



Helping advance science,
one protein at a time.

June
2014



this issue

Rho GTPases & Reactive Oxygen Species: Crosstalk & Feedback

Related Publications

Research Tools

Meetings

Muscle & Molecular Motors GRC 2014

July 6-11, West Dover, VT
Sponsor

Intrinsically Disordered Proteins GRC 2014

July 6-11, Easton, MA
Sponsor

Society for Neuroscience 2014

November 15-19,
Washington, D.C.
Booth #1917

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
W: cytoskeleton.com

Distributors

www.cytoskeleton.com/distributors/

Rho GTPases & Reactive Oxygen Species: Crosstalk & Feedback

Redox agents, such as reactive oxygen species (ROS) and reactive nitrogen species (RNS), are key regulators in a variety of signal transduction pathways, including integrin signaling, extracellular matrix adhesion, and inflammation¹⁻³. Rho GTPases are also key regulators of many cellular processes, including cell growth, motility, and adhesion⁴. While redox agents and Rho GTPases operate through a wide array of regulatory mechanisms, crosstalk between ROS/RNS and Rho GTPases is thought to play a pivotal role in many of their physiological functions^{5,6} and in a growing number of pathological processes such as acute lung injury and cancer^{7,8}. This newsletter broadly summarizes some established and emerging concepts in this rapidly developing field (Fig. 1).

Regulation of ROS by Rho GTPases

Major sources of cellular ROS/RNS production are the primary redox-modulating enzymes such as NADPH oxidases (NOX), nitric oxide synthase (NOS), and superoxide dismutase (SOD)^{5,9,10} (Fig. 1). Rho GTPases, in particular Rac1 and 2, are a necessary component of activated NOX complexes and the subsequent generation of ROS (superoxide anion, O₂⁻) from molecular oxygen (O₂)¹¹. Upon activation by extracellular signals, the NOX complex assembles at the cell membrane where Rac (GTP-bound and possibly GDI-bound) is required for ROS generation^{12,13} (Fig. 1). It is not yet clear if Rac acts as an adaptor or mediates electron transfer during O₂⁻ production¹². Rac, specifically Rac1, has also been shown to directly interact with NOS enzymes to regulate the generation of nitric oxide (NO)¹³. The authors found that NOS preferentially binds nucleotide-free GTPase, suggesting the involvement of a guanine nucleotide exchange factor (GEF) and GEF-mediated cellular compartmentalization of Rac-mediated NO generation¹³. Finally, Rac has been shown to interact directly, in a redox-dependent manner, with SOD, an enzyme that catalyzes the conversion of O₂⁻ to hydrogen peroxide (H₂O₂)¹⁴. Under reducing conditions, similar to those in a cell, Rac-GTP binds to SOD, which allows Rac to remain in the active state and propagate the production of O₂⁻ by NOX. Under oxidizing conditions, like those produced locally by O₂⁻ / H₂O₂ generation, SOD was found to dissociate from Rac which was then converted to its inactive GDP form, possibly by a yet-to-be-identified GTPase activating protein (GAP)¹⁴. These reports clearly indicate that Rho GTPases can generate a multitude of reactive oxygen intermediates (RIs) which

can, in turn, result in a diverse array of protein modifications. Which species are generated and the cellular responses elicited will depend upon a number of factors, including the given stimuli, the local cell redox potential, and cellular location.

Direct and Indirect Regulation of Rho GTPases by ROS

Recent evidence points towards the interesting possibility that ROS/RNS can directly regulate Rho GTPases through redox-mediated post-translational modifications (PTMs). For example, Heo and Campbell¹⁵ demonstrated that the conserved redox

