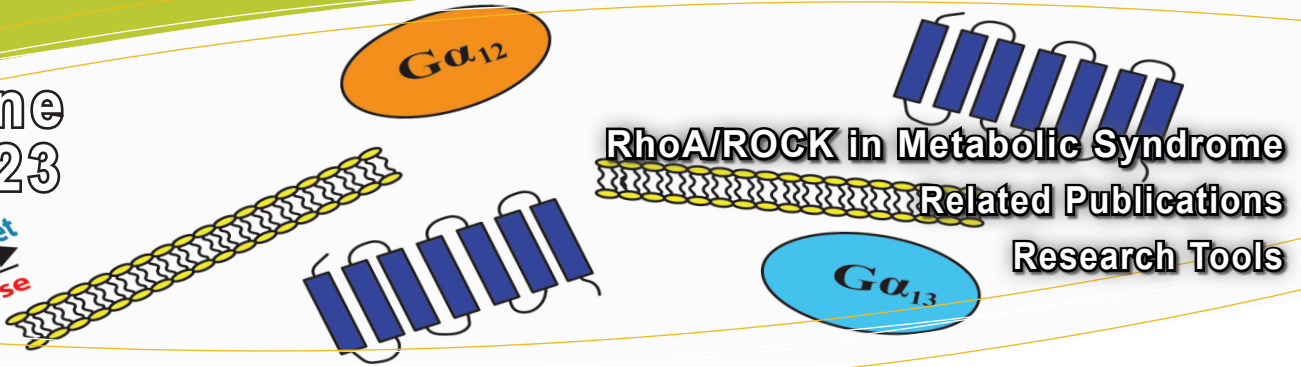




Helping advance science,  
one protein at a time.

June  
2023

High Fat Diet  
↔  
Exercise



## Sponsored Conferences

Gordon Scientific  
Inhibition in the CNS  
Les Diablerets,  
Switzerland  
July 16<sup>th</sup>-21<sup>st</sup>

Kern Lipid Conference  
Vail, CO  
August 14<sup>th</sup>-16<sup>th</sup>

## Cytoskeleton Products

- Actin Proteins
- Activation Assays
- Antibodies
- ECM Proteins
- ELISA Kits
- G-LISA® Kits
- Live Cell Imaging
- 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)

## For More News

W: [cytoskeleton.com /blog/](http://cytoskeleton.com/blog/)

## RhoA/ROCK in Metabolic Syndrome

### Introduction

Metabolic syndrome (MetS) or insulin resistance syndrome, is a cluster of conditions that raises the risk of heart disease, type II diabetes, and stroke. These conditions include central obesity (i.e., excess body fat at the waist), hypertension, hyperglycemia, high blood triglycerides, and low HDL cholesterol. Those with three or more of these conditions may be diagnosed with MetS. MetS affects approximately one-third of adults in the United States. Pathogenesis of MetS is still not fully understood, though a variety of genetic, environmental, and lifestyle factors are thought to be involved.

Rho GTPases are a family of signaling proteins that play a key role in regulating the cytoskeleton. Activation of RhoA by GTP binding results in the activation of downstream effector proteins that mediate actin polymerization and stabilization and myosin II function<sup>1</sup>. RhoA is involved in essential cellular processes, such as cytokinesis, stress fiber formation, and smooth muscle cell contractility, as well as several pathological processes, including tumor cell migration and invasion and cardiovascular diseases<sup>2-5</sup>. RhoA and its downstream effector proteins, Rho-associated protein kinases (ROCK1 and ROCK2), are also critical in MetS pathogenesis<sup>6</sup>. Elevated ROCK activity has been demonstrated in models of MetS conditions, including obesity, insulin resistance, elevated triglycerides, and hypertension<sup>4,7</sup>. In the PBMCs of patients with type 2 diabetes mellitus, ROCK activity is elevated 420-570% compared to healthy controls<sup>8</sup>. Serum ROCK activity is also elevated in those diagnosed with MetS and correlated with waist circumference, fasting glucose, and triglyceride levels<sup>9</sup>. As a result, RhoA and ROCK have become attractive therapeutic targets for MetS treatment<sup>10,11</sup>. One randomized controlled trial revealed that atorvastatin inhibits ROCK activity, which they suggest contributes to the anti-atherosclerotic benefit of statins<sup>12</sup>. This newsletter will focus on recently identified roles for RhoA and ROCK signaling in the biological processes involved in MetS conditions.

