

# in vivo-jetRNA®+

A Ready-to-Use Lipid-Based Nanoparticle Solution for In Vivo and Ex Vivo mRNA Delivery

## Product Information

in vivo-jetRNA®+ is a non-viral delivery solution composed of pre-formed lipid-based nanoparticles. It has been specifically developed for in vivo mRNA delivery. Its high delivery efficiency combined with its safety make in vivo-jetRNA®+ an innovative alternative to lipid nanoparticles (LNPs), with the added advantage of being user-friendly with a two-step protocol.

in vivo-jetRNA®+ can be used to deliver mRNA, self-amplifying RNA (saRNA), and guide RNA (gRNA) via systemic or local administration routes in various animal models, making it an ideal tool for scientists developing vaccines, protein replacement therapies, oncology treatment, or working on genome editing. It has also proven highly effective for ex vivo applications, such as the generation of mRNA CAR-T cells.



## Features and Benefits

- **Efficient:** Delivery efficiency comparable to LNPs
- **Time-Saving:** Ready-to-use, no equipment or formulation expertise needed
- **Safe:** Low pro-inflammatory effects in animal models
- **Stable:** Efficiency maintained for one month at 4 °C
- **Universal:** mRNA, saRNA, and gRNA delivery via any administration route

# Introduction

## Relevant Applications

- mRNA vaccines
- mRNA therapeutics
- Protein replacement therapy
- Gene therapy
- Genome editing
- Ex vivo cell therapy
- Proof of concept studies
- Academic research

## Relevant Process Steps

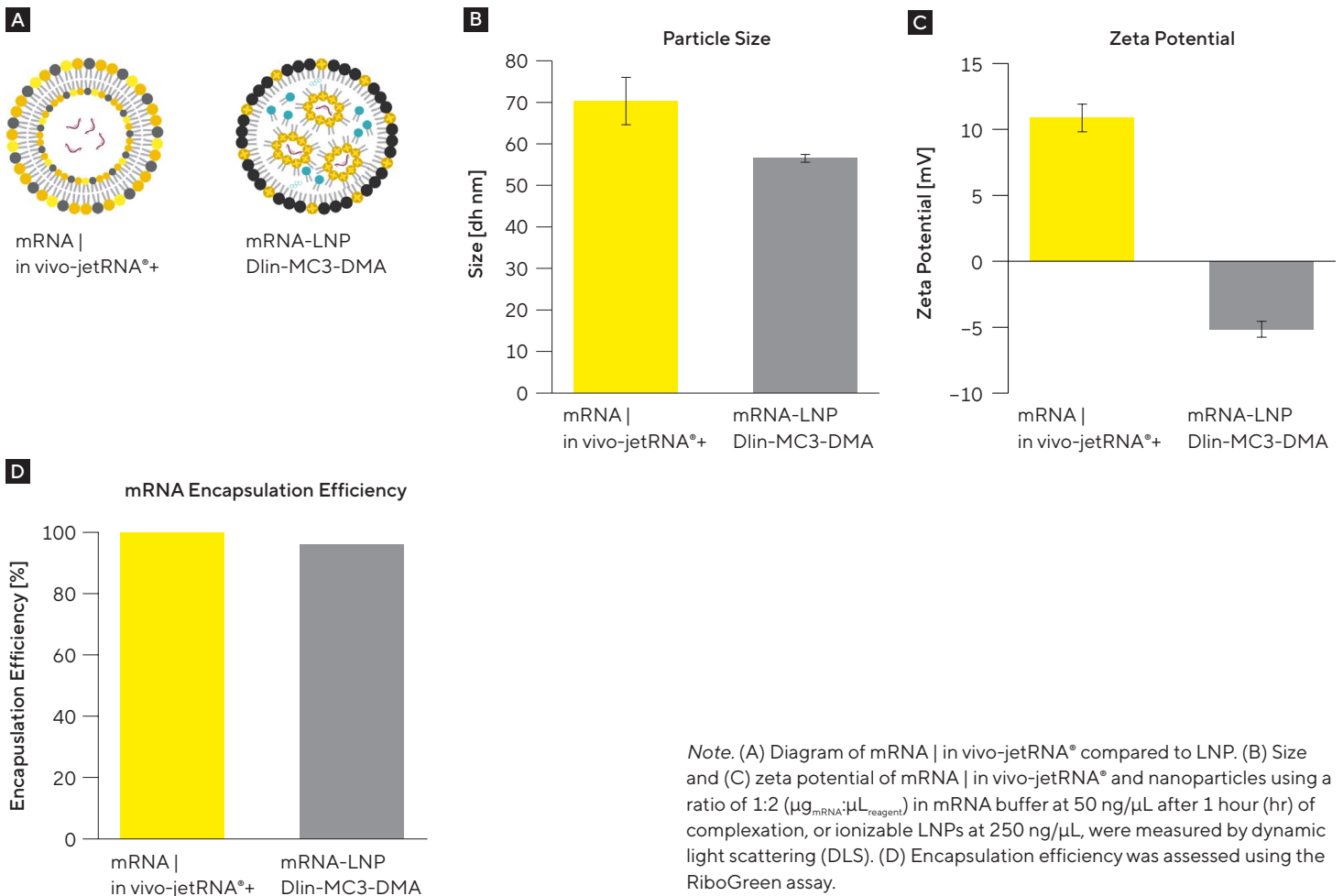
- mRNA nanoparticles formulation
- Drug product formulation
- Delivery | transfection

