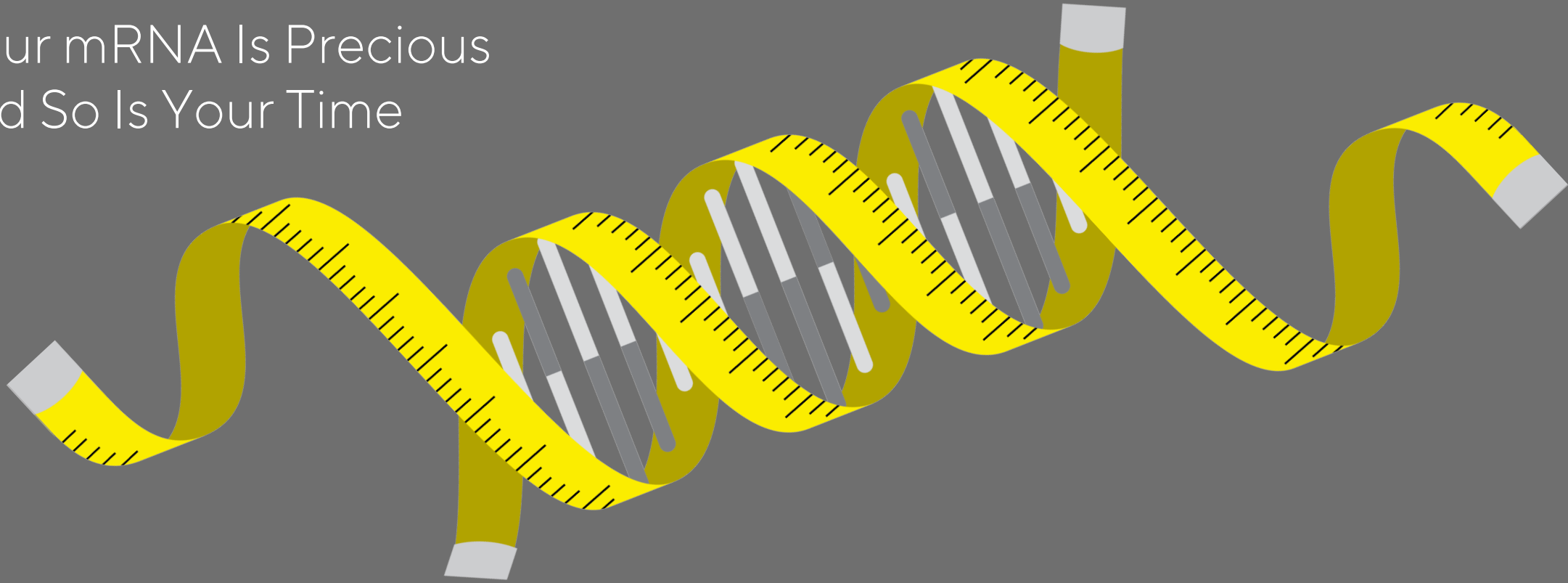
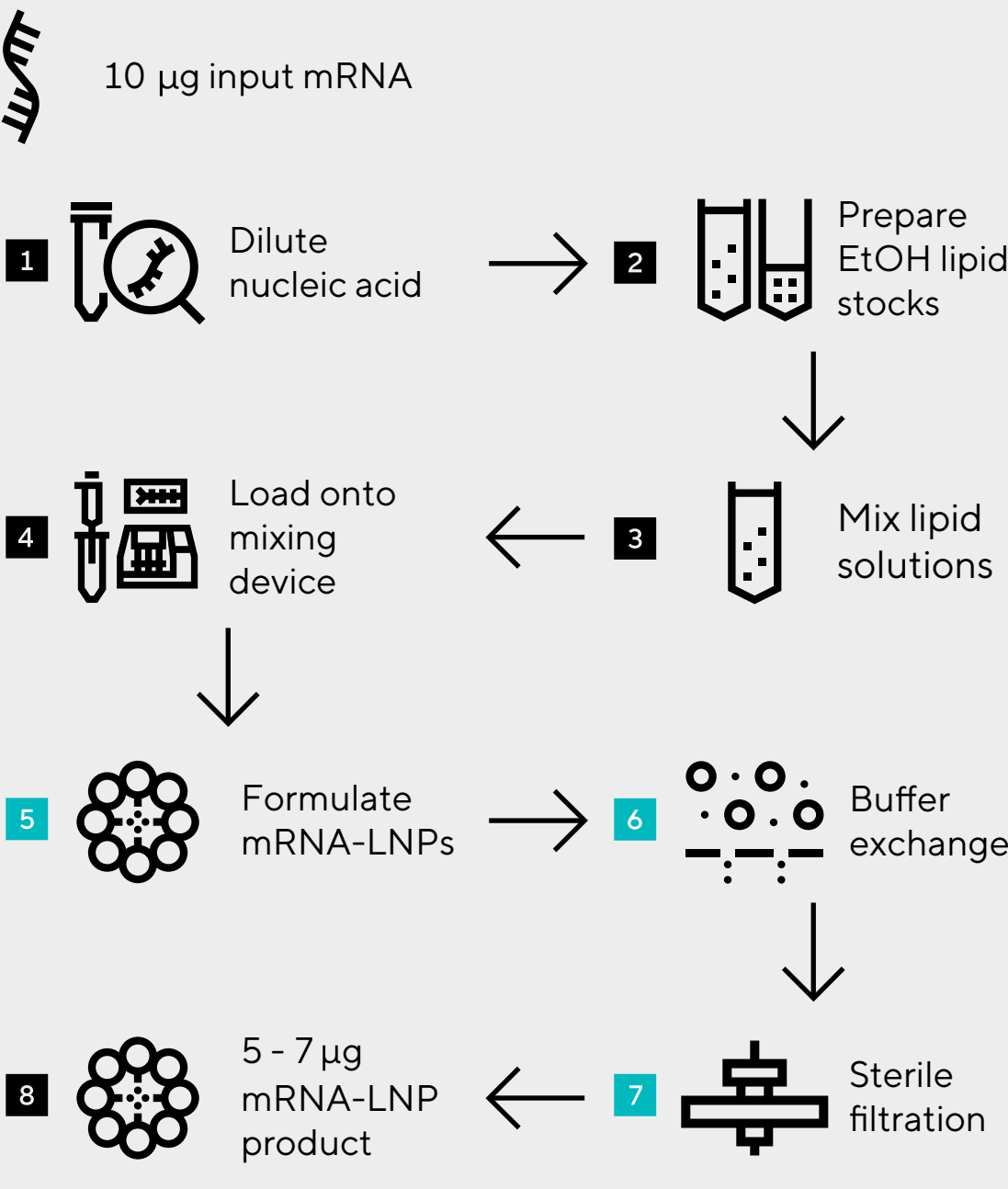


