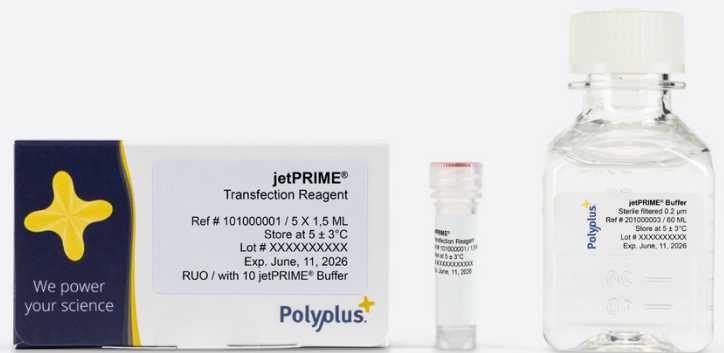


# jetPRIME®

## Versatile DNA | siRNA Transfection Reagent



### Product Information

jetPRIME® is a polyvalent transfection reagent that ensures effective and reproducible DNA and siRNA transfection in several adherent cell lines. This versatile reagent is also perfect for co-transfection of DNA and siRNA. Transfection with jetPRIME® requires only low amounts of nucleic acid, hence resulting in very low cytotoxicity.

jetPRIME® is suited for different applications, such as plasmid transfection, gene silencing, co-transfection of different nucleic acids, and CRISPR/Cas9-mediated gene knock-out.

### Benefits

- **Versatile:** One reagent for DNA and siRNA transfection
- **Timesaving:** Easy-to-use transfection reagent
- **Gentle on cells:** Low cytotoxicity
- **Cost-effective:** Requires less DNA and less reagent for efficient transfection

# Introduction

## Relevant Applications

- Plasmid transfection
- Genome editing
- siRNA transfection
- Co-delivery of different nucleic acids

## Relevant Process Steps

- Upstream Process
- Transfection

# Technical Specifications

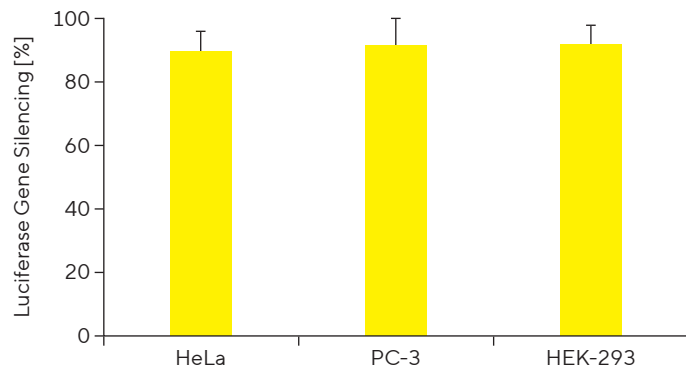
Molecule delivered	sDNA, siRNA, DNA & siRNA (co-transfection)
Cell Types	Adherent cells
Number of transfections	1.5 mL of jetPRIME® transfection reagent is sufficient to perform up to 1500 transfections in 24-well plates or 375 transfections in 6-well plates
Storage	Store jetPRIME® at 5 °C ± 3°C. Expiry date is indicated in the certificate of analysis and on the product
Provided with	jetPRIME® buffer

# Capabilities

## ▪ One Reagent for DNA and | or siRNA Transfection

jetPRIME® is a polyvalent transfection reagent well suited for DNA and siRNA delivery. jetPRIME® is also ideal for DNA | siRNA co-transfection, ensuring high DNA Transfection and efficient gene silencing in a variety of cell lines with exceptionally low toxicity.