### RhoA/ROCK Signaling Mediates Actin Cytoskeleton Dynamics in Adipogenesis

There are three main types of adipocytes. White adipocytes store excess lipids as triglycerides and release free fatty acids during times of high energy expenditure<sup>13</sup>. Brown adipocytes are thermogenic, that is, they convert energy into heat<sup>14</sup>. Brown adipocytes decrease with age, while overall adiposity, particularly white adipocyte accumulation, increases with age. Adipocytes with intermediate characteristics are referred to as beige adipocytes and can switch between the functionality of white and brown adipocytes based on

current energy conditions<sup>13,14</sup>. Brown and beige adipocytes are critical for maintaining energy homeostasis and their decline is associated with an increased risk of obesity, diabetes, and other metabolic disorders<sup>15</sup>. For this reason, identifying signaling components that can modulate the adipogenesis program to favor thermogenic adipocytes is a promising therapeutic strategy for MetS treatment.

RhoA and ROCK play a key role in regulating actin cytoskeleton dynamics in all metabolic tissues and during adipogenesis, the process of creating new fat cells<sup>6,16</sup>. RhoA/ROCK activity inhibits the differentiation of white and beige adipocytes as well as the transdifferentiation of white adipocytes to beige<sup>6</sup>. Rho and ROCK activity further promote obesity by shifting away from thermogenic brown and beige adipogenesis in favor of myogenesis and cardiomyocyte differentiation, respectively<sup>6</sup>. Thus, targeting RhoA/ROCK activity represents a means to shift toward brown and beige adipogenesis by affecting multiple points of the adipogenesis program. Indeed, ROCK2 inhibition in mice enhances the thermogenic program in both white and brown adipocytes<sup>17</sup>. Another study demonstrated that mice deficient in the transcription factor MRTFA, which lies downstream of ROCK signaling, were protected from diet-induced obesity and demonstrated "beiging" of white adipose tissue<sup>18</sup>.

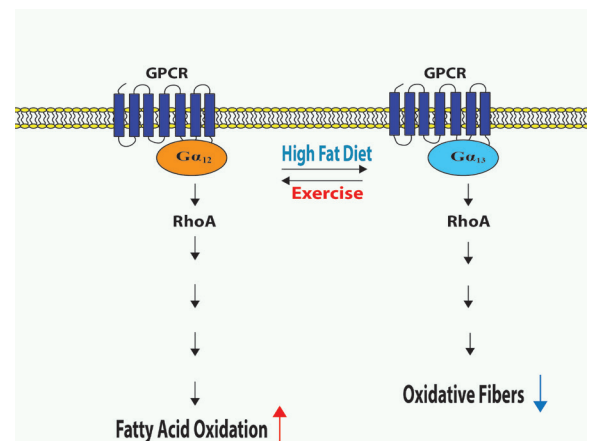


Figure 1. Schematic of how distinct GPCRs induce RhoA activation resulting in opposing metabolic regulatory mechanisms.



Helping advance science,  
one protein at a time.

## Continued from Page 1

### G<sub>α<sub>12,13</sub></sub>/RhoA Affect Obesity, Whole-Body Energy Metabolism, Insulin Sensitivity, and Liver Steatosis

Upstream of RhoA, heterotrimeric G-proteins containing G<sub>α<sub>12/13</sub></sub> subunits transduce signals from a wide range of G protein-coupled receptors (GPCRs), including sphingosine 1-phosphate (S1P) receptors, thrombin receptors, and purinergic receptors<sup>19</sup>. Significant evidence indicates that GPCR signaling through G<sub>α<sub>12/13</sub></sub> influences a wide-range of metabolic disorders<sup>19-21</sup>. Though both G<sub>α<sub>12</sub></sub> and G<sub>α<sub>13</sub></sub> lead to the activation of RhoA, they appear to exert their effects in a tissue-specific manner and serve opposing roles in the regulation of pathways involved in energy homeostasis in certain tissues<sup>19</sup> (Figure 1).

