

LipidBrick®: Innovative Cationic Lipids Fuel Next-Gen LNPs for Targeted Oncology and Advanced Cell Therapies

Pauline Urquiza¹, Quentin Huauilmé¹, Mélodie Seiler¹, Marine Fenat¹, Margaux Briand¹, Kassandra Renaud¹, Julie Chevrier¹, Thibaut Benchimol¹, Malik Hellal², Claire Guéguen¹, Patrick Erbacher¹

¹Sartorius Polyplus, Vectura, 75 rue Marguerite Perey, 67400 Illkirch, France

²Sartorius Stedim FMT S.A.S. Zone Industrielle les Paluds, Avenue de Jouques, 13400 Aubagne, France

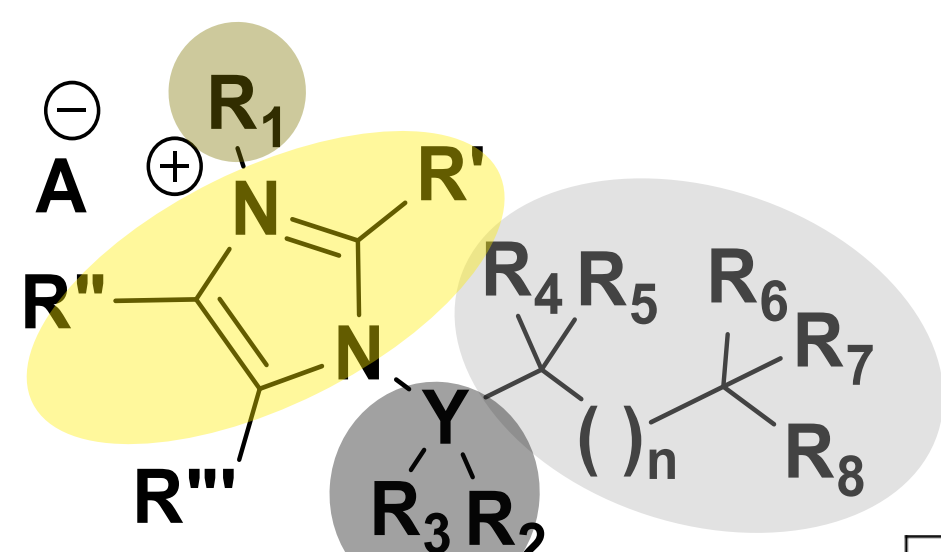
Lipid nanoparticles (LNPs) have shown promising potential in organ selective targeted -Oncology and Gene Modified Cell Therapies; however, the critical challenge lies in fine-tuning their formulation to enhance specificity, efficiency, and safety, to maximize therapeutic outcomes.

Here, we present a novel approach through the development of lipid nanoparticles (LNPs) derived from our library of Innovative cationic lipids. These LNPs demonstrate expanded extrahepatic biodistribution, improving drug delivery and cellular targeting. This addresses major challenges in cancer treatment for the development of therapeutic vaccines via targeting antigen presenting cells (APCs) and T cells in the spleen, and organ selective oncology treatments via local injection. Optimization of LNP formulations around our cationic lipids leads to localized expression without off-target effects unlike commonly used ionizable lipid-based LNPs.

In addition, we investigate ex vivo cell engineering, employing LNPs. We highlight the efficacy and versatility of LNPs based on our library of lipids, which enable the transfection of mRNA in different types of immune cells, leading to transient engineering of CAR-T and NK cells. This data suggests that LNP formulations based on our cationic lipids, with the addition of targeting ligands, could also prove effective for the generation of CAR-Ts and NKs in vivo. The data underpinning these applications were collected through a blend of proof-of-concept studies conducted both in vitro and in vivo, aimed at evaluating the efficacy and safety of imidazolium lipid-based LNP systems. Molecular assays were employed to monitor biodistribution and cellular uptake, offering real-world insights with an emphasis on safety and dosage optimization

