}]^{-1} \text{cm}^{-1}$) and the multiplier of 50 is the dilution factor for the lysate in Precision Red reagent (20 μ l lysate in 1 ml Precision Red reagent).

Thus, for a 20 μ l sample in 1 ml of Precision Red, the equation becomes $C = A \times 5$

For a 10 μ l sample in 1 ml Precision Red, the equation becomes $C = A \times 10$

Section II: Preparation of Lysates for G-LISA

Quick Protein Concentration Method for 96 Well Plate

- Add 10 μ l of each lysate or Lysis Buffer into the well of a 96 well plate.
- Add 290 μ l of Precision Red™ Advanced Protein Assay Reagent to each well.
- Incubate for 1 min at room temperature.
- Blank spectrophotometer with 290 μ l of Precision Red plus 10 μ l of Lysis Buffer at 600 nm.
- Read absorbance of lysate samples.
- Multiply the absorbance by 3.75 to obtain the protein concentration in mg/ml

96 Well Plate Method

The linear range of this assay is 0.05 - 0.4 and is recommended when lysates are below the linear range of the 1 ml cuvette method. The pathlength for 96 well plate readings is 0.8 cm, hence the equation is modified as shown in the example below:

Example Calculation for 96 Well Plate Measurement

Assume a 10 μ l sample of cell lysate added to 290 μ l of Precision Red reagent gives an absorbance reading of 0.1

$$C = \frac{A}{\epsilon l} = \frac{0.1}{10 \times 0.8} \times 30 = 0.375 \text{ mg/ml}$$

Where c = protein concentration (mg/ml), A = absorbance reading, l = pathlength (cm), ϵ = extinction coefficient ($[\text{mg/ml}]^{-1} \text{ cm}^{-1}$) and the multiplier of 30 is the dilution factor for the lysate in Precision Red reagent (10 μ l lysate in 290 μ l Precision Red reagent).

Thus, for a 10 μ l sample in 290 μ l Precision Red, the equation becomes $C = A \times 3.75$

For a 5 μ l sample in 295 μ l Precision Red, the equation becomes $C = A \times 7.5$

NOTE: The protein concentrations generated by the multipliers stated in the method will result in approximate lysate concentrations. Data will be highly reproducible from lysate to lysate and will generate excellent values for relative concentrations of a series of lysates. It should be noted for activation assays, the relative protein concentration between experimental extracts is far more important than the absolute protein quantitation. However, if desired, one can create a standard curve using BSA or IgG protein standards for each experiment. The standard curve should be performed prior to lysate preparations due to the labile nature of the lysates.

Section III: Assay Technical Tips

Use of a multichannel pipettor

When processing more than 16 wells, it is recommended to use a multi-channel or multi-dispensing pipettor with a range of 25 to 200 μ l per dispense. Critical steps such as lysate addition, post-binding wash step and the Antigen Presenting Buffer step all have requirements for accurate and timely additions. Attempting to perform >16 assays with a single channel pipettor will also increase the likelihood of allowing wells to dry out before reagent addition can be completed, resulting in variable signals. Therefore, use a multi-channel or a multi-dispensing pipettor wherever possible.

Plate shaker recommendations

It is recommended to use an orbital plate shaker at 400 rpm. As a back-up you can use a 200 rpm orbital shaking culture incubator or a normal orbital rotating platform set to 200 rpm. Signals may be lower with the 200 rpm option. Rocking or tilting plate shakers will not be sufficient for this assay. Below are some examples of acceptable plate shakers:

Model # 4625, Titer Plate Shaker, Lab-Line Instruments, Barnstead Intl.

Model # RF7854, Digital Microplate Shaker, ML Market Lab, researchml.com

Model # RF7855, Incubating Microplate Shaker, ML Market Lab, researchml.com

Vortex of samples after binding buffer addition

Binding buffer is a viscous solution and requires thorough mixing after addition to cell lysates. The best way to do this is a brief 3-5 second vortex on high. This step will greatly reduce variability between duplicate samples and will increase overall assay accuracy.

Spectrophotometer / Luminometer settings

Spectrophotometer: The majority of the work in the design of the colorimetric G-LISA assays has been based on the Molecular Devices SpectraMax 250. The spectrophotometer settings are given in Table 4 as a reference.