In mice, exercise reduces G<sub>α<sub>13</sub></sub> levels in skeletal muscle while high-fat diet (HFD)-fed mice have elevated G<sub>α<sub>13</sub></sub> levels. Similarly, G<sub>α<sub>13</sub></sub> is elevated in the skeletal muscle of patients with type II diabetes mellitus<sup>19</sup>. Skeletal muscle-specific ablation of G<sub>α<sub>13</sub></sub> in mice leads to higher levels of whole-body metabolism and increased insulin sensitivity<sup>22</sup>. This effect is dependent on G<sub>α<sub>13</sub></sub>/RhoA/ROCK2-mediated suppression of nuclear factor of activated T cells 1 (NFATc1) by phosphorylation at Ser243<sup>22</sup>. Phosphorylation of NFATc1 at this site results in its inactivation and is elevated in obese mice and reduced following exercise<sup>23</sup>. While G<sub>α<sub>13</sub></sub>-mediated activation of RhoA in skeletal muscle leads to ROCK-mediated diet-induced adiposity, G<sub>α<sub>12</sub></sub>-mediated activation of RhoA leads to HIF-1α-dependent fatty acid oxidation<sup>19,22,23</sup>. Mice fed a HFD and patients with liver steatosis have lower levels of liver G<sub>α<sub>12</sub></sub> while expression is upregulated under fasting conditions<sup>23</sup>. Ablation of G<sub>α<sub>12</sub></sub> in mice leads to higher rates of lipid accumulation<sup>23</sup>. Mechanistically, this role for G<sub>α<sub>12</sub></sub> in the regulation of fatty acid oxidation and mitochondrial respiration is mediated by the SIRT1/PPARα pathway<sup>23,24</sup>. Thus, in skeletal muscle, G<sub>α<sub>12</sub></sub>/RhoA and G<sub>α<sub>13</sub></sub>/RhoA serve opposing roles and together maintain energy homeostasis.

This dual regulatory program is absent in the liver where NFATc1 is not expressed. In contrast to skeletal muscle expression levels, hepatic G<sub>α<sub>13</sub></sub> levels are lower in HFD-fed or genetically obese mice and patients with diabetes<sup>19</sup>. Additionally, hyperglycemia in mice leads to decreased hepatic G<sub>α<sub>13</sub></sub> levels, contributing to glucose intolerance and insulin resistance<sup>25</sup>. In the liver, G<sub>α<sub>12</sub></sub> contributes to hepatic fibrosis by promoting JNK-dependent autophagy in hepatic stellate cells (HSCs)<sup>19</sup>.

### Summary and Future Perspective: G<sub>α<sub>12/13</sub></sub>/RhoA/ROCK as Therapeutic Targets in MetS

Though clinical trials to evaluate the pharmacological modulation of these signaling components for MetS are not yet underway, many preclinical animal studies have underscored the opportunities that exist. In one such example, the inhibition of ROCK suppressed obesity, hypercholesterolemia, and glucose intolerance in mice through the activation of the LKB1/AMPK pathway<sup>26</sup>. Because RhoA downstream effectors and subsequent cellular processes are tissue-specific, targeted therapeutic approaches such as liver-specific ROCK1 inhibition, have been evaluated in preclinical mouse models, resulting in increased energy expenditure, greater insulin sensitivity, and reduced lipid accumulation<sup>27</sup>. The multifaceted roles of G<sub>α<sub>12/13</sub></sub>/RhoA/ROCK signaling in metabolic syndrome continue to be uncovered. Recent evidence suggests this signaling axis represents a novel therapeutic target for MetS treatment based on several preclinical studies that have investigated RhoA signaling components as therapeutic targets.