LipidBrick® Library: A Range of Proprietary Cationic Lipids

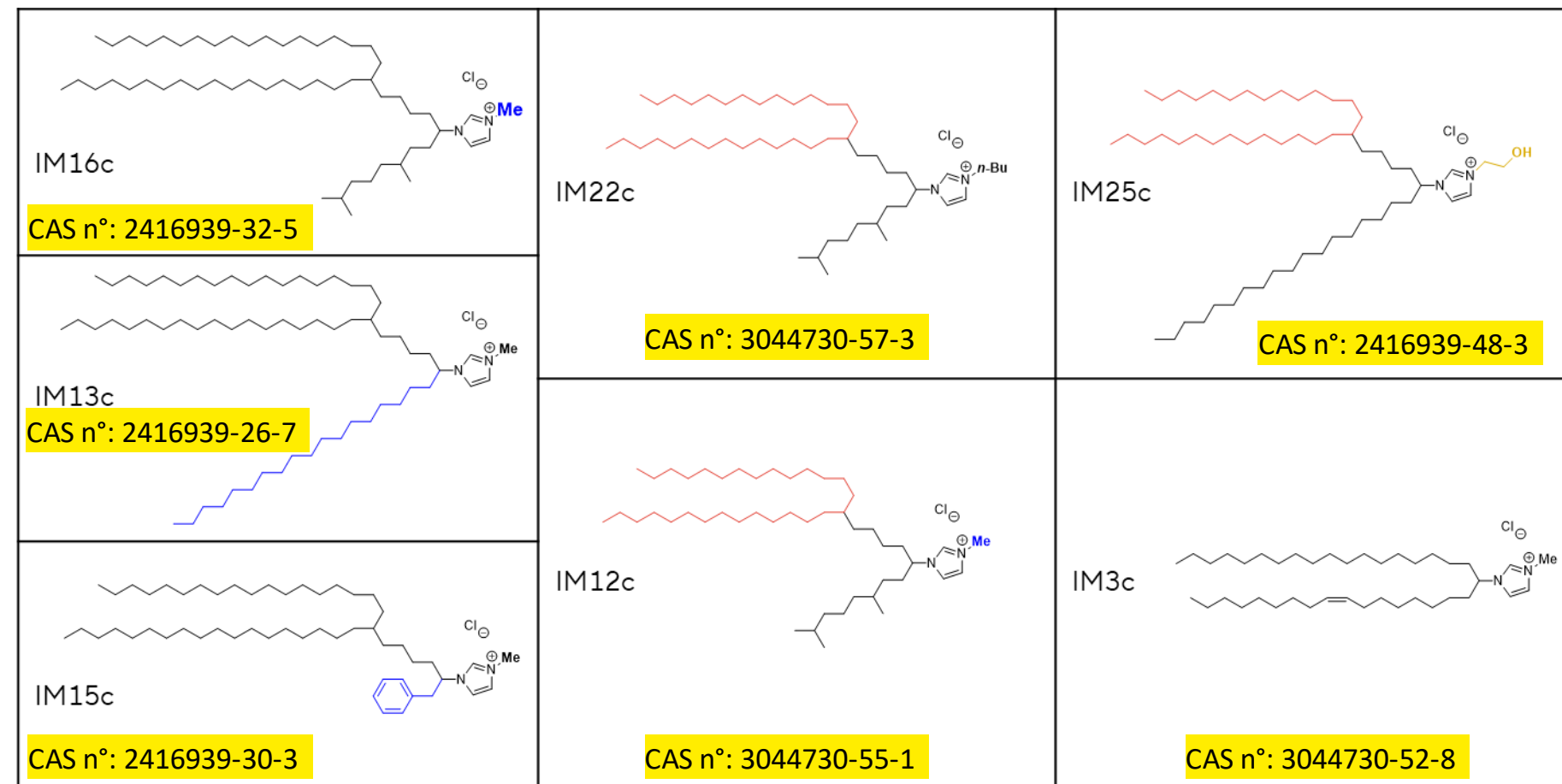
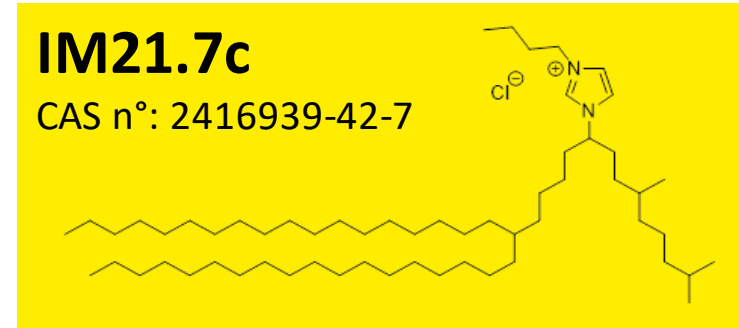
Imidazolium-Based Cationic Lipids: Offering New Targeting Options and Improving Safety



