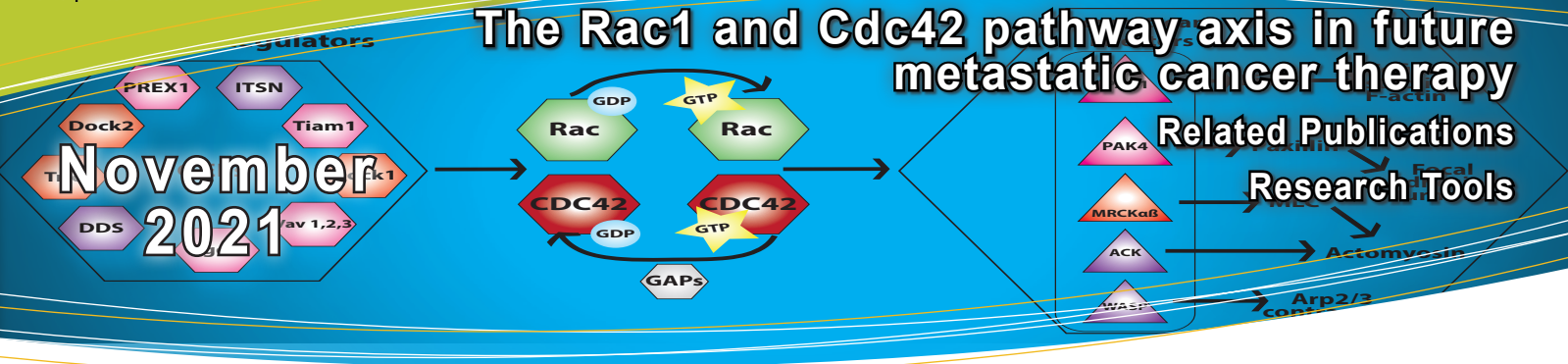




Helping advance science,  
one protein at a time.



**Sponsored Conferences**

Society for Neuroscience 2021  
Virtual Meeting  
November 8th-11th 2021

ASCB/EMBO Cell Bio 2021  
Virtual Meeting  
December 1st -10th, 2021

ASN: American Society for Neurochemistry  
Roanoke, VA  
April 10th - 14th, 2022

Journal of Cell Science Meeting 2022  
Wotton House, Surrey, UK  
May 8th-11th 2022

**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

### Rac1 and Cdc42 pathways as pivotal axis in future metastatic cancer therapy

The appearance of Rac and Cdc42 proteins on the radar of the medical scientific community started in the 1990's with seminal papers by Professor Alan Hall's laboratory at MRC, UK. They identified Rac and Cdc42 small G-proteins (R&C) as central regulators of cell motility, shape and protrusions (lamellipodia and filopodia)<sup>1,2</sup>. Now several decades later, the concept has been fully proven in a multitude of publications (see refs. in these reviews 3 & 4), and the realization that these proteins are central to cancer cell motility and invasion has led the field to explore them as candidates for therapeutic intervention. Targeting metastatic cancer has become mainstream due to an average of 58% of cancer-related deaths, 10-90% depending on cancer type, are due to consumption by metastasis<sup>5,6</sup>. It must be noted, therapies based on the R&C axis are cell stasis interventions rather than the more typical cytotoxic cancer therapies such as cisplatin and paclitaxel; however, these different mechanisms could be more interdependent than previously thought.

of metastatic breast cancer<sup>10</sup>. These effects are highly dominant, but when considering other genes, which also dominate a metastatic phenotype, the frequency of their occurrence is very low i.e. mutations occur at 0-1%, and altered levels at 0.5-4.0%<sup>11,12</sup>. On the other hand, these proteins are ubiquitous in all mammalian cell types, so therapeutic intervention has to circumvent the normal cell's functional requirements. As a result, most of the early R&C direct binding drug candidates have faltered due to toxic side effects (reviewed in 13). In an effort to find a good therapeutic window, the field is turning its focus to the upstream and downstream regulators of the R&C axis as shown in Fig1.

### The Rac & Cdc42 central axis

As shown in Figure 1, the Rho family small G-proteins hold a central regulator focal point between upstream receptor signaling pathways and down stream effector pathways. In some instances, this centrality is highlighted with mutated isoforms or altered levels of protein which directly link to cancer development and prognosis. For example, the human Rac1 mutation P29S/L is present in melanomas<sup>7</sup> where haptotaxis (directionality) and velocity of cells are reduced<sup>8</sup>, and also Programmed Cell Death Ligand (PD-L1) is up-regulated<sup>9</sup>. In contrast, elevated levels of Cdc42 expression in mammary gland epithelial cells increased motility and invasion and poor prognosis

### Upstream Activators

Upstream activators are dominated by GTP Exchange Factors (GEFs), which number almost 100 different genes in the human genome. GEFs are activated by cell surface receptors e.g. GPCRs, RTKs, integrins. In turn, GEFs activate small G-proteins by increasing GTP/GDP exchange rate at the nucleotide-binding site. e.g. Tiam, Vav1 and Vav2 are GEFs of Rac, and Dbl, Dbp, and ITSN are GEFs of Cdc42 and some, Tiam1, Vav1,2 activate both GTPases (Fig.1). Due to Mg<sup>2+</sup>:GTP having a higher binding affinity than GDP, and a ten-fold higher level than GDP in cells, the exchange of nucleotides favors GTP in a homeostatic cancer cell. GEFs can be upregulated by mutations that increase binding affinity to the small G-protein, or mutations that simulate its phosphorylated form, or simply increased expression via demethylation of their promoter regions (reviewed in 14).

