



Helping advance science,
one protein at a time.

March 2021

Small G-proteins Related Publications Research Tools

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

Recent Citation

Spotlights

SARS-CoV-2 infection alters the transcriptomic landscape of human airway cells

A Live Look At Membrane Tension At The Plasma Membrane And Beyond

RHOA-mediated mechanical force generation through Dectin-1

Small G Proteins and Microvesicle-Mediated Cancer Biogenesis

Under control of guanine nucleotide exchange factors (GEFs), small G-proteins cycle between the inactive GDP-bound state and the active GTP-bound state; thus, acting as molecular switches in a diverse multitude of cellular processes¹. Small G-proteins have been implicated as central modulators of biological processes including cell growth, cell differentiation, and cell movement^{2,3}. It is now widely recognized that aberrant activity of small G-proteins belonging to the Ras superfamily, which were initially characterized as tumor oncoproteins, contribute heavily to the role of human diseases such as cancer⁴. Their importance is so integral that scientists continue to link small G-proteins to a wide range of cellular processes involved in tumor progression and metastasis.

Microvesicles (MVs), a type of extracellular vesicle (EV) shed from the outward blebbing of cellular plasma membranes with a diameter of 100-1000 nm, play a role in oncogenesis⁵. Although the underlying mechanisms of MV generation and release are not completely understood, the impact of microvesicular cargo is well known. MVs alter cellular biology through delivery of their diverse cargo that is known to include proteins, lipids, microRNAs (miRNAs), messenger RNAs (mRNAs), and long non-coding RNAs (lncRNAs)^{6,7}. MVs are known to participate in several critical events in cancer biogenesis such as tumor pathogenesis and metastasis⁸, which seems quite logical given that EVs in circulation can influence distant tissues⁹. One interesting development has been the discovery that small G-protein activation states tune EV content and secretion, and may be another mechanism by which small G-proteins influence cancer progression.

Cdc42 and EGFR regulate microvesicle shedding in MDA-MB-231 breast cancer cells

It has previously been shown that GTP-bound Cdc42 binds Ras GTPase-activating-like protein 1 (IQGAP1)¹⁰. A recent study by Wang et al. further revealed that RNAi-mediated inhibition of IQGAP1 in MDA-MB-231 triple-negative breast cancer cell line significantly reduced MV shedding concomitant with reduction of MV biogenesis biomarkers VEGF90k and flotillin-2¹¹. In this study, counts of MV release were quantified with fluorescent dyes and revealed that cells expressing plasmids for constitutively active or inhibited Cdc42 (Cdc42Q61L and Cdc42T17N, respectively) were not significantly different from controls; conversely, fast GTP-GDP cycling-expressing cells (Cdc42F28L) exhibited enhanced MV release and biomarker expression. Furthermore, Cdc42Q61L-expressing cells capable of binding with IQGAP exhibited MV shedding, while binding-incapable Cdc42T17N cells displayed reduced shedding ability. Cdc42's control of MV shedding was linked

to epidermal growth factor receptor (EGFR) regulation of cell surface protein internalization of VEGF90k, and this process was reversed with EGFR-inhibitor AG1478. These results suggest that Cdc42 complexed with IQGAP signaling is required for MV biogenesis, and Cdc42-mediated stimulation of EGFR facilitates MV biogenesis.

ARF6 and Exportin-5 regulate EV trafficking in a GEF-dependent manner

miRNAs that target tumor suppressor genes contribute to the development of tumor microenvironments; therefore, understanding how miRNAs are selected for packaging into microvesicles may be critically important⁵. A study by Clancy et al. revealed a link between ARF6 activation and the packaging of pre-miRNA and other miRNA-related cargo into tumor derived microvesicles¹². Cells expressing constitutively active ARF6-Q67L, which exhibit a 5-fold increase in ARF6, shed 3-fold greater MVs than controls, and miRNA sequencing of TMVs reveals a global increase of TMV miRNA cargo. Increased TMV miRNA content