## RhoA Small G-Proteins

Labeled Actin	Amount	Cat. #
<b>RhoA Protein: His Tagged Human Wild Type</b>	1 x 100 µg	RH01-A
	3 x 100 µg	RH01-C
	1 x 1 mg	RH01-XL
<b>RhoA Protein: His Tagged: Human Constitutively Active</b>	1 x 10 µg	R6301-A

## References

- Lessey EC, Guilluy C, Burridge K. From mechanical force to RhoA activation. *Biochemistry*. 2012;51(38):7420-7432. doi:10.1021/bi300758e
- Basant A, Glotzer M. Spatiotemporal Regulation of RhoA during Cytokinesis. *Curr Biol*. 2018;28(9):R570-R580. doi:10.1016/j.cub.2018.03.045
- Ridley AJ, Hall A. The small GTP-binding protein rho regulates the assembly of focal adhesions and actin stress fibers in response to growth factors. *Cell*. 1992;70(3):389-399. doi:10.1016/0092-8674(92)90163-7
- Duong-Quy S, Bei Y, Liu Z, Dinh-Xuan AT. Role of Rho-kinase and its inhibitors in pulmonary hypertension. *Pharmacol Ther*. 2013;137(3):352-364. doi:10.1016/j.pharmthera.2012.12.003
- O'Connor K, Chen M. Dynamic functions of RhoA in tumor cell migration and invasion. *Small GTPases*. 2013;4(3):141-147. doi:10.4161/sqtp.25131
- Wei L, Shi J. Insight Into Rho Kinase Isoforms in Obesity and Energy Homeostasis. *Front Endocrinol (Lausanne)*. 2022 Jun 13;13:886534. doi:10.3389/fendo.2022.886534
- Lee SH, Huang H, Choi K, et al. ROCK1 isoform-specific deletion reveals a role for diet-induced insulin resistance. *Am J Physiol Endocrinol Metab*. 2014;306(3):E332-E343. doi:10.1152/ajpendo.00619.2013
- Ocaranza, M.P, Valderas, P, Moya, J, et al. Rho kinase cascade activation in circulating leukocytes in patients with diabetes mellitus type 2. *Cardiovasc Diabetol* 19, 56 (2020). <https://doi.org/10.1186/s12933-020-01027-2>
- Liu PY, Chen JH, Lin LJ, Liao JK. Increased Rho kinase activity in a Taiwanese population with metabolic syndrome. *J Am Coll Cardiol*. 2007;49(15):1619-1624. doi:10.1016/j.jacc.2006.12.043
- Jahani V, Kavousi A, Mehri S, Karimi G. Rho kinase, a potential target in the treatment of metabolic syndrome. *Biomed Pharmacother*. 2018 Oct;106:1024-1030. doi:10.1016/j.biopha.2018.07.060. Epub 2018 Jul 14.
- Hahmann C, Schroeter T. Rho-kinase inhibitors as therapeutics: from pan inhibition to isoform selectivity. *Cell Mol Life Sci*. 2010;67(2):171-177. doi:10.1007/s00018-009-0189-x
- Nohria A, Prsic A, Liu PY, et al. Statins inhibit Rho kinase activity in patients with atherosclerosis. *Atherosclerosis*. 2009;205(2):517-521. doi:10.1016/j.atherosclerosis.2008.12.023
- Zoico E, Rubel S, De Caro A, et al. Brown and Beige Adipose Tissue and Aging. *Front Endocrinol (Lausanne)*. 2019;10:368. Published 2019 Jun 20. doi:10.3389/fendo.2019.00368
- Ikeda K, Maretich P, Kajimura S. The Common and Distinct Features of Brown and Beige Adipocytes. *Trends Endocrinol Metab*. 2018;29(3):191-200. doi:10.1016/j.tem.2018.01.001
- Mancini C, Gohlke S, Garcia-Carrizo F, Zagoriv V, Stephanowitz H, Schulz TJ. Identification of biomarkers of brown adipose tissue aging highlights the role of dysfunctional energy and nucleotide metabolism pathways. *Sci Rep*. 2021;11(1):19928. Published 2021 Oct 7. doi:10.1038/s41598-021-99362-1
- Hara Y, Wakino S, Tanabe Y, et al. Rho and Rho-kinase activity in adipocytes contributes to a vicious cycle in obesity that may involve mechanical stretch. *Sci Signal*. 2011;4(157):ra3. Published 2011 Jan 25. doi:10.1126/scisignal.2001227