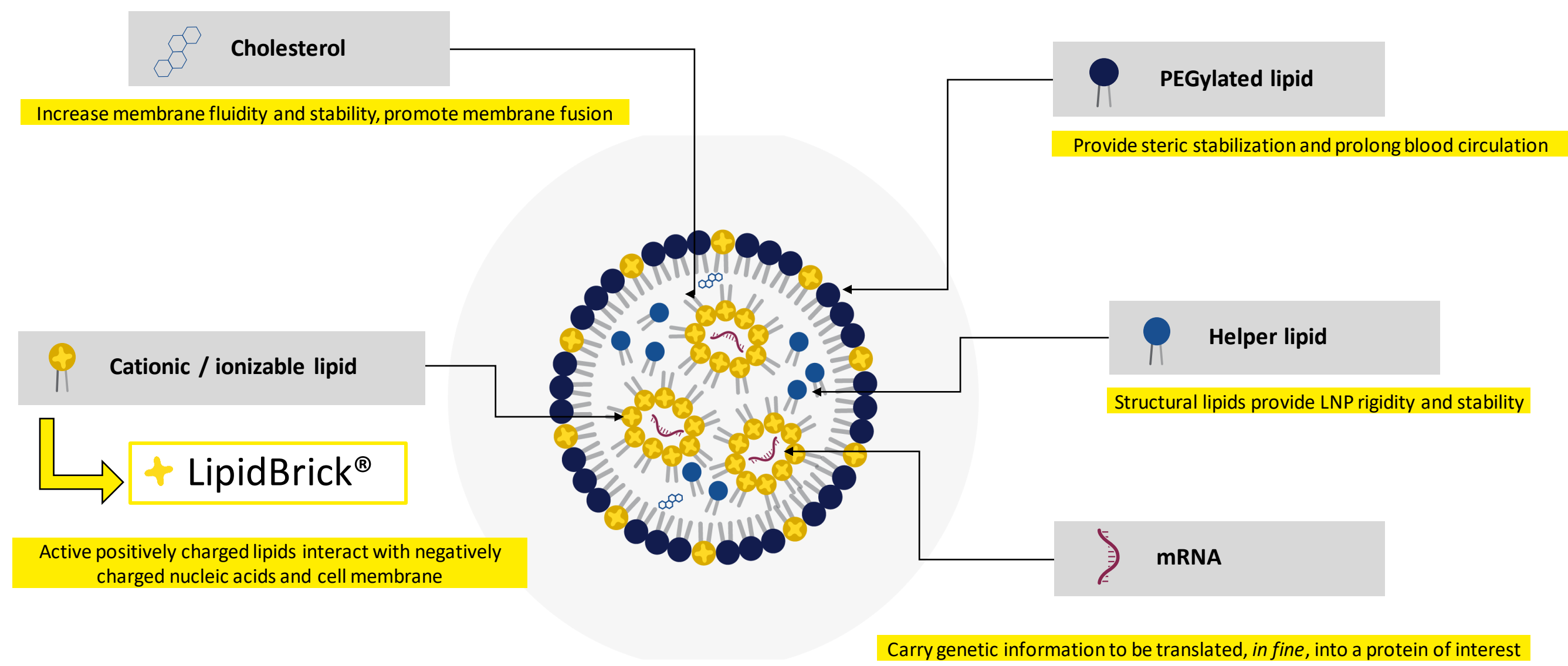
➤ New generation of cationic lipids dedicated to LNPs or liposomes formulation.