Only in the last 10 years, has the prevalence of GEF oncogenes in metastatic cancer been found to be broad

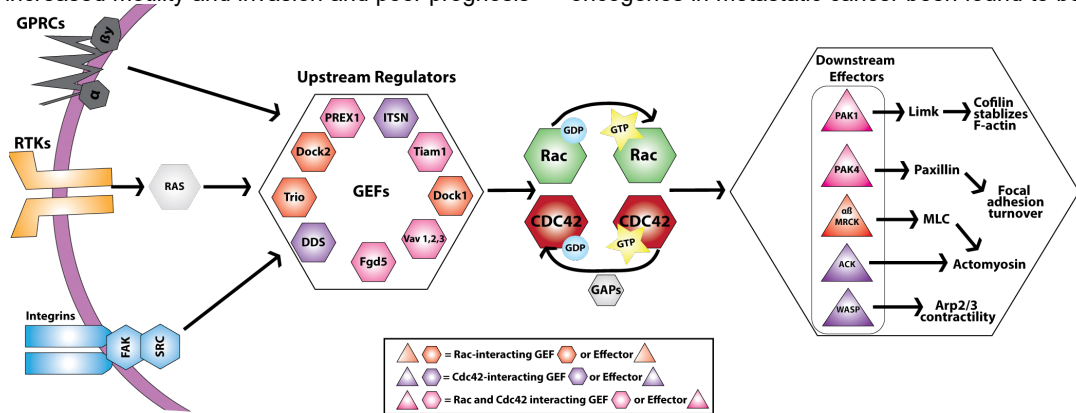


Figure legend: Upstream and downstream pathways converging on the Rac and Cdc42 axis (adapted from refs. 14 & 17).



Helping advance science,  
one protein at a time.

# Rac1 and Cdc42 Detection PRODUCTS

## Continued from Page 1

due to the advent of genomic databases<sup>15</sup>. Although GEFs are broadly represented as oncogenes, individually mutated GEF genes make up a small portion of the total pool of players. For example, in melanoma metastasis, Tiam GEF mutations have a 9.0% prevalence while Vav2 mutations have a 3.6% prevalence, whereas all mutated RC GEFs are present in 80.7% of melanomas<sup>14</sup>. Under these circumstances, many GEF targets are necessary for a comprehensive drug development effort, and tumor screening is necessary to delineate causative GEF genes in individual patients. There are many options for drug development in R&C GEFs, for example small molecules could bind the protein interaction site, or an allosteric site, or mask the phosphorylation site thus inhibiting its activity<sup>16</sup>.

### Downstream Activators

Fewer players are currently known on the downstream side of the R&C axis. Most are kinases that are activated by GTP-bound forms of R&C. In particular, p21-activated kinase (PAK), activated Cdc42-associated kinase (ACK), and more recently myotonic dystrophy-related Cdc42 binding kinase (MRCK) have been the focus of drug development efforts<sup>17</sup>. For example, Unbekandt et al. (2018) present a tour de force of MRCK inhibitor development for melanoma<sup>18</sup>.

As a side note, many of the oncogenic GEFs' promoters are demethylated during transition to metastasis<sup>14</sup>, which could conceivably be targeted with demethylase inhibitors, but the question of specificity comes into the picture when using those enzymes with a broad, sometimes undefined, substrate range.

### Interdependence of the Ras and R&C axes in metastatic cancer

Critically, it has been reported that primary cancer oncogenes, such as K-Ras and N-Ras, work in concert with the R&C pathway when transitioning from primary to metastatic phenotype<sup>19,20</sup>, which makes primary cancer progression to metastasis highly dependent on R&C axis activation/dysfunction. Some GEFs, e.g. P-Rex, have been shown to contribute at a later stage metastatic cancer progression<sup>21,22</sup>, which opens the possibility to develop later stage-specific therapeutics. The multitude of RC axis modulators make it essential to develop diagnostics to determine which GEFs and effectors are differentially expressed or mutated, this information will allow oncologists to pick the matched drug for a particular patient.