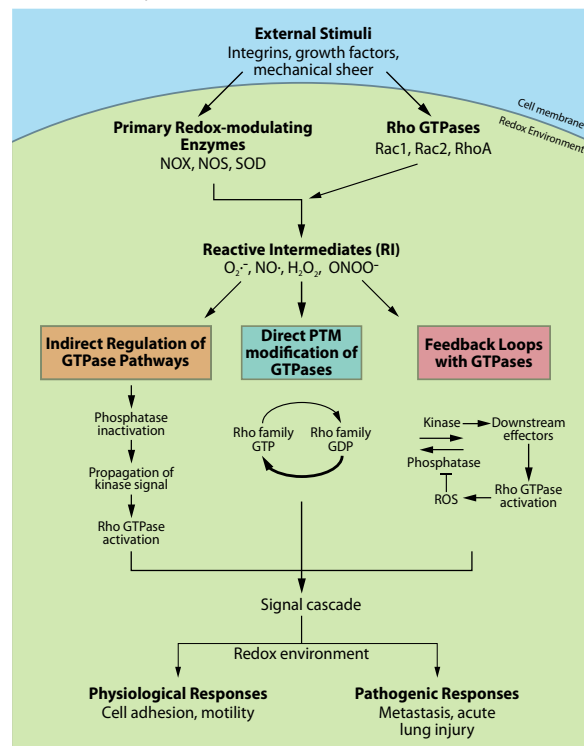


Figure 1: Schematic Representation of Crosstalk Between Rho GTPases and ROS in Signal Transduction. The diagram is a schematic representation of possible crosstalk between Rho GTPases and ROS/RNS in a stylized cell. Reactions given under indirect (orange), direct (blue) and feedback (pink) sections are representations of the most commonly reported mechanisms that fall under these headings. For specific pathways, the reader is directed to the references and examples given in this newsletter.



Continued from Page 1

sensitive motif Cys¹⁸ (Rac numbering) located at the end of the conserved sequence (GXXXGK(S/T)), found in the p-loop of several Rho GTPases, including RhoA, Rac, and Cdc42, is the site of redox-mediated reversible cysteine oxidation *in vitro*. This oxidation results in nucleotide displacement and, under reducing conditions, GTP nucleotide exchange and GTPase activation¹⁵. This mechanism was later verified *in vivo* when it was shown that H₂O₂-mediated activation of RhoA and stress fiber formation is abolished by mutation of the cysteines Cys¹⁶ and Cys²⁰ (Rho numbering)¹⁶. Importantly, the cysteine mutations do not inhibit the ability of RhoA to be activated by GEFs¹⁶. The authors suggested that GTPase activation likely occurs via two parallel mechanisms, classical enzymatic GEFs/GAPs and a redox-mediated exchange (Fig. 1). They postulate that the cells' redox potential may help determine which mechanism is operational¹⁶. Further evidence for a redox-driven exchange mechanism has come from the finding that RhoA is nitrated at Tyr³⁴, within the switch one region of the protein, via enhanced NO production due to acute lung injury⁷. The authors went on to demonstrate that this PTM produced a GEF-like activation of RhoA. Finally, redox-mediated PTMs have also been proposed to inactivate Rho GTPases¹⁷.

In addition to direct regulation, indirect redox-dependent regulation of Rho GTPases involving phosphatases has been reported. For example, Rac-activated ROS production leads to inactivation of RhoA via a mechanism involving redox-mediated inactivation of the low molecular weight protein tyrosine phosphatase (LMW-PTP), subsequent elevation of p190 RhoGAP activity, and deactivation of RhoA¹⁸. Interestingly, inactivation of phosphatases by ROS appears to be a common mechanism for ROS regulation of diverse cell pathways¹⁹ (Fig. 1).

Feedback Loops between ROS and Rho GTPases

The hypothesis that Rho GTPases participate in redox feedback loops⁵ has recently been demonstrated for regulation of leukocyte chemotaxis²⁰ (Fig. 1). In a familiar theme, Rac mediates the production of ROS that results in a redox-mediated inactivation of a phosphatase. This in turn allows upregulation of a kinase that mediates GEF activation of Rac²⁰ (Fig. 1). A feedback mechanism has also been proposed for SOD regulation of Rac-mediated generation of ROS¹⁴.

Summary

In conclusion, the relationship between Rho GTPases and redox agents is highly complex. Interesting areas for future work might examine the relationship between GAP/GEF-regulated and redox-mediated exchange processes, addressing whether they work in parallel or synergistically, and would they make good druggable targets.