➤ Featuring an Imidazolium polar head, resulting in Improved Safety compared to traditional cationic lipids.

➤ Expanded into an 8-lipid library for greater flexibility and screening options.



Enabling Formulation of Suitable 4- and 5-lipid Lipid Nanoparticles



➤ LipidBrick® Library is composed of permanently charged cationic lipids that play a key role in LNP composition, functioning as the active lipid that electrostatically binds with the genetic payload to enable improved encapsulation efficacy, cellular uptake and a broadened biodistribution beyond the liver.

Powering Tailored LNP Formulations Across Cell Types

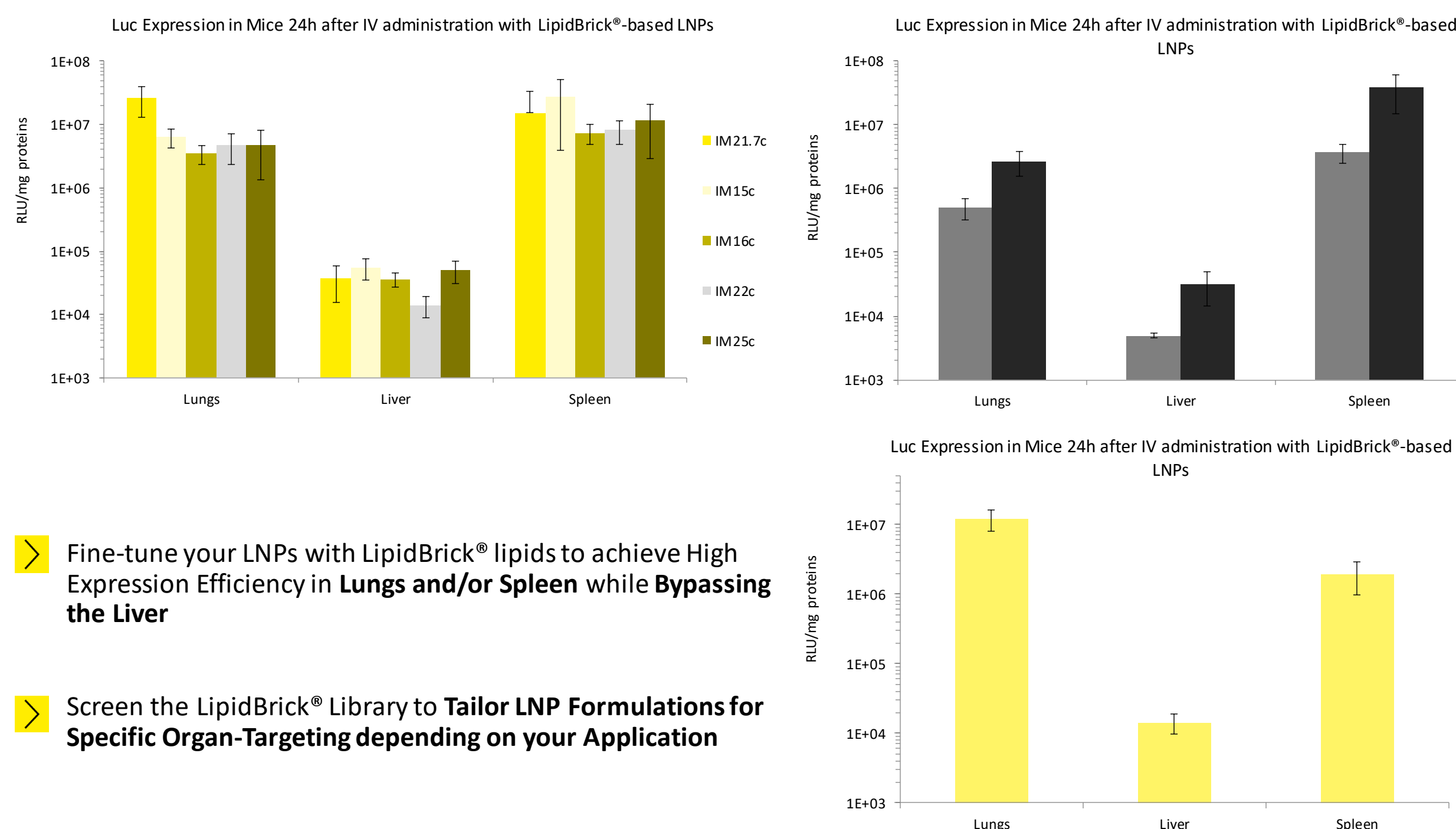
Screening conditions:		Performance of best LNP Formulation for each LipidBrick® lipid:							
LNP Composition	Molar ratio range (%)	LipidBrick®	IM3c	IM12c	IM13c	IM15c	IM16c	IM21.7c	IM22c
LipidBrick® IMXXc	20 – 50	HepG2							
Phospholipid (DOPE or DSPC)	10 – 40	HEK-293							
Cholesterol	28.5 – 48.5	C2C12							
DMG-PEG2k	1.5	Huh7							
		A-498							
		Caco-2							
		Primary T cells							

