



Helping advance science,  
one protein at a time.

MAR  
2012

this issue

Cardiac Research Using G-LISA Technology  
G-LISA Publications  
Rho Protein Research Tools

## Upcoming Meetings

American Association for  
Cancer Research (AACR)  
March 31st - April 4th, 2012  
Booth # 4207

## Cytoskeleton Products

Actin Proteins  
Activation Assays  
Antibodies  
ECM Proteins  
ELISA Kits  
G-LISA® Kits  
Pull-down Assays  
Motor Proteins  
Small G-Proteins  
Tubulin & FtsZ Proteins

## Contact Us

P: 1 (303) 322.2254  
F: 1 (303) 322.2257  
E: [cserve@cytoskeleton.com](mailto:cserve@cytoskeleton.com)  
W: [cytoskeleton.com](http://cytoskeleton.com)

## Distributors

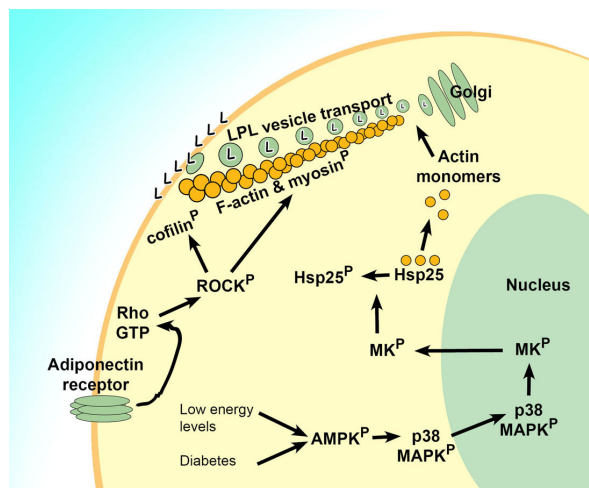
[www.cytoskeleton.com/distributors/](http://www.cytoskeleton.com/distributors/)

## Cardiac research using G-LISA® technology: Studying the Rho pathway in diabetic cardiomyopathy.

The Rho/ROCK pathway is understood to regulate cardiac hypertension in cardiovascular disease<sup>1</sup>. The main manifestation is to regulate the tension of the actomyosin cytoskeleton<sup>2,3,4,5</sup>. It's also well documented that inducing RhoA activation causes rearrangement of the actin cytoskeleton in many cell types. In cardiac tissue, one consequence of this re-organization leads to increased localization of lipoprotein lipase (LPL) to the plasma membrane<sup>6,7</sup>, which causes increased reliance on fatty acid metabolism (breakdown of chylomicrons and very low-density lipoproteins) instead of glucose oxidation<sup>8</sup>. Similar physiological effects occur in the hyperglycemic state of type 2 diabetes, which exacerbates the poor metabolic state of cells in this disease<sup>7,8</sup>. Rho activation is a focal point in the pathway from extracellular signals to intracellular mobilization of the cytoskeleton, and hence, re-localization of LPL depends on Rho activation<sup>6</sup>. There are believed to be two pathways that cause LPL mobility to the plasma membrane these are the Adiponectin Receptor - RhoA - Rock pathway and the AMPK - p38 - MK pathway (see Figure 1). The normal physiological cause of LPL re-localization is in response to low energy levels which are detected by AMP-activated protein kinase. This kinase phosphorylates p38 MAPK which translocates to the nucleus and increases transcription of MAPK-activated protein kinase 2 (MK2). MK2 phosphorylates HSP25 which uncouples from actin monomers creating a larger pool of free actin for assembly<sup>8</sup>. In concert with this cascade, circulating in the blood, adiponectin binds to its extracellular receptor activating the RhoA/ROCK pathway, which causes F-actin fiber formation with an appropriate orientation towards the plasma membrane<sup>6</sup>.

The RhoA G-LISA format offered by Cytoskeleton, Inc. has enabled the measurement of Rho activity in cardiology studies despite the small amounts of primary cardiac myocytes or aorta tissues that are available<sup>4,6,9</sup>. The benefits of this format are 1) improved accuracy (cv = 13%) over the traditional Western blot approach<sup>10</sup>, 2) rapid processing, 3) small sample size (10-50 µg total protein),

Figure 1. Signalling pathways regulating LPL re-localization.