Your mRNA Is Precious
and So Is Your Time



8-Step Protocol With Standard LNPs





Lengthy



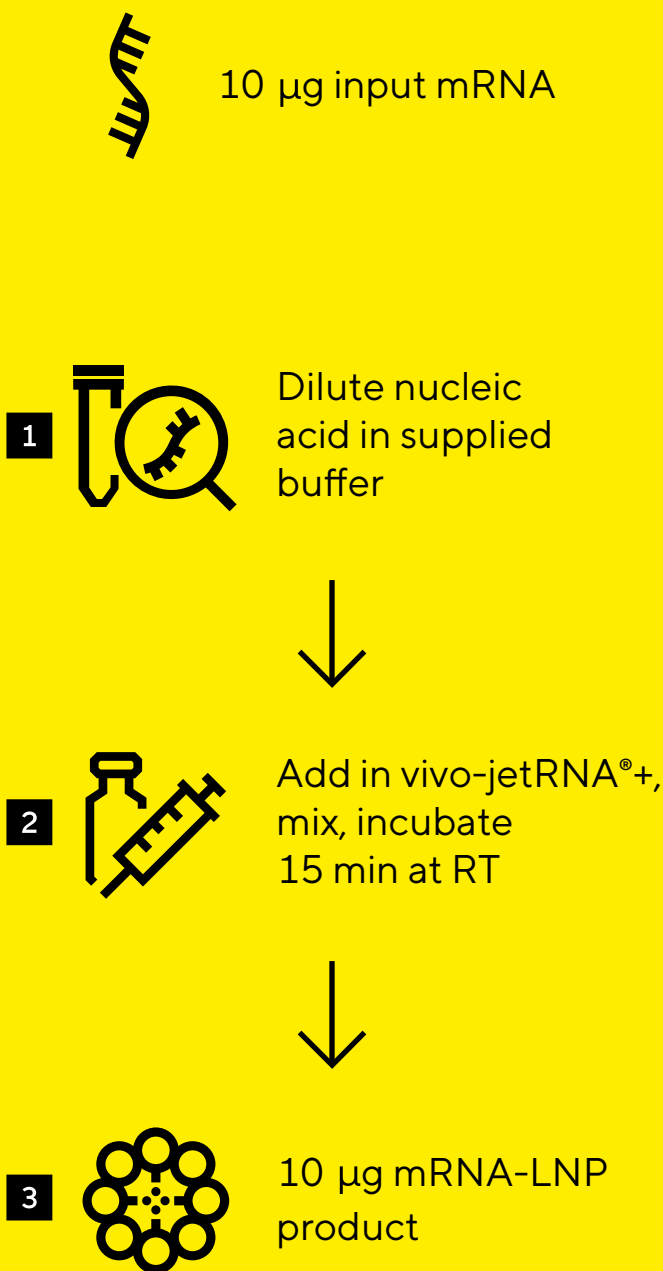
Equipment



≥30% mRNA Loss

567

3-Step Protocol With in vivo-jetRNA®+