**Figure 1:** Exogenous Luciferase Silencing in Several Cell Lines After DNA and siRNA Cotransfection Using jetPRIME®

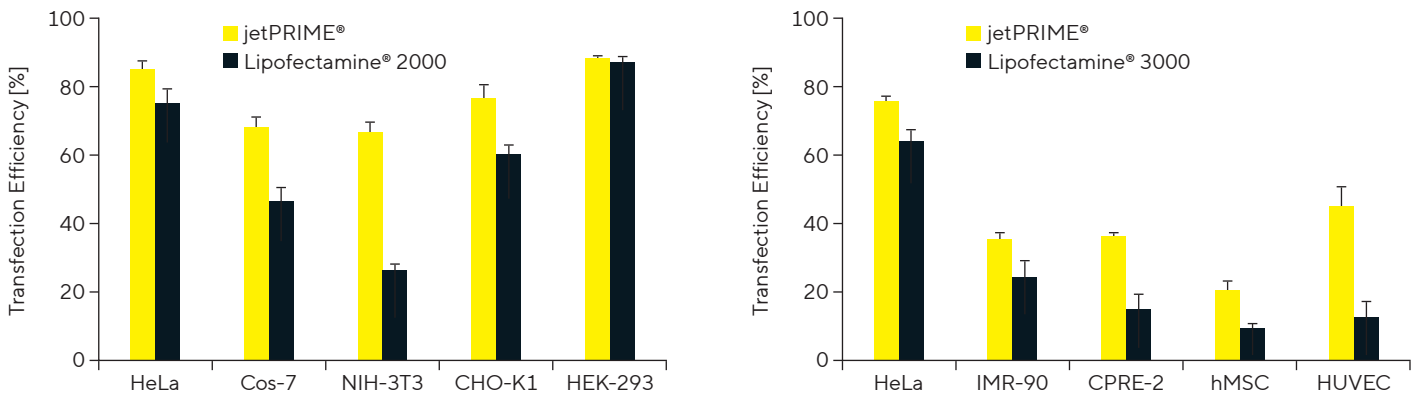


Note. Experiments were performed with 400 ng p4CMV-Luc and 10 nM of luciferase siRNA per well in 6-well plates.

▪ **High Transfection Efficiency in a Wide Variety of Adherent Cell Lines**

jetPRIME® is an easy-to-use transfection reagent for day-to-day experiments that leads to efficient and reliable scientific results. It leads to outstanding high percentage of transfected adherent cell lines of various origins as well as primary cells

**Figure 2:** Comparative Transfection Efficiency of jetPRIME® Versus Its Main Competitors

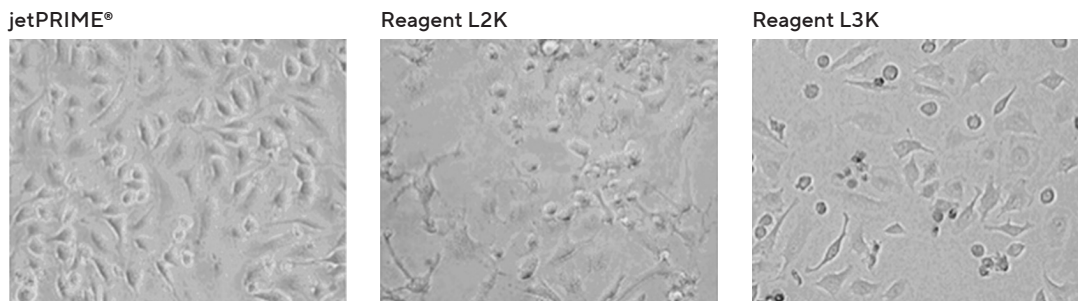


Note. Transfection efficiency was assessed by FACS analysis in various cell lines 24 hours after transfection in 24-well plates. Conditions were used according to the manufacturer's recommendation.

▪ **Extremely Gentle on Cells**

Transfection with jetPRIME® requires low amounts of reagent and nucleic acids, leading to increased cell viability while keeping high transfection efficiency.

**Figure 3:** Cells Transfected With jetPRIME® Are Healthy, While Major Cytotoxicity Is Observed With Competitors: Cell Viability 24 Hours After Transfection



Note. Phase contrast microscopy HeLa Cells 24 hours after transfections performed according to the manufacturer's recommendations for each reagent.

### ■ Requires Less DNA and Less Reagent

Transfection with jetPRIME® requires a small amount of reagent and DNA | siRNA, making its use very cost-effective.

**Table 1:** Conditions for Transfection Given per 6-Well Plates According to Manufacturer's Recommendations

Reagent	Volume of reagent per well	Amount of DNA per well	Number of transfection per 1.5 mL vial
jetPRIME®	2 - 4 µL	1 - 2 µg	375 - 750
Lipofectamine® 2000	5 - 12.5 µL	2.5 µg	120 - 300
Lipofectamine® 3000	3.75 - 7.5 µL	2.5 µg	200 - 400

## Ordering Information

Item	Description	Package   Volume   Quantity   Size	Order Number
jetPRIME® 0.1 mL	Versatile DNA   siRNA transfection reagent provided with associated Buffer	0.1 mL Vial (+ 5 mL Buffer Bottle)	101000027
jetPRIME® 0.75 mL		0.75 mL Vial (+ 60 mL Buffer Bottle)	101000015
jetPRIME® 1.5 mL		1.5 mL Vial (+ 2x60 mL Buffer Bottle)	101000046
jetPRIME® 5x 1.5 mL		5x 1.5 mL Vial (+ 10x60 mL Buffer Bottle)	101000001
jetPRIME® buffer 60 mL	jetPRIME® associated Buffer	60 mL Bottle	201000003

Bulk quantities are available upon request.

### Germany

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

### USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

 For further information, visit  
[sartorius.com](https://www.sartorius.com)