High-Throughput Screening of Diverse LipidBrick® Lipid-Based LNPs for Encapsulation and Delivery of Fluc mRNA: 25-100 ng of mRNA-LNPs were applied per well in 96-well plates, with luciferase expression evaluated 24- or 48-hours post-transfection across various cell types in vitro.

➤ Each application requires tailored LNP formulations to achieve optimal results (one size ≠ fit all!). LipidBrick® Library offers diverse options with varied chemical structures, facilitating high-throughput screening (HTS).

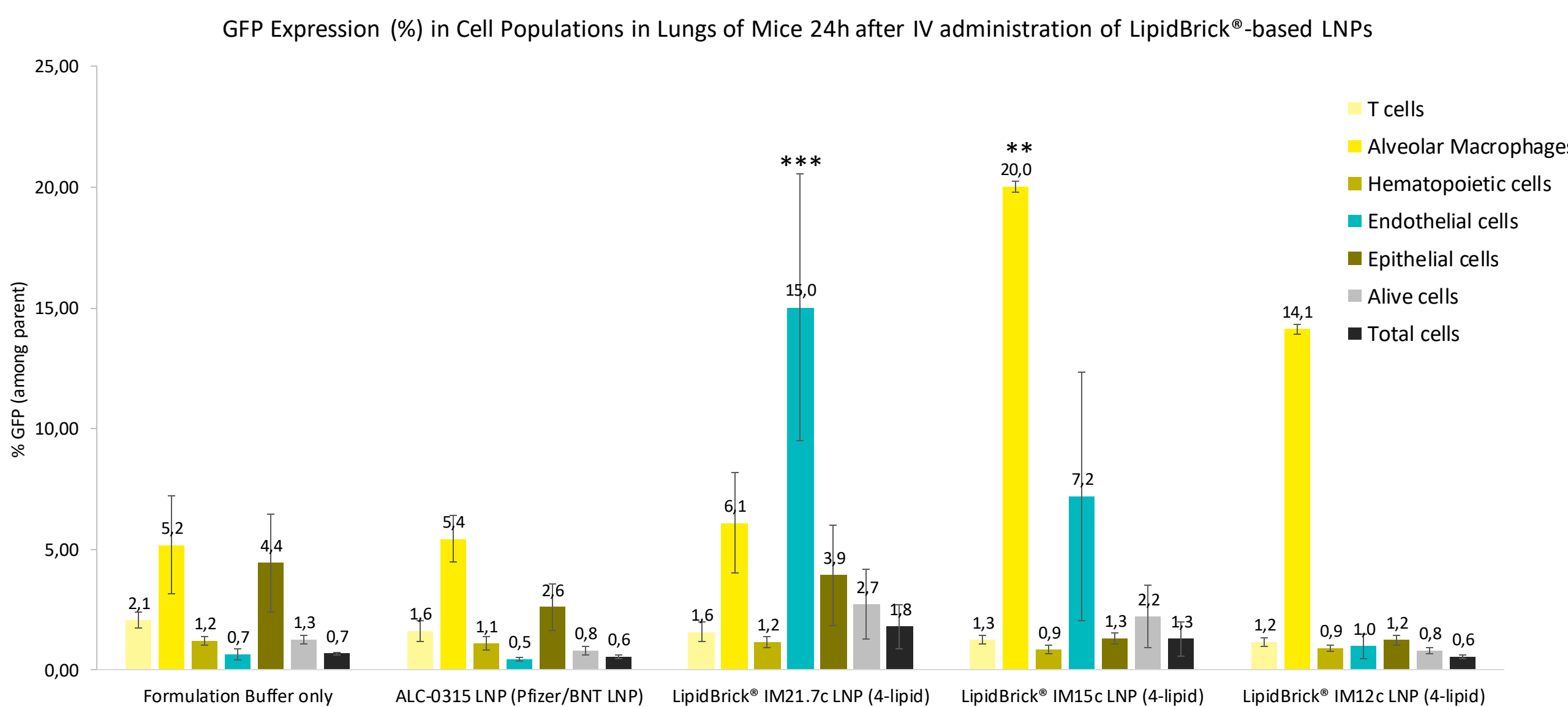
Targeted Oncology: LipidBrick® Library for Targeted Cancer Therapies