*Legend: Extracellular factors such as adiponectin and intracellular cues such as low energy levels initiate the signalling pathways through AMPK and Rho/Rock to create a cytoskeletal framework to transport LPL from the Golgi to the plasma membrane.*

and 4) improved economy per assay. These strengths have allowed researchers to measure Rho activity in smaller pieces of tissue such as mouse aorta, which is 2 mm wide and 8 mm long and too small to be measured by previous methods<sup>4,6,9</sup>. In addition, studying RhoA in aorta homogenates has been a challenge due to high levels of contaminating phosphatases which reduce the GTP-Rho signals to indistinguishable levels in some circumstances. Thus, aorta tissue homogenates require phosphatase inhibitors in order to stabilize the GTP-RhoA signal long enough to measure with G-LISA. Appropriate phosphatase inhibitors are a mixture of false substrates and inhibitors of serine/threonine phosphatases. We recommend 50 mM NaF,

## References

20 mM Na-pyrophosphate, 1  $\mu$ M microcystin and 1 mM ortho-vanadate. Without these inhibitors, some researchers have found samples produce very similar and low (degraded) GTP-RhoA signals.

Finally, Rho specific inhibitors and activators are an emerging trend in the clinic, such as Bioaxone's BA-210 (Cethrin®) which is indicated for neuro-regeneration. Cytoskeleton offers a Rho inhibitor (Cat. # CT04) and a Rho Activator (Cat. # CN03) which are highly potent cell permeable reagents for Rho pathway research.

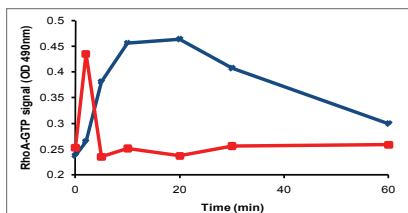
### References:

- Hong Z and Yong-Jun L. 2010. RhoA/Rho kinase: a novel therapeutic target in diabetic complications. *Chinese Medical Journal*. 123: 2461-2466.
- Hirata et al., 1992. Involvement of Rho p21 in the GTP-enhanced calcium ion sensitivity of smooth muscle contraction. *J. Biol. Chem.*, **267**, p.8719-8722.

- Gong et al., 1995. Arachidonic acid and diacylglycerol release associated with inhibition of myosin light chain dephosphorylation... *J. Physiol.*, **486**, p. 113-122.
- Matsumoto et al., 2010. Enhancement of mesenteric artery contraction to 5-HT depends on Rho kinase and Src kinase pathways in the ob/ob mouse model of type 2 diabetes. *Br J Pharmacol.*, **160**(5): 1092-1104.
- Yang et al., 2012. Mechanism of fibrotic cardiomyopathy in mice expressing truncated Rho-associated coiled-coil protein kinase 1. *FASEB J.* Jan 25th e-pub ahead of print.
- Ganguly et al., 2011. Adiponectin increases LPL activity via RhoA/ROCK-mediated actin remodeling in adult rat cardiomyocytes. *Endocrinology*, **152**, p.247-254.
- Punlinilkunnit and Rodrigues, 2006. Cardiac lipoprotein lipase: Metabolic basis for diabetic heart disease. *Cardiovasc. Res.*, **69**, p.329-340.
- Kim et al., 2007. Acute diabetes moderates trafficking of cardiac lipoprotein lipase through p38 mitogen-activated protein kinase-dependent actin cytoskeleton organization. *Diabetes*, **57**, p.64-76.
- Seok et al., 2008. Isoflavone Attenuates Vascular Contraction through Inhibition of the RhoA/Rho-Kinase Signaling Pathway. *J. Pharmacol. Exp. Ther.*, **326**, p.991-998.
- Benard and Bokoch, 2002. Assay of Cdc42, Rac, and Rho GTPase activation by affinity methods. *Methods Enzymol.*, **345**, p.349-359.

## Rho Protein Research Tools

### RhoA G-LISA and ELISA



**Time course of activation of RhoA in Swiss 3T3 cells by CN01 and LPA.** Serum-starved Swiss 3T3 cells were treated with Rho Activator, cat. # CN01 (blue diamonds) or LPA (magenta squares). RhoA activity was measured by reading signals at OD<sub>490nm</sub>.