Table 4: Spectrophotometer settings for colorimetric G-LISAs

Parameters	Character	Contents
Wavelength	490 nm	Bandwidth 2 nm (can be \pm 20 nm for filter based machines)
Protocol	End point	Standard end point assay
Shaking	Medium, orbital	5 s
Temperature	24°C	Room temperature is also fine for readings

Section III: Assay Technical Tips

Luminometer: Luminometers vary widely in their sensitivity and absolute readings. It is therefore recommended to run a G-LISA assay with blank and positive control to determine that you are in the linear range of the assay. When in the linear range of the assay the positive control should read 4-10 fold higher than the blank wells. Table 5 gives guidelines for luminometer settings;

Table 5: Luminometer settings for luminescence based G-LISAs

Parameters	Description and Recommendations
Gain	<p>Gain controls the sensitivity of the machine. Most luminometers do not allow manual alteration of gain and use an auto-calibration or limited calibration function. It is important to contact the luminometer manufacturer or consult the users manual to determine the best way to alter the machine sensitivity.</p> <p>If gain can be altered one should read at low, medium and high gains to determine the reading within the linear range of the assay (positive control should be 4-7X higher than blank). Gain range varies with instrument, for example gain in the Tecan GmbH SpectroFluor Plus ranges from 0 - 150 (where 150 is the highest).</p>
Integration Time	<p>This parameter can be varied on most machines. It is a good idea to set the machine as the lowest integration time (usually 10 – 100 ms). Integration times greater than 200 ms are likely to read out of the linear range of the assay and may require lowering of gain or dilution of primary and/or secondary antibodies (see below).</p>
Shaking	<p>Most machines give the shaking option. The recommended setting is 5 s shake, medium orbital speed before read. This option is not essential to the assay.</p>
Temperature	<p>Room temperature</p>
Plate type	<p>Any setting that specifies 96 well flat, white will be sufficient</p>
Filters	<p>Luminescence does not require excitation or emission filters the filter spaces should be left blank.</p> <p>If this is not an option excitation can be set at any value and emission should be set between 400-500nm, with 430-445 as optimal setting.</p>

