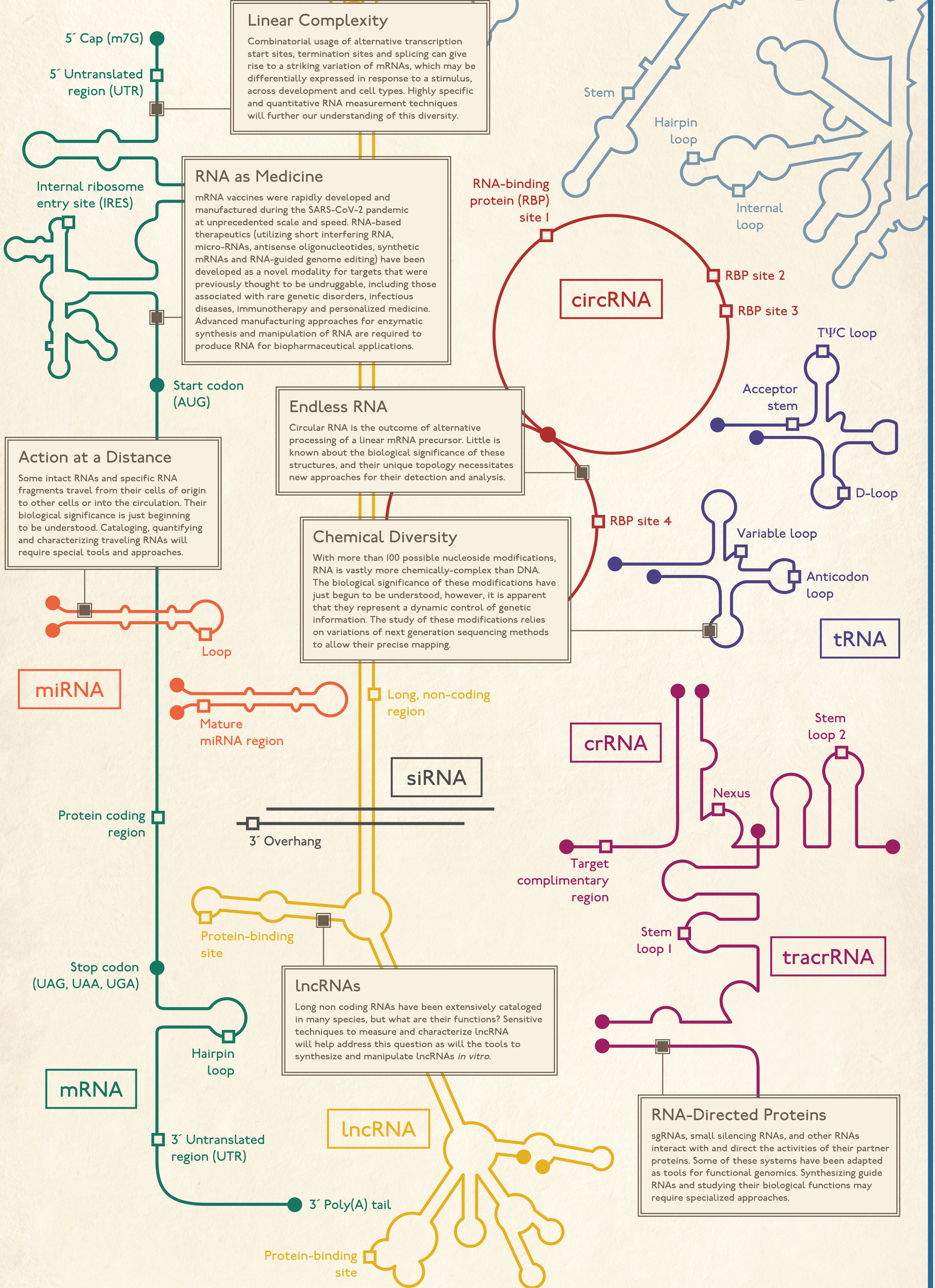


RNA METRO MAP



LET PRODUCTS FROM NEB HELP YOU NAVIGATE THE COMPLEXITY OF RNA

RNases

ShortCut® RNase III
RNase H & RNase III
XRN-I
Monarch® RNase A

RNase Inhibitors

RNase Inhibitor, Human Placenta & Murine
Ribonucleoside Vanadyl Complex

RNA Modifying Enzymes

mRNA Cap 2'-O-Methyltransferase
5' Deadenylase
RNA 5' Pyrophosphohydrolase (RppH)
Vaccinia Capping System
Faustovirus Capping Enzyme
Sce PUS1
NudC Pyrophosphatase
mRNA Decapping Enzyme