## Performance

### Efficient

*in vivo*-jetRNA<sup>®</sup>+ is an optimal reagent for delivering mRNA *in vivo* due to its ability to protect its payload from ubiquitous endonucleases, prevent non-specific interactions with proteins, and promote efficient cell entry. *in vivo*-jetRNA<sup>®</sup>+ has been demonstrated to encapsulate 100% of the mRNA, leading to gene expression comparable to LNPs currently used as the standard for non-viral delivery (Figure 1).

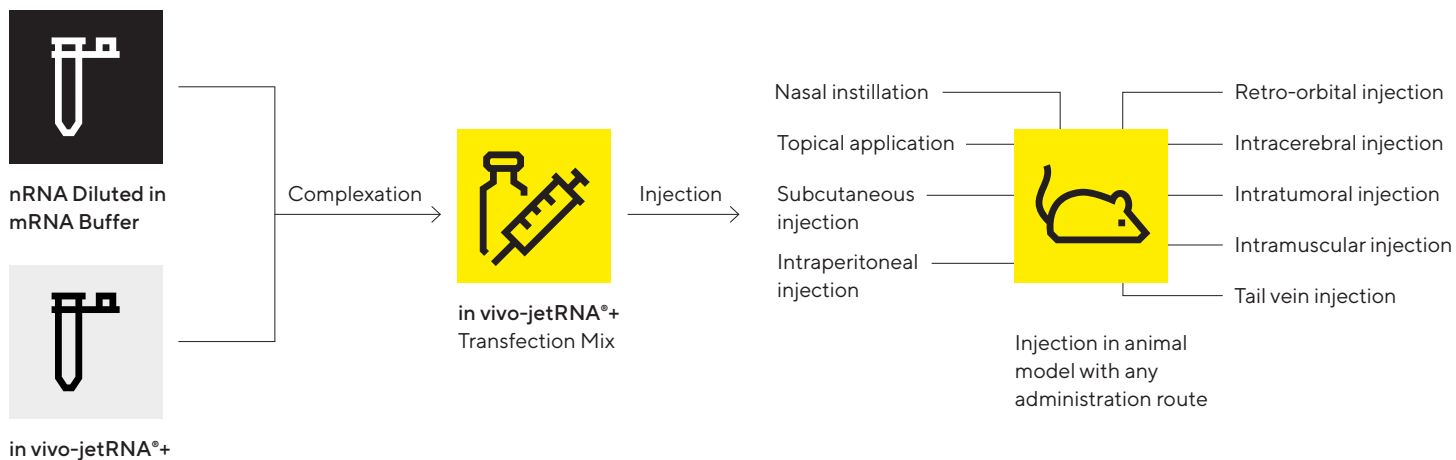
**Figure 1:** *in vivo*-jetRNA<sup>®</sup>+ Leads to 100% mRNA Encapsulation, Enabling Efficient mRNA Delivery



## Time-Saving

LNPs are widely used for mRNA delivery, but their optimal formulation requires expertise, months of work, and specific equipment and consumables. With *in vivo-jetRNA*<sup>®+</sup>, there is no need to spend time and budget on formulation, as it is a ready-to-use optimized formulation with a simple protocol (Figure 2). mRNA | *in vivo-jetRNA*<sup>®+</sup> lipid-based nanoparticles are prepared in just two steps and ready to use in 15 minutes, with no need for formulation equipment.