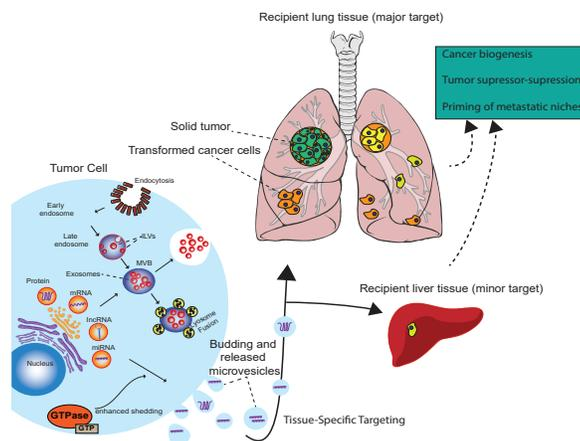


Figure 1. Diagram of microvesicle trafficking to tissues and outcomes.

led investigators to suspect, and through interrogation of public databases, implicate pre-miRNA transporter Exportin-5 as an ARF6 binding partner. ARF6-Exportin-5 binding was confirmed through protein-protein interaction mapping with mass spectrometry and immunoprecipitation. ARF6-Exportin-5 binding was reversed in cells through SecinH3 treatments, which broadly inhibits cytohesins (ARF6-GEFs), resulting in the loss of Exportin-5 from TMVs. Immunofluorescence analysis also confirmed ARF6-Exportin-5 co-localization in budding TMVs with



The Protein
Experts

Helping advance science,
one protein at a time.

Small G-Protein PRODUCTS

Continued from Page 1

continued Exportin-5 detection within isolated tumor derived microvesicles. Investigators utilized public databases to deduce cytohesin 3 (GRP1) as the most likely binding partner for the ARF6-GTP-Exportin-5 complex. siRNA-mediated knockdown confirmed that GRP1 acts in concert with ARF6 and Exportin-5 to traffic pre-miRNA cargo into nascent TMVs. This study highlights the importance of the ARF6-Exportin-5 complex as a regulator of miRNA packaging into TMVs.

RalA and RalB control tumor volume and package EVs for tissue-specific metastatic niche formation

Ghoroghi et al. sought to further clarify how RalA and RalB GTPases can tune both the content of TMVs and rate of secretion. Interestingly, it was shown that shRNA-mediated knockdown of RalA or RalB, could significantly reduce the amount of EVs secreted in both human cancer cells lines and the nematode *C. elegans*; illustrating that EV secretion under control of GTPases is evolutionarily conserved⁵. Confocal quantification confirmed a concomitant and significant reduction in the number of multivesicular bodies (MVBs), which contain exosomes prior to release; thereby, suggesting that MVBs are under the control of RalA and RalB expression. To determine how GTPases can influence MVB homeostasis investigators focused on phospholipase Ds (PLDs) which are known to impact cancer progression, exosome secretion, and are targets of RalA and RalB. Suppression of PLD1 through chemical inhibitor CAY10594 in PLD expressing 4T1 cells led to a recapitulation of the suppressed EV secretion and MVB counts seen by RalA/B knockdown. Automated examination of RalA/B protein levels by immunohistochemistry in breast cancer patient-tumors revealed significant overexpression. Xenografting small hairpin RNA (shRNA)-mediated RalA/B-depleted cells into Balb/c mice mammary ducts revealed that while shRalA depletion significantly increased tumor volume, shRalB depletion resulted in tumors with smaller volume than shControl. Intriguingly, although RalA/B depletion increased tumor volume it appeared to reduce metastasis, with shControl-grafted mice exhibiting the highest metastatic foci in serially-sectioned lung tissue.

Functional assays determined that EV content had vascular permeability-enhancing properties that disrupted endothelial cell adherent and tight junctions that could be reversed by small hairpin depletion. EVs isolated from control 4T1 cells were serially injected intra-orbitally into mice to prime metastatic niches and resulted in predominant localization in the lungs that was not observed when EVs were isolated from shRalA/B cells indicating that RalA/B prime EVs to target lung tissues. Interrogation of the EV content with mass spectroscopy revealed differentially expressed proteins found in RalA/B EVs but not shRalA/B EVs, including CD146 and MCAM. Treatment of RalA/B isolated EVs with anti-CD146 and anti-MCAM antibodies abolished lung targeting and pre-metastatic niche formation suggesting that RalB-mediated tuning predisposes EVs to enhance metastatic niche formation specifically in lung tissues.

Summary

Altogether these data present an exciting new perspective for small G-protein research. Perturbation of GTPases belonging to the RAS superfamily can modulate the content and quantity of cancer-propagating EVs. Importantly, these works also identified small G-protein accomplices, implicating GEFs and scaffolding proteins, required for microvesicle packaging complexes or downstream effectors which may be important for the development of future druggable targets and improving our understanding of cancer biogenesis. Cytoskeleton has an array of small G-protein tools and membrane probes to aide investigators as they work to define the critical role that small GTPases play in regulating TMVs.