15 min

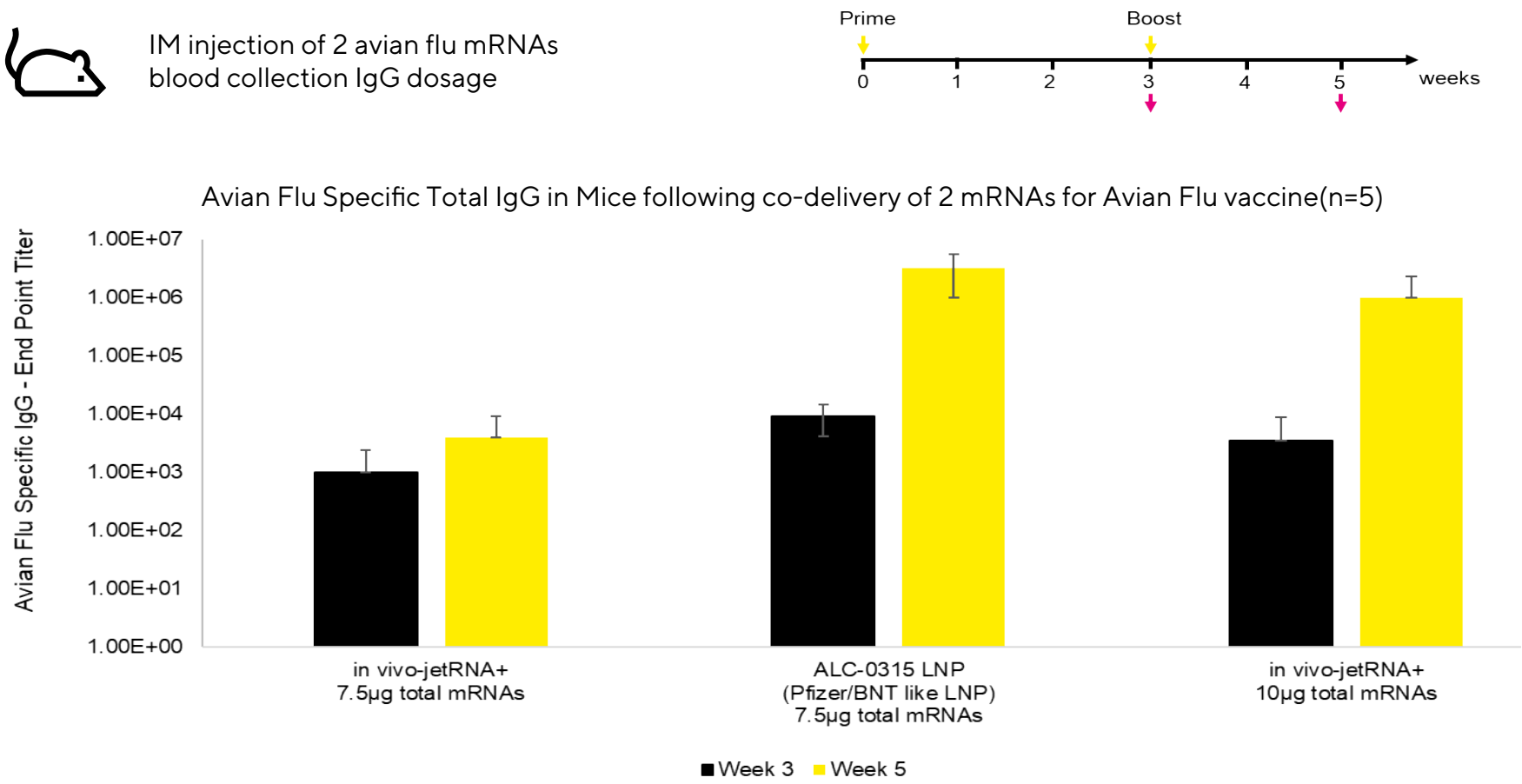


No equipment



No mRNA loss

Achieve Cost-Efficient LNP-Level Immunization with in vivo-jetRNA®+



Scan the
QR code to
learn more
about RNA

