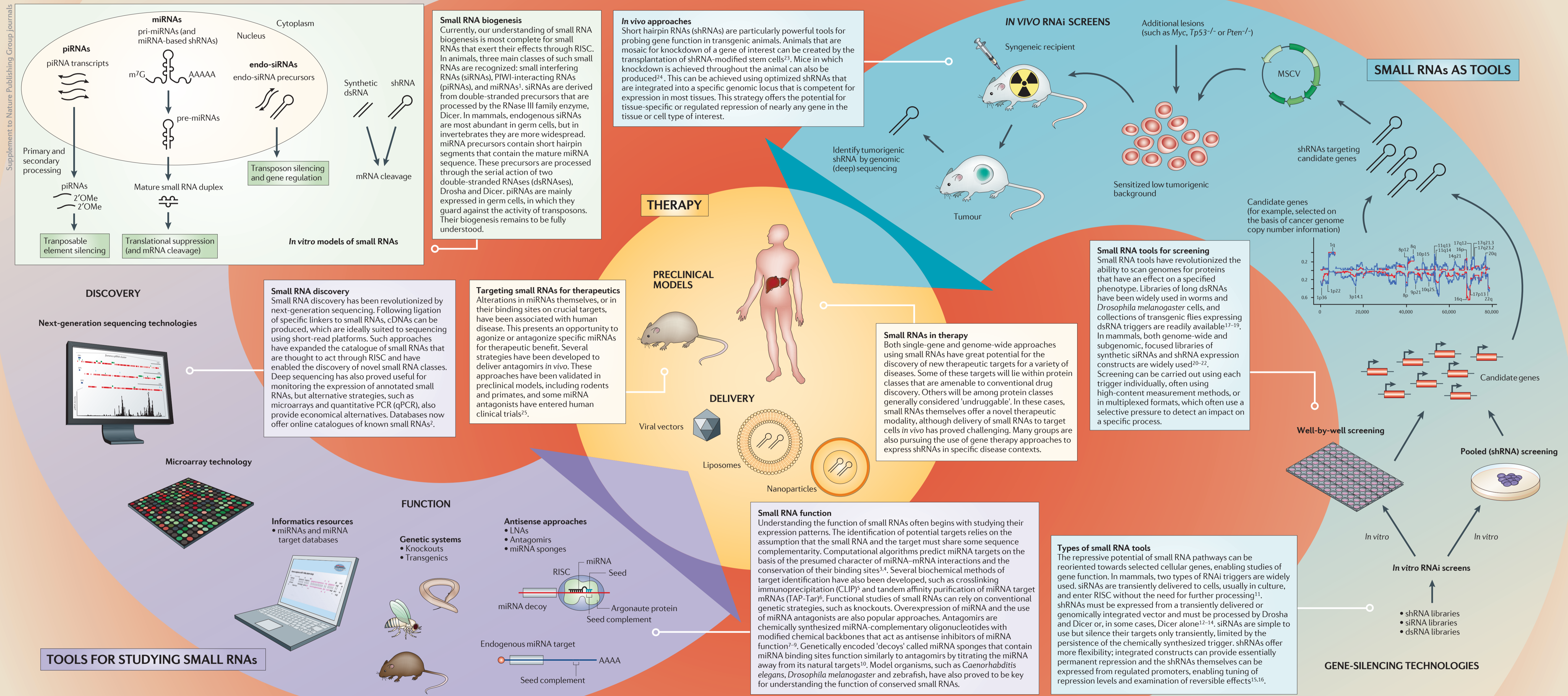


During the past decade, small RNAs have emerged as crucial regulators of gene expression and genome function, having roles in almost every aspect of biology¹. Many small RNAs act through RNA interference (RNAi)-related mechanisms, which involve programming the RNA-induced silencing complex (RISC) to recognize and repress targets. One class of small RNA, the microRNAs (miRNAs), naturally regulates programmes of gene expression. Altered miRNA function contributes to human disease, and manipulation of specific miRNAs is now being pursued as a novel therapeutic modality. Small RNAs have

also been adapted for use as tools based on reprogramming the RNAi machinery to silence specific coding or non-coding RNAs. These tools have been exploited to investigate gene function in cultured cells and in living animals. Genome-scale collections of silencing triggers permit phenotype-based genetic screens to be carried out easily in organisms in which they were previously difficult or impossible. Such strategies are being used to discover and validate new therapeutic targets, and small RNAs themselves may offer a mechanism for inhibiting targets that are currently viewed as 'undruggable'.



Horizon Discovery RNAi Solutions from Dharmacon™
Access the largest and most complete portfolio of innovative and technologically advanced RNAi tools for gene silencing. Since discovery of the endogenous RNAi pathway, our scientists have made key contributions to the field of RNAi that have been foundational to the development of potent, specific and highly functional siRNA, microRNA-adapted shRNA and microRNA reagents. Today, Horizon's RNAi products from Dharmacon facilitate a variety of applications including transient, long-term, inducible and *in vivo* RNAi strategies that extend from basic research in biology to improved therapeutic strategies in medicine.

- siRNA — Patented dual-strand modifications in ON-TARGETplus™ siRNA provides unrivaled specificity and potency; Accell™ siRNA is the only siRNA reagent that can be delivered to difficult-to-transfect cells WITHOUT a transfection reagent
- SMARTvector™ shRNA — microRNA-adapted shRNA for constitutive and inducible RNAi provide specific and potent, long-term silencing
- miRIDIAN™ microRNA — Up-regulate or suppress endogenous mature microRNA function with rationally-designed synthetic and expressed microRNA mimic and hairpin inhibitors

In addition, our portfolio of molecular biology tools includes large collections of cDNAs and ORFs for gene overexpression and RNAi rescue, PCR and qPCR reagents, delivery reagents as well as validated protocols to support the entire RNAi workflow. Horizon Discovery continues to be the industry leader in the field of RNA chemistry, RNAi biology and high-throughput screening, and partners with the RNAi screening community through technical support and screening services. From single gene knockdown to genome-scale RNAi screens, find your RNAi solutions at <https://horizondiscovery.com/en/navigation/gene-modulation/knockdown>

Abbreviations
m⁷G, 7-methylguanosine cap; 2' OMe, 2'-O-methyl; MSCV, murine stem cell virus.
Affiliations
Kenneth Chang and Gregory J. Hannon are at Cold Spring Harbor Laboratory, Watson School of Biological Sciences, 1 Bungtown Road, Cold Spring Harbor, New York 11724, USA.
The authors declare competing financial interests: G.J.H. is a consultant for GE Healthcare Dharmacon, Inc. K.C. declares no competing interests.

Acknowledgements
K.C. and G.J.H. are supported by grants from the US National Institutes of Health, the US Department of Defense Breast Cancer Research Program and a kind gift from K. W. Davis.
Edited by Louisa Flintoft; copy-edited by Matthew Smyllie; designed by Claudia Bentley.
© 2020 Nature Publishing Group.
<http://www.nature.com/nrg/posters/>
References are available online.