Kits	Cat #	Amount
RhoA G-LISA® Activation Assay, colorimetric	BK124	96 assays
RhoA G-LISA® Activation Assay, luminescence	BK121	96 assays
Total RhoA ELISA	BK150	96 assays

### GEF and GAP Assays

Products	Cat. #	Amount
RhoGAP Assay Biochem Kit™	BK105	80-160 assays
RhoGEF Exchange Assay Biochem Kit™	BK100	60-300 assays

### Rhotekin-RBD Beads

Products	Purity	Cat. #	Amount
Rhotekin-RBD Protein	>90%	RT01-A RT01-B	1 x 500 $\mu$ g 3 x 500 $\mu$ g
Rhotekin-RBD Beads	>85%	RT02-A RT02-B	2 x 2 mg 6 x 2 mg

### RhoA Antibody

Product	Host	Type	Species Reactivity	Cat. #	Amount
RhoA Specific Antibody	Mouse	mAb	Hu, Ms, Rt, other extracts	ARH03-A ARH03-B	1 x 100 $\mu$ g 3 x 100 $\mu$ g
Human RhoA Peptide (a.a. 120-150)					

### Rho Activators & Inhibitors

G-protein Modulator	Cell Entry Mechanism	Protein Modulation	Cat. #	Amount
<b>Rho Activator II</b> Deamidation of Rho Gln-63	Cell permeable	Direct	CN03-A CN03-B	3 x 20 $\mu$ g 9 x 20 $\mu$ g
<b>Rho Inhibitor I</b> ADP ribosylation of Rho Asn-41	Cell permeable	Direct	CT04-A CT04-B	1 x 20 $\mu$ g 5 x 20 $\mu$ g
<b>Rho/Rac/Cdc42 Activator I</b> Deamidation of Rho Gln-63 & Rac/Cdc42 Gln-61	Cell permeable	Direct	CN04-A CN04-B	3 x 20 $\mu$ g 9 x 20 $\mu$ g
<b>Rho Pathway Inhibitor I</b> Rho kinase (ROCK) inhibitor Y-27632	Cell permeable	Direct	CN06-A CN06-B	5 x 10 units 20 x 10 units
<b>Rho Activator I</b> SHP-2 phosphatase-mediated Rho activation	Cell permeable	Indirect	CN01-A CN01-B	5 x 10 units 20 x 10 units

### Rho Modulator & Effector Proteins

Products	Purity	Cat. #	Amount
<b>C3 Transferase Protein</b> Specific inhibitor of Rho activity (not cell permeable, but also see CT04 above).	>90%	CT03-A CT03-B CT03-C	1 x 25 $\mu$ g 2 x 25 $\mu$ g 4 x 25 $\mu$ g
<b>Dbs His Protein, RhoGEF domain (DH/PH)</b> GEF for Cdc42 and RhoA	>80%	GE01-A GE01-B	2 x 50 $\mu$ g 8 x 50 $\mu$ g
<b>p50RhoGAP GST Protein, full length</b> GAP for Cdc42, Rac and Rho	>90%	GAP01-A GAP01-B	1 x 50 $\mu$ g 4 x 50 $\mu$ g
<b>p50RhoGAP GST Protein, GAP domain</b> GAP for Cdc42, Rac and Rho	>90%	GAS01-A GAS01-B	1 x 50 $\mu$ g 4 x 50 $\mu$ g
<b>RhoGDI GST Protein</b> Inhibitor of Cdc42, Rac and Rho	>90%	GDI01-A GDI01-B	1 x 25 $\mu$ g 4 x 25 $\mu$ g

### Rho Purified Proteins

Purified G-proteins	Purity	Cat. #	Amount
<b>RhoA His Protein, constitutively-active (Q63L)</b>	>90%	R6301-A R6301-C	1 x 10 $\mu$ g 4 x 10 $\mu$ g
<b>RhoA GST Protein, wild-type</b>	>90%	RHG01-C	8 x 25 $\mu$ g
<b>RhoA His Protein, wild-type</b>	>80%	RH01-A RH01-C RH01-XL	1 x 100 $\mu$ g 3 x 100 $\mu$ g 1 x 1 mg
<b>RhoC His Protein, wild-type</b>	>90%	RH03-A RH03-C	1 x 100 $\mu$ g 3 x 100 $\mu$ g