References

1. Nestler, E. J. & Duman, R. S. Small G Proteins. (1999).
2. Bos, J. L., Rehmann, H. & Wittinghofer, A. GEFs and GAPs: Critical Elements in the Control of Small G Proteins. *Cell* vol. 129 865–877 (2007).
3. Csépanyi-Kömi, R., Lévy, M. & Ligeti, E. Small G proteins and their regulators in cellular signalling. *Molecular and Cellular Endocrinology* vol. 353 10–20 (2012).
4. Colicelli, J. Human RAS superfamily proteins and related GTPases. *Science's*

5. Ghoroghi, S. et al. Ral GTPases promote breast cancer metastasis by controlling biogenesis and organ targeting of exosomes. *Elife* 10, 1–29 (2021).
6. Kelemen, E., Danis, J., Göblös, A., Bata-Csörgő, Z. & Széll, M. Exosomal long non-coding RNAs as biomarkers in human diseases. *Electron. J. Int. Fed. Clin. Chem. Lab. Med.* 30, 224–236 (2019).
7. D'Asti, E. et al. Oncogenic extracellular vesicles in brain tumor progression. *Front. Physiol.* 3 JUL, 1–15 (2012).
8. Wang, T. et al. Hypoxia-inducible factors and RAB22A mediate formation of microvesicles that stimulate breast cancer invasion and metastasis. *Proc. Natl. Acad. Sci. U. S. A.* 111, (2014).
9. D'Souza-Schorey Crislyn, C. & Clancy, J. W. Tumor-derived microvesicles: Shedding light on novel microenvironment modulators and prospective cancer biomarkers. *Genes Dev.* 26, 1287–1299 (2012).
10. Ratajczak, M. Z. & Ratajczak, J. Extracellular microvesicles/exosomes: discovery, disbelief, acceptance, and the future? *Leukemia* vol. 34 3126–3135 (2020).
11. Wang, J. et al. Cdc42 functions as a regulatory node for tumour-derived microvesicle biogenesis. *J. Extracell. Vesicles* 10, (2021).
12. Clancy, J. W., Zhang, Y., Sheehan, C. & D'Souza-Schorey, C. An ARF6–Exportin-5 axis delivers pre-miRNA cargo to tumour microvesicles. *Nat. Cell Biol.* 21, 856–866 (2019).

Bead pull-down Activation Assays

Product	Assays	Cat. #
Arf1 Activation Assay Biochem Kit	20	BK032-S
Arf6 Activation Assay Biochem Kit	20	BK033-S
GGA3-PBD Beads (Arf1 + Arf6)	20-25	GGA07
Cdc42 Activation Assay Biochem Kit	50	BK034
Rac + Cdc42 Activation Assay Biochem Kit	50	PAK02
Rac1 Activation Assay Biochem Kit	50	BK035
Ras Activation Assay Biochem Kit	50	BK008
Rhotekin RBD beads (RhoA/B/C)	30	RT02
RhoA Activation Assay Biochem Kit	80	BK036
RhoA/Rac1/Cdc42 Activation Assay Combo Biochem Kit	3 x 10	BK030
RalA Activation Assay Biochem Kit	50	BK040

G-LISA Activation Assay Kits

Product	Assays	Cat. #
Arf1 G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK132
Arf6 G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK133
Cdc42 G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK127
Rac1 G-LISA™ Activation Assay Kit (Luminescence format)	96	BK126
Rac1 G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK128
RalA G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK129
RhoA G-LISA™ Activation Assay (Luminescence format)	96	BK121
RhoA G-LISA™ Activation Assay Kit (Colorimetric format)	96	BK124
RhoA / Rac1/ Cdc42 G-LISA™ Activation Assay Bundle 3 kits (24 assays per kit)	96	BK135
Rac1,2,3 G-LISA™ Activation Assay (Colorimetric format)	96	BK125
Rac1 G-LISA™ Activation Assay (Luminescence format)	96	BK126
Rac1 G-LISA™ Activation Assay Kit (Colorimetric Based)	96	BK128
Ras G-LISA™ Activation Assay Kit (Colorimetric Based)	96	BK131
Total RhoA ELISA	96	BK150