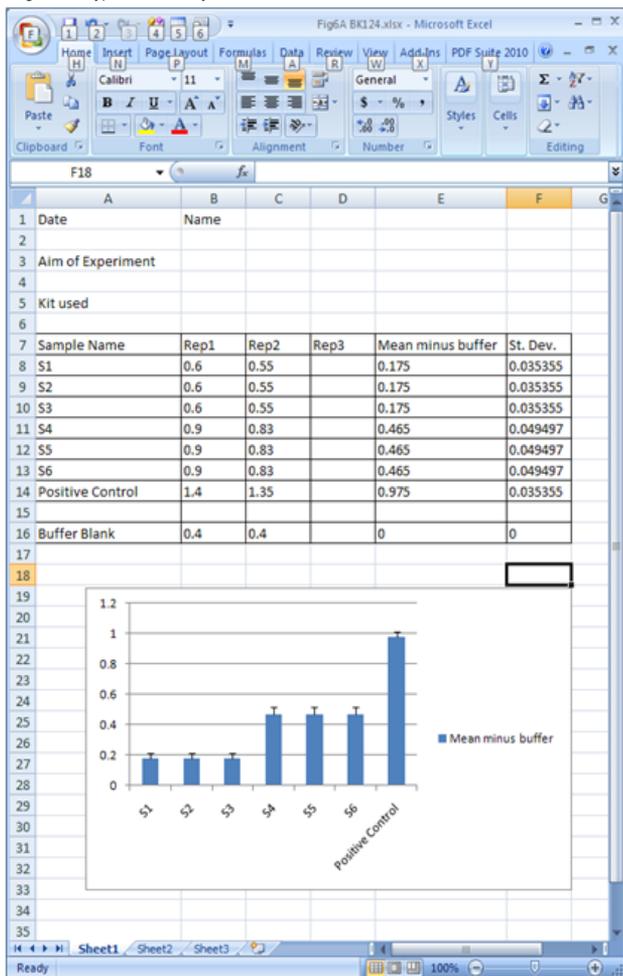
Section IV: Data Analysis

Data Analysis: Excel format

1. It is recommended to use the Lysis Buffer wells as reference blanks in all studies with this kit. Based on the operator designating the appropriate wells, most machines have associated protocols that perform this operation automatically. Call Technical Help for the company supplying the plate reader for information on how to perform this function. When the data are “Lysis Buffer subtracted” (Lysis Buffer only samples have been allocated as Blanks in the assay), then you can import them into a simple graph software such as Excel. Alternatively, the Lysis Buffer background can be subtracted manually or in the spreadsheet application.
2. Data should be arranged in columns where the headings are “Sample”, “Mean”, “Standard Deviation”, “rep1”, “rep2”, “rep3” and “rep4” for the number of replicates performed on each sample. E-mail tservice@cytoskeleton.com or visit www.cytoskeleton.com to download a free Excel Template.
3. List your samples under the “Sample” column in the same order that they were assayed in the plate.
4. Enter the following formula into the first sector under “Mean”, “=average(Xn:Yn)” where X = the column designator for “rep1”, Y = column designator for “rep4”, and n= row designator of the row that you are working on. Repeat for each sector under the “Mean” header until there are sufficient rows to cover the number of samples in your experiment.
5. Enter the following formula into the first sector under “Standard Deviation”, “=stdev(Xn:Yn)” where X = the column designator for “rep1”, Y = column designator for “rep4”, and n= row designator of the row that you are working on. Repeat for each sector under the “Standard Deviation” header until there are sufficient rows to cover the number of samples in your experiment.
6. Enter your replicate data into rep1, rep2 etc. It doesn’t matter if you only have duplicates because the program will ignore any sectors that do not contain data. The program will calculate the Mean and Standard Deviation of your replicates.
7. When the data has been entered select the Sample, Mean and Standard Deviation data sectors by the click and drag method. Then select the chart making function, in Excel this looks like a clickable square with a mini-bar chart inside. This will guide you through the chart making process with the data you have selected. Choose “column chart” initially, designate the Mean numbers for input values. The Standard Deviation column for the y-axis error bars needs to be designated after the Mean numbers chart is made. This is achieved by double clicking on the graph bars, and selecting the “Y-axis error” tab, then entering the location of the Standard Deviation data by clicking the “Custom” option and selecting the area in the worksheet. E-mail tservice@cytoskeleton.com or visit www.cytoskeleton.com to download a free Excel Template.
8. An example of a typical Excel layout and data plot is shown in Figure 1.

Section IV: Data Analysis

Figure 1: Typical Excel Layout



Section IV: Data Analysis

Experiment Record Sheet

Scientist Name

Contact Tel. #

e-mail

Kit Cat. # / Lot #

STEP

Comments or Changes

- 1 Type of cells or tissue
- 2 How were the cells prepared prior to lysis? days in culture
.....% confluency
..... inducer
.....mg/ml of protein in lysate
- 3 How long were the ice-cold solutions on ice before lysis? Min
- 4 Time that cultures were removed from incubator? am or pm
- 5 Was Binding Buffer added? Y or N
- 6 Time that binding reactions were placed on the shaker? am or pm
- 7 Did you add 50 μ l of extract per well? Y or N
- 8 What locations are the 50 μ l Lysis Buffer controls? Wells
- 9 What speed and time was the shaking for the binding reaction? rpmmin
- 10 How long did you wait after the post-binding wash step? s or min
- 11 What was the time when the primary antibody reaction was started?am or pm
- 12 What was the time when the Secondary antibody reaction was started?..... am or pm
- 13 What was the time when detection reagent was added? am or pm
- 14 What was the time when the plate was read? am or pm

Section IV: Data Analysis

Plate Record Template

Name of experiment:

Date of experiment:

	H	G	F	E	D	C	B	A	
1									
2									
3									
4									
5									
6									
7									
8									
9									
10									
11									
12									

Technical Assistance: call either 303-322-2254 or e-mail tservice@cytoskeleton.com

Section V: References

1. Niu T.-K. et al. 2005. Dynamics of GBF1, a brefeldin A-sensitive Arf1 exchange factor at the Golgi. *Mol. Biol. Cell.* **16**, 1213-1222.
2. Matto M. et al. 2011. Role of ADP ribosylation factor 1 in the regulation of Hepatitis C virus replication. *J. Virol.* **85**, 946-956.
3. Tanaka M. et al. 2007. The C-terminus of ephrin-B1 regulates metalloproteinase secretion and invasion of cancer cells. *J. Cell Sci.* **120**, 2179-2189.
4. Wessels E. et al. 2006. A viral protein that blocks Arf1-mediated COP-1 assembly by inhibiting the guanine nucleotide exchange factor GBF1. *Dev. Cell.* **11**, 191-201.
5. El Azreq M.-A. et al. 2010. Cytohesin-1 regulates the Arf6-phospholipase D signaling axis in human neutrophils: impact on superoxide anion production and secretion. *J. Immunol.* **184**, 637-649.
6. Zhang Q. et al. 2007. Small-molecule synergist of the Wnt/b-catenin signaling pathway. *Proc. Natl. Acad. Sci. U.S.A.* **104**, 7444-7448.
7. Belov G.A. et al. 2007. Hijacking components of the cellular secretory pathway for replication of poliovirus RNA. *J. Virol.* **81**, 558-567.
8. Balasubramanian N. et al. 2007. Arf6 and microtubules in adhesion-dependent trafficking of lipid rafts. *Nat. Cell Biol.* **9**, 1381-1391.
9. Goldfinger L. E. et al. 2006. RLIP76(RalBP1) is an R-Ras effector that mediates adhesion-dependent Rac activation and cell migration. *J. Cell Biol.* **174**, 877-888.
10. Choi W. et al. 2006. Arf6 plays an early role in platelet activation by collagen and convulxin. *Blood.* **107**, 3145-3152.
11. Hafner M. et al. 2006. Inhibition of cytohesins by SecinH3 leads to hepatic insulin resistance. *Nature.* **444**, 941-944.
12. Palacios F. & D'Souza-Schorey C. 2003. Modulation of Rac1 and ARF6 activation during epithelial cell scattering. *J. Biol. Chem.* **278**, 17395-17400.
13. Sakurai, A. et al. 2010. Semaphorin 3E initiates antiangiogenic signaling through Plexin D1 by regulating Arf6 and R-Ras. *Mol. Cell. Biol.* **30**, 3086-3098.
14. Yao et al., 2013. Nonmuscle myosin light-chain kinase mediates microglial migration induced by HIV Tat: Involvement of $\beta 1$ integrins. *FASEB J.* doi: 10.1096/fj-219600
15. Chen et al., 2012. Inhibition of tumor cell growth, proliferation and migration by X-387, a novel active-site inhibitor of mTOR. *Biochem. Pharmacol.* **83**, 1183-1194.
16. Romero et al., 2010. Chronic Ethanol Exposure Alters the Levels, Assembly, and Cellular Organization of the Actin Cytoskeleton and Microtubules in Hippocampal Neurons in Primary Culture. *Toxicol. Sci.* **118**, 602-612.