Tailored Formulations: Providing to modulate Biodistribution



10µg of mRNA encoding Fluc were injected through retro-orbital (RO) injection in mice. Luciferase expression was assessed 24 hours post-injection.

Empowering RNA Therapies: Cell-Specific Targeting with LipidBrick®-based LNPs

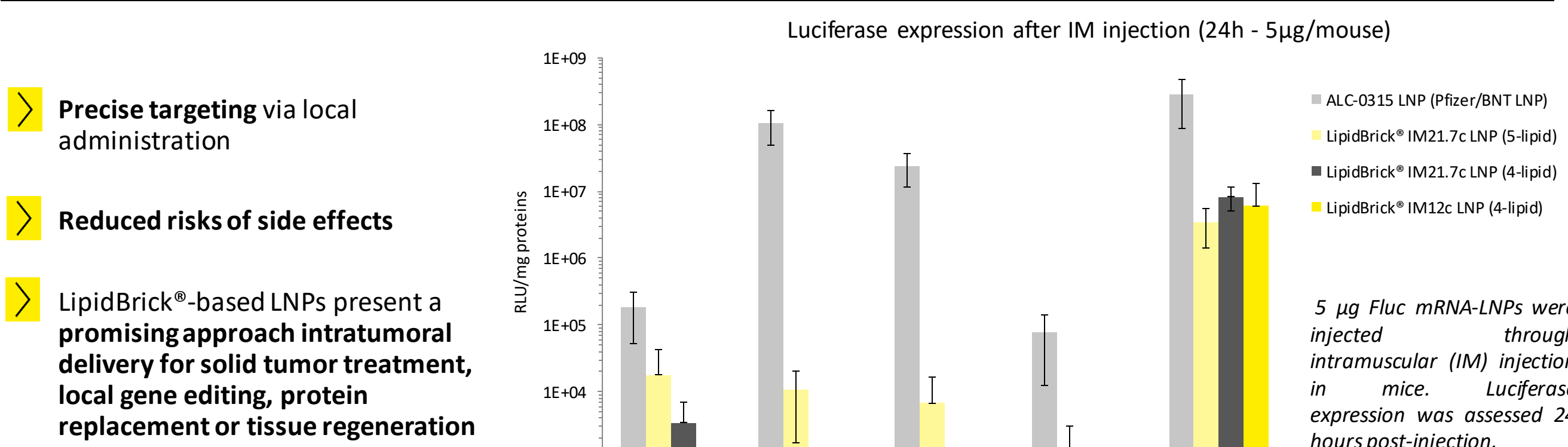


10 µg CleanCap EGFP 5mU mRNA-LNPs at N/P ratio of 6.8 were injected through retro-orbital (RO) injection in mice. Lungs were harvested 24 hours post-injection, dissected and labelled before being analyzed by flow cytometry. One-way ANOVA, Tukey's multiple comparison vs. Formulation buffer only, **p<0.01 and ***p<0.001.