Although metastatic cancer is destined to kill on average 58% of cancer patients<sup>5,6</sup>, this realization is not equated to the NIH & FDA's focus to encourage therapeutics in this area. In fact, only 16% of active cancer clinical trials are for metastatic disease<sup>23</sup>. Clinical trials targeting prevention and treatment of metastasis are a difficult sell when evidence of metastasis is not apparent early in the disease. And hence, its imperative for scientists and clinical oncologists to communicate this aspect at every opportunity so the barriers are decreased as early as possible. Cytoskeleton is proud to provide highly dependable and accurate Pull-down and GLISA Activation Assay Kits which have been used by researchers in this field over the past 20 years.

## Bead pull-down Activation Assays

Product	Assays	Cat. #
Cdc42 Activation Assay Biochem Kit	50	BK034
Rac1 Activation Assay Biochem Kit	50	BK035
Ras Activation Assay Biochem Kit	50	BK008
RhoA Activation Assay Biochem Kit	80	BK036
RhoA/Rac1/Cdc42 Activation Assay Combo Biochem Kit	3 x 10	BK030

## References

- Ridley AJ, & 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, p389-399.
- Ridley AJ, et al. 1992. The small GTP-binding protein rac regulates growth factor-induced membrane ruffling. *Cell*, 70, p401-410.
- Murali A, and Rajalingam K. 2014. Small Rho GTPases in the control of cell shape and mobility. *Cell Mol Life Sci*,71(9):1703-21. doi: 10.1007/s00018-013-1519-6. Epub 2013 Nov 26.
- Warner H, Wilson BJ and Caswell PT. 2019. Control of adhesion and protrusion in cell migration by Rho GTPases. *Curr. Op. Cell Biol*. 56, p64-70.
- Chaffer C, and Weinberg RA. 2011. A perspective on cancer cell metastasis. *Science*, 331, p1559-64.
- Dillekás H, Rogers MS, Straume O. 2019. Are 90% of deaths from cancer caused by metastases? *Cancer Medicine*, 8, p5574–5576.
- Mar VJ et al. 2014. Clinical and pathological associations of the Rac1 P29S mutations in primary cutaneous melanoma. *Pigment Cell Melanoma Res.*, 27, p1117-1125.
- King SJ, et al. 2016. Lamellipodia are crucial for haptotactic sensing and response. *J. Cell Sci.* 129 p2329-2342.
- Vu HL, et al. 2015. Rac1 P29S regulates PD-L1 expression in melanoma. *Pigment Cell melanoma Res.* 28, p590-598.
- Bray K, et al. 2013. Cdc42 overexpression induces hyperbranching in the developing mammary gland by enhancing cell migration. *Breast Cancer Research*, 15, R91.
- Svensmark JH & Brakebusch C. 2019. Rho GTPases in cancer: friend or foe? *Oncogene*, 38, 7447-7456.
- Aspenstrom P. 2018. Activated Rho GTPases in cancer - The beginning of a new paradigm. *Intl. J. Mol. Sci.* 19, p3949-63.
- Maldonado MM, & Dharmawardhane S. 2018. Targeting Rac and Cdc42 GTPases in Cancer. *Cancer Res.* 78 (12), p3101–3111.
- Maldonado MM, et al. 2020. Targeting Rac and Cdc42 GEFs in metastatic cancer. *Front. Cell Dev. Biol.* 8, Article 201.
- Gao et al. 2013. Integrative analysis of complex cancer genomics and clinical profiles using the cBioPortal. *Sci. Signal.* 6, Article 269.
- Murphy N. et al. 2021. Progress in the therapeutic inhibition of Cdc42 signalling. *Biochem. Soc. Trans.* 49, p 1443-1456.
- Clayton NS, & Ridley AJ. 2020. Targeting Rho GTPase signalling networks in cancer. *Front. Cell Dev. Biol.* 8, Article 222.
- Unbekandt M. et al. 2018. Discovery of potent and selective MRCK inhibitors with therapeutic effect on skin cancer. *Cancer Res.*, 78(8), p2096-2114.
- Cooke M. et al 2020. Rac-GEF/Rac Signaling and Metastatic Dissemination in Lung Cancer. *Front. Cell Dev. Biol.* 8, Article 118.
- Kissil JL. 2007. Requirement for Rac1 in a K-ras induced lung cancer in the mouse. *Cancer Res.* 67(17), 8089-94.
- Cardama et al. 2014(a). Preclinical Development of Novel Rac1-GEF Signaling Inhibitors using a Rational Design Approach in Highly Aggressive Breast Cancer Cell Lines. *Anticancer Agents Med Chem.* 14(6), p840–851.
- Cardama et al. 2014(b). Proapoptotic and antiinvasive activity of Rac1 small molecule inhibitors on malignant glioma cells. *Onc. Targets Ther.* 7, p2021–2033.
- Clinicaltrials.gov as of 2021-09-16, and search for "cancer" and limit to "active, not recruiting". Then calculate percent of titles containing the word "metastatic". Result: 1271 active, not recruiting, cancer trials, wherein 203 contained the word "metastatic" in the title.

## G-LISA Activation Assay Kits

Product	Assays	Cat. #
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