**Figure 2:** Protocol of *in vivo-jetRNA*<sup>®+</sup> for In Vivo mRNA Delivery

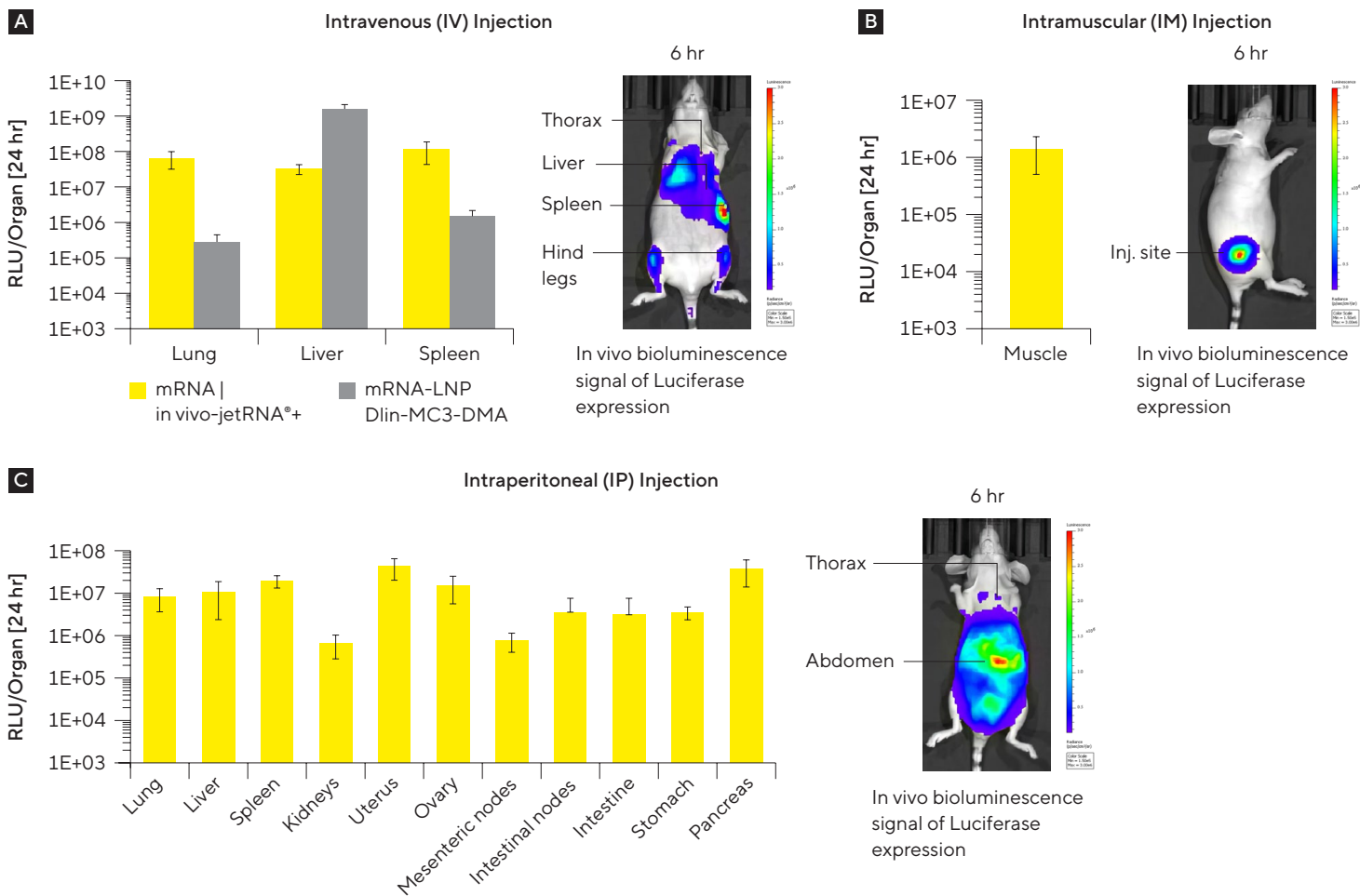


*Note.* This two-step protocol is suitable for direct injection of mRNA | *in vivo-jetRNA*<sup>®+</sup> nanoparticles by any administration route, local or systemic.