➤ Screen the LipidBrick® Library to Fine-Tune your LNP Formulations for Specific Cell-Targeting depending on your application.

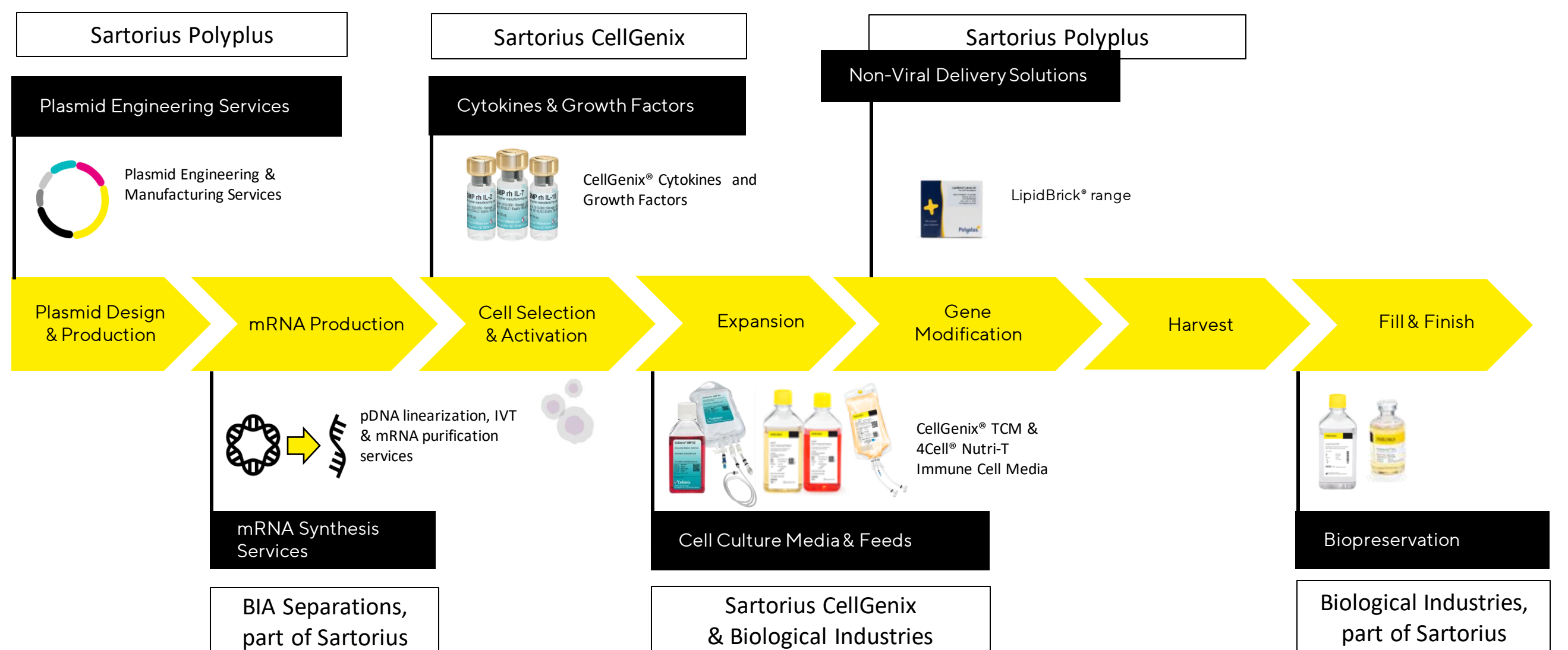
➤ LipidBrick®-based LNPs present a promising approach for high transfection efficacy in endothelial cells and/or alveolar macrophages, key targets for innovative lung disease therapies, including lung cancer.