G-Switch™ Activators and Inhibitors

Product	Cat. #	Amount
Rho inhibitor I: ADP ribosylation of Rho Asn-41	CT04-A	1 x 20µg
	CT04-B	5 x 20µg
Rho Activator II: Deamidation of Rho Gln-63	CN03-A	3 x 20µg
	CN03-B	9 x 20µg
Rho/Rac/Cdc42 Activator I	CN04-A	3 x 20µg
	CN04-B	9 x 20µg
Rho Pathway Inhibitor I: Rho kinase (ROCK) inhibitor Y-27632	CN06-A	5 x 10 units
	CN06-B	20 x 10 units

References

1. Yan B. & Smith J.W. 2000. A redox site involved in integrin activation. *J. Biol. Chem.* **275**, 39964-39972.
2. Chiarugi P. et al. 2003. Reactive oxygen species as essential mediators of cell adhesion: the oxidative inhibition of a FAK tyrosine phosphatase is required for cell adhesion. *J. Cell Biol.* **161**, 933-944.
3. Leto T.L. & Geiszt M. 2006. Role of Nox family NADPH oxidases in host defense. *Antioxid. Redox Signal.* **8**, 1549-1561.
4. Jaffe A.B. & Hall A. 2005. Rho GTPases: biochemistry and biology. *Annu. Rev. Cell Dev. Biol.* **21**, 247-269.
5. Mitchell L. et al. 2013. Redox regulation of Ras and Rho GTPases: mechanism and function. *Antioxid. Redox Signal.* **18**, 250-258.
6. Ferro E. et al. 2014. Ras GTPases are both regulators and effectors of redox agents. *Ras Signaling: Methods Mol. Biol.* **1120**, 55-74.
7. Rafikov R. et al. 2014. Lipopolysaccharide-induced lung injury involves the nitration-mediated activation of RhoA. *J. Biol. Chem.* **289**, 4710-4722.
8. Myant K.B. et al. 2013. ROS production and NF-kappaB activation triggered by Rac1 facilitate WNT-driven intestinal stem cell proliferation and colorectal cancer initiation. *Cell Stem Cell.* **12**, 761-773.
9. Ueyama T. et al. 2006. Involvement of Rac1 in activation of multicomponent Nox-1 and Nox-3 based NADPH oxidases. *Mol. Cell Biol.* **26**, 2160-2174.
10. Werner E. 2004. GTPases and reactive oxygen species: switches for killing and signaling. *J. Cell Sci.* **117**, 143-153.
11. Hordijk P.L. 2006. Regulation of NADPH oxidases: the role of Rac proteins. *Circ. Res.* **98**, 453-462.
12. Bokoch G.M. 1995. Regulation of the phagocyte respiratory burst by small GTP-binding proteins. *Trends Cell Biol.* **5**, 109-113.
13. Di-Poi N. et al. 2001. Mechanism of NADPH oxidase activation by Rac/Rho-GDI complex. *Biochemistry.* **40**, 10014-10022.
14. Harraz M.M. et al. 2008. SOD1 mutations disrupt redox-sensitive Rac regulation of NADPH oxidase in a familial ALS model. *J. Clin. Invest.* **118**, 659-670.
15. Heo J. and Campbell S.L. 2005. Mechanism of redox-mediated guanine nucleotide exchange on redox-active Rho GTPases. *J. Biol. Chem.* **280**, 31003-31010.
16. Aghajanian A. et al. 2009. Direct activation of RhoA by reactive oxygen species requires a redox-sensitive motif. *PLoS ONE.* **4**, e8045.
17. Zuckerbraun B.S. et al. 2007. Nitric oxide-induced inhibition of smooth muscle cell proliferation involves S-nitrosation and inactivation of RhoA. *Am. J. Physiol. Cell Physiol.* **292**, C824-C831.
18. Nimnual A.S. et al. 2003. Redox-dependent downregulation of Rho by Rac. *Nat. Cell Biol.* **5**, 236-241.
19. Meng T.C. et al. 2002. Reversible oxidation and inactivation of protein tyrosine phosphatases *in vivo*. *Mol. Cell.* **9**, 387-399.
20. Kuiper J.W. et al. 2011. Rac regulates PtdInsP3 signaling and the chemotactic compass through a redox-mediated feedback loop. *Blood.* **118**, 6164-6171.

Small G-Protein Activation Assay Kits

Product	Cat. #	Amount
RhoA Activation Assay Biochem Kit (bead pull down format)	BK036	80 assays
Rac1 Activation Assay Biochem Kit (bead pull down format)	BK035	50 assays
Cdc42 Activation Assay Biochem Kit (bead pull down format)	BK034	50 assays
G-LISA™ RhoA Activation Assay Biochem Kit (colorimetric format)	BK124	96 assays
G-LISA™ Rac 1 Activation Assay Biochem Kit (colorimetric format)	BK128	96 assays
G-LISA Cdc42 Activation Assay Biochem Kit (colorimetric format)	BK127	96 assays