Nucleotides & Cap Analogs

Pyrophosphatase
3'-Biotin-GTP
3'-Desthiobiotin-GTP
m7G(5')ppp(5')G RNA Cap Structure Analog
m7G(5')ppp(5')A RNA Cap Structure Analog
G(5')ppp(5')A RNA Cap Structure Analog
G(5')ppp(5')G RNA Cap Structure Analog
3'-O-Me-m7G(5')ppp(5')G RNA Cap Structure Analog

In Vitro Transcription Kits

HiScribe® T7 High Yield RNA Synthesis Kit
HiScribe T7 Quick High Yield RNA Synthesis Kit
HiScribe T7 ARCA mRNA Kit with and without tailing
HiScribe SP6 RNA Synthesis Kit
HiScribe T7 mRNA Kit with CleanCap® Reagent AG
EnGen® sgRNA Synthesis Kit, *S. pyogenes*

RNA Polymerases

RNA Polymerases (T7, SP6, T3)
E. coli RNA Polymerase, Core & Holo Enzyme
E. coli Poly(A) Polymerase
Hi-T7® RNA Polymerase

Ligases

T4 RNA Ligase I (ssRNA Ligase)
T4 RNA Ligase 2
T4 RNA Ligase 2, truncated
T4 RNA Ligase 2, truncated KQ
5' DNA Adenylation Kit
SplintR® Ligase
Thermostable 5' AppDNA/RNA Ligase
T4 RNA Ligase 2, truncated K227Q
RtcB Ligase
Poly(U) Polymerase

Ladders

ssRNA & dsRNA Ladders
Low Range ssRNA Ladder
microRNA Marker

RT-qPCR & RT-PCR

Luna® Universal One-Step RT-qPCR Kit
Luna Universal Probe One-Step RT-qPCR Kit
Luna Probe One-Step RT-qPCR 4X Mix with UDG
LyoPrime Luna™ Probe One-Step RT-qPCR Mix with UDG
LunaScript® RT SuperMix Kit
LunaScript RT Master Mix Kit (Primer-free)
LunaScript Multiplex One-Step RT-PCR Kit
OneTaq® One-Step RT-PCR Kit
OneTaq RT-PCR Kit

Purification & Isolation

Monarch Total RNA Miniprep Kit
Monarch RNA Cleanup Kits (10 µg, 50 µg, 500 µg)
Epimark® N6-Methyladenosine Enrichment Kit
Oligo (dT)₂₅-Cellulose Beads
Magnetic mRNA Isolation Kit

NEBNext® Reagents for RNA-seq

Ultra™ II RNA Library Prep Kits (Directional & Non-directional)
rRNA Depletion Kit v2 (Human/Mouse/Rat)
Globin & rRNA Depletion Kit (Human/Mouse/Rat)
rRNA Depletion Kit (Bacteria)
RNA Depletion Core Reagent Set, for customized depletion
Immune Sequencing Kit (Human & Mouse)
Poly(A) mRNA Magnetic Isolation Module
Multiplex Small RNA Prep Sets
Single Cell/Low Input RNA Library Prep Kit

Get started today! Next stop: NEBna.com

Printed in the USA on recycled paper (50% PCF).
One or more of these products are covered by patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc. The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read, rather, it indicates where the document was originally developed. For more information, please email us at usabio@neb.com. The use of these products may require you to obtain additional third party intellectual property rights for certain applications.
Your purchase, acceptance, and/or payment of and for NEB's products is pursuant to NEB's Terms of Sale at www.neb.com/support/terms-of-sale. NEB does not agree to and is not bound by any other terms or conditions, unless those terms and conditions have been expressly agreed to in writing by a duly authorized officer of NEB.
CleanCap® and products incorporating it are sold under license from TruLink BioTechnologies, LLC and may be used for research use only, not for use in diagnostics, therapeutic procedures or for use in humans. For additional information, please see the product's Limited Use License at <https://www.trulinkbiotech.com/cleancap-research-license>.
© Copyright 2022, New England Biolabs, Inc., all rights reserved.

NEW ENGLAND
BioLabs Inc.

be INSPIRED
drive DISCOVERY
stay GENUINE