- Wei L, Surma M, Yang Y, Tersey S, Shi J. ROCK2 inhibition enhances the thermogenic program in white and brown fat tissue in mice. *FASEB J*. 2020;34(1):474-493. doi:10.1096/fj.201901174RR
- McDonald ME, Li C, Bian H, Smith BD, Layne MD, Farmer SR. Myocardin-related transcription factor A regulates conversion of progenitors to beige adipocytes. *Cell*. 2015;160(1-2):105-118. doi:10.1016/j.cell.2014.12.005
- Yang YM, Kuen DS, Chung Y, Kurose H, Kim SG. Ga12/13 signaling in metabolic diseases. *Exp Mol Med*. 2020 Jun;52(6):896-910. doi:10.1038/s12276-020-0454-5. Epub 2020 Jun 23.
- Barella LF, Jain S, Kimura T, Pydi SP. Metabolic roles of G protein-coupled receptor signaling in obesity and type 2 diabetes. *FEBS J*. 2021;288(8):2622-2644. doi:10.1111/febs.15800
- Yanagida K, Igarashi H, Yasuda D, et al. The Ga12/13-coupled receptor LPA4 limits proper adipose tissue expansion and remodeling in diet-induced obesity. *JCI Insight*. 2018;3(24):e97293. Published 2018 Dec 20. doi:10.1172/jci.insight.97293
- Koo JH, Kim TH, Park SY, et al. Ga13 ablation reprograms myofibers to oxidative phenotype and enhances whole-body metabolism. *J Clin Invest*. 2017;127(10):3845-3860. doi:10.1172/JCI92067
- Kim TH, Yang YM, Han CY, et al. Ga12 ablation exacerbates liver steatosis and obesity by suppressing USP22/SIRT1-regulated mitochondrial respiration. *J Clin Invest*. 2018;128(12):5587-5602. doi:10.1172/JCI97831
- Purushotham A, Schug TT, Xu Q, Surapureddi S, Guo X, Li X. Hepatocyte-specific deletion of SIRT1 alters fatty acid metabolism and results in hepatic steatosis and inflammation. *Cell Metab*. 2009;9(4):327-338. doi:10.1016/j.cmet.2009.02.006
- Kim TH, Koo JH, Heo MJ, et al. Overproduction of inter-α-trypsin inhibitor heavy chain 1 after loss of Ga13 in liver exacerbates systemic insulin resistance in mice. *Sci Transl Med*. 2019;11(513):eaan4735. doi:10.1126/scitranslmed.aan4735
- Noda K, Nakajima S, Godo S, Saito H, Ikeda S, Shimizu T, Enkhjargal B, Fukumoto Y, Tsukita S, Yamada T, Katagiri H, Shimokawa H. Rho-kinase inhibition ameliorates metabolic disorders through activation of AMPK pathway in mice. *PLoS One*. 2014 Nov 3;9(11):e110446. doi:10.1371/journal.pone.0110446
- Landry T, Shookster D, Huang H. Tissue-Specific Approaches Reveal Diverse Metabolic Functions of Rho-Kinase 1. *Front Endocrinol (Lausanne)*. 2021;11:622581. Published 2021 Feb 9. doi:10.3389/fendo.2020.622581

## Rho, Rac, and Cdc42 Activation Assays

Product	Amount	Cat #
<b>RhoA / Rac1 / Cdc42 Activation Assay Combo Biochem Kit</b>	3 x 10 Assays	BK030
<b>RhoA Pull-down Activation Assay Biochem Kit</b>	80 Assays 20 Assays	BK036 BK036-S
<b>RhoA G-LISA™ Activation Assay (Luminescence format)</b>	96 Assays	BK121
<b>RhoA G-LISA™ Activation Assay Kit (Colorimetric format)</b>	96 Assays 24 Assays	BK124 BK124-S
<b>Total RhoA ELISA</b>	96 Assays	BK150
<b>Rac1,2,3 G-LISA™ Activation Assay (Colorimetric format)</b>	96 Assays	BK125
<b>Rac1 G-LISA™ Activation Assay (Luminescence format)</b>	96 Assays	BK126
<b>Rac1 G-LISA™ Activation Assay Kit (Colorimetric Based)</b>	96 Assays 24 Assays	BK128 BK128-S