Fine-Tuning LNPs with LipidBrick® Library to Enhance Targeting Specificity

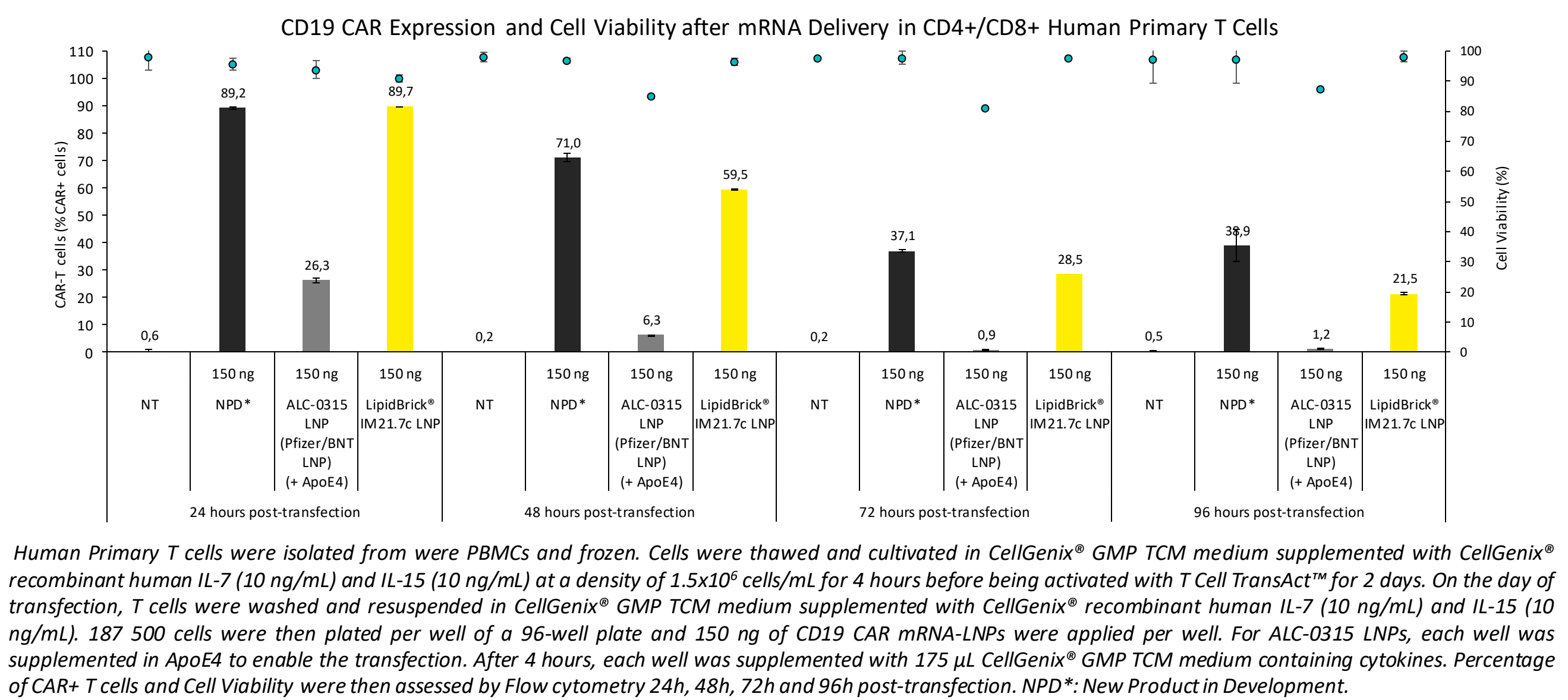


Advanced T Cell Therapy: LipidBrick® for Non-Viral Delivery

Comprehensive Portfolio for Non-Viral Gene-Modified Cell Therapy Solutions



Enhanced & Durable CD19 CAR mRNA Expression in T Cells with LipidBrick®-based LNPs



➤ LipidBrick®-based LNPs present a promising Non-Viral Delivery approach for ex vivo Gene-Modified Cell Therapy applications, such as CAR-T Cell Therapies.

➤ Data suggests LipidBrick®-based LNPs with targeting ligands could be effective for the generation of in vivo CAR-Ts and NKs

Achieve Promising Transfection Efficacy in Primary Human NK Cells