Section V: References

17. Kim H.D. et al. 2008. Epidermal growth factor-induced enhancement of glioblastoma cell migration in 3D arises from an intrinsic increase in speed but an extrinsic matrix- and proteolysis-dependent increase in persistence. *Mol. Biol. Cell.* **19**, 4249-4259.
18. Tanaka T. et al. 2010. Monocyte chemoattractant proein-1/CC chemokine ligand 2 enhances apoptotic cell removal by macrophages through Rac1 activation. *Biochem. Biophys. Res. Commun.* **399**, 677-682.
19. Yang C. et al. 2006. Essential role for Rac in heregulin beta1 mitogenic signaling: a mechanism that involves epidermal growth factor receptor and is independent of ErbB4. *Mol. Cell. Biol.* **26**, 831-842.
20. Grill B. & Schrader J.W. 2002. Activation of Rac-1, Rac-2, and Cdc42 by hemopoietic growth factors or cross-linking of the B-lymphocyte receptor for antigen. *Blood.* **100**, 3183-3192.
21. Takaya A. et al. 2004. RalA action at nascent lamellipodia of epidermal growth factor stimulated Cos7 cells and migrating madin-darby kidney cells. *Mol. Biol. Cell* **15**, 2549-2557.
22. Omerovic J. et al. 2008. Ras isoform abundance and signaling in human cancer cell lines. *Oncogene.* **27**, 2754-2762.
23. Satoh T. et al. 1993. Platelet-derived growth factor receptor mediates activation of ras through different signaling pathways in different cell types. *Mol. Cell Biol.* **13**, 3706-3713.
24. Perez de Castro I. et al. 2004. Ras activation in Jurkat T cells following low-grade stimulation of the T-cell receptor is specific to N-Ras and occurs only on the Golgi apparatus. *Mol. Cell Biol.* **24**, 3485-3496.
25. Scott G.A. et al. 2007. Lysophosphatidylcholine mediates melanocyte dendricity through PKC ζ activation. *J. Invest. Dermatol.* **127**, 668-675.
26. Aguilar H.N. et al. 2011. Phos-tag-based analysis of myosin regulatory light chain phosphorylation in human uterine myocytes. *PLoS ONE.* **6**, e20903.
27. Ridley A.J. & Hall A. 1992. The small GTP-binding protein rho regulates the assembly of focal adhesions and actin stress fibers in response to growth factors. *Cell.* **70**, 389-399.
28. Schoenwaelder S.M. & Burridge K. 1999. Evidence for a calpeptin-sensitive protein-tyrosine phosphatase upstream of the small GTPase Rho. *J. Biol. Chem.* **274**, 14359-14367.
29. Ren X.D. et al. 1999. Regulation of the small GTP-binding protein Rho by cell adhesion and the cytoskeleton. *EMBO J.* **18**, 578-585.
30. Kranenburg, O. et al. 1999. Activation of RhoA by lysophosphatidic acid and Ga12/13 subunits in neuronal cells: induction of neurite retraction. *Mol. Biol. Cell.* **10**, 1851-1857.

Section V: References

31. Zhang, Q. et al. 1997. Lysophosphatidic acid and microtubule-destabilizing agents stimulate fibronectin matrix assembly through Rho-dependent actin stress fiber formation and cell contraction. *Mol. Biol. Cell.* **8**, 1415-1425.
32. Vouret-Craviari, V. et al. 2002. Distinct signals via Rho GTPases and Src drive shape changes by thrombin and spingosine-1-phosphate in endothelial cells. *J. Cell Sci.* **115**, 2475-2484.