## Universal

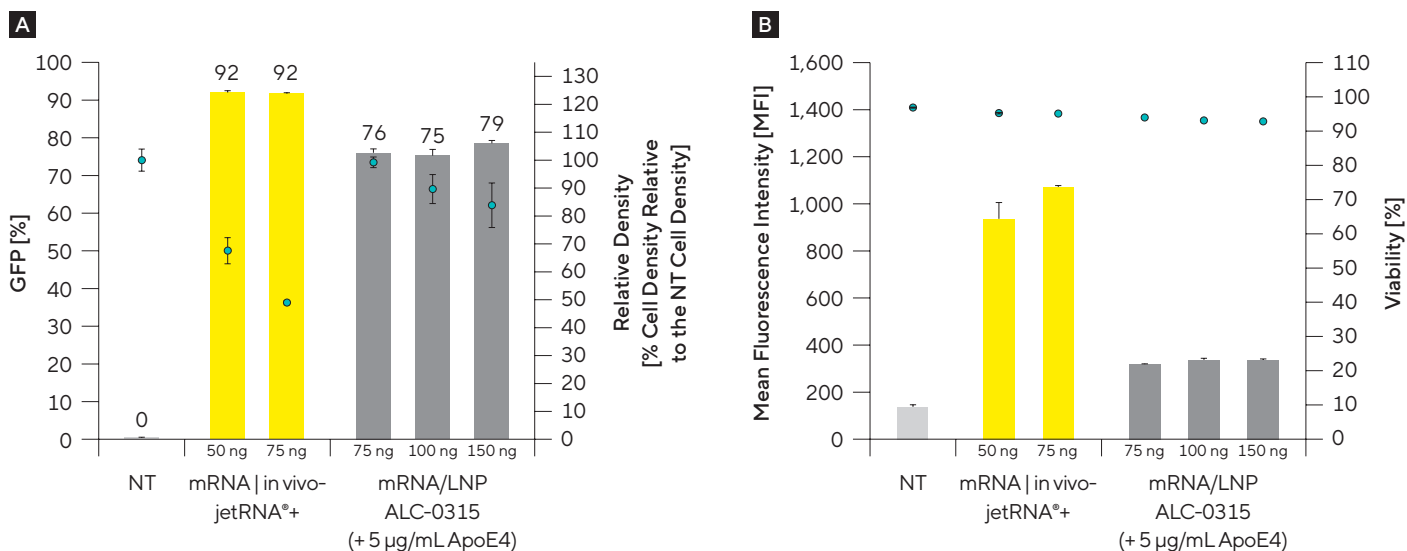
*in vivo-jetRNA*<sup>®</sup>+ has been specifically developed for the delivery of mRNA *in vivo* and is also highly efficient with saRNA and gRNA. It can be administered systemically or locally, allowing scientists to tailor biodistribution to their needs and objectives (Figure 3). In addition, *in vivo-jetRNA*<sup>®</sup>+ has also proved highly effective for *ex vivo* applications, such as the generation of transient mRNA CAR-T cells (Figure 4).

**Figure 3:** *in vivo-jetRNA*<sup>®</sup>+ Supports to Efficient mRNA Delivery in Different Organs Depending on the Administration Route



Note. mRNA encoding Luciferase was injected into mice using *in vivo-jetRNA*<sup>®</sup>+ through different administration routes: (A) intravenous (IV), (B) intramuscular (IM), and (C) intraperitoneal (IP). Complexes were formed with a mRNA | *in vivo-jetRNA*<sup>®</sup>+ ratio of 1:2 ( $\mu\text{g}_{\text{mRNA}};\mu\text{L}_{\text{reagent}}$ ) in mRNA Buffer using 10  $\mu\text{g}$  mRNA for IV, 5  $\mu\text{g}$  mRNA for IM, and 20  $\mu\text{g}$  mRNA for IP injection. Luciferase expression was assessed 6 and 24 hours post-injection.

**Figure 4:** *in vivo-jetRNA*<sup>®</sup>+ Supports Very High Ex Vivo Transfection Efficiency of Primary Human T Cells



Note. mRNA encoding GFP was transfected into primary human T cells using *in vivo-jetRNA*<sup>®</sup>+ and ionizable LNPs (based on ALC-0315). Complexes were formed with a mRNA | *in vivo-jetRNA*<sup>®</sup>+ ratio of 1:2 ( $\mu\text{g}_{\text{mRNA}}:\mu\text{L}_{\text{reagent}}$ ) in mRNA Buffer using 50 or 75 ng mRNA or with ALC-0315-LNPs (coated with 5 µg/mL ApoE4) using 75, 100, or 150 ng mRNA for 187,500 activated primary T cells in 67.5 µL of X-VIVO™ 15 (Lonza) in a 96-well plate. A further 175 µL was added after 4 hours. GFP expression, relative cell density, mean fluorescence intensity and viability were assessed 24 hours post-transfection .

## Technical Specifications

Attribute	<i>in vivo-jetRNA</i> <sup>®</sup> + 1 mL
Quality grade	Research grade
Type	Pre-formed lipid-based nanoparticle solution
Volume	1 mL
Container	Polypropylene vial
Storage	5 ± 3°C. Do not freeze
Expiry date	Indicated in the certificate of analysis and on the product
Provided with	mRNA buffer (60 mL bottle)
Number of injections	Sufficient for 25 to 50 intravenous injections or 50 to 100 intramuscular injections in mice

## Ordering Information

Item	Description	Volume	Order Number
<i>in vivo-jetRNA</i> <sup>®</sup> + 1 mL	One 1 mL vial of <i>in vivo-jetRNA</i> <sup>®</sup> + reagent supplied with one 60 mL bottle of mRNA buffer	1 mL vial (+ 60 mL mRNA buffer bottle)	101000122

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

**France**

Polyplus®—Now part of Sartorius  
75 Rue Marguerite Perey  
67400 Illkirch  
Phone +33 390 406 180



**For more information, visit**

[sartorius.com/transfection-reagents-plasmids](https://sartorius.com/transfection-reagents-plasmids)