

## jetPRIME<sup>®</sup> Transfection Reagent

### Short Protocol – DNA Transfection

#### Day 0: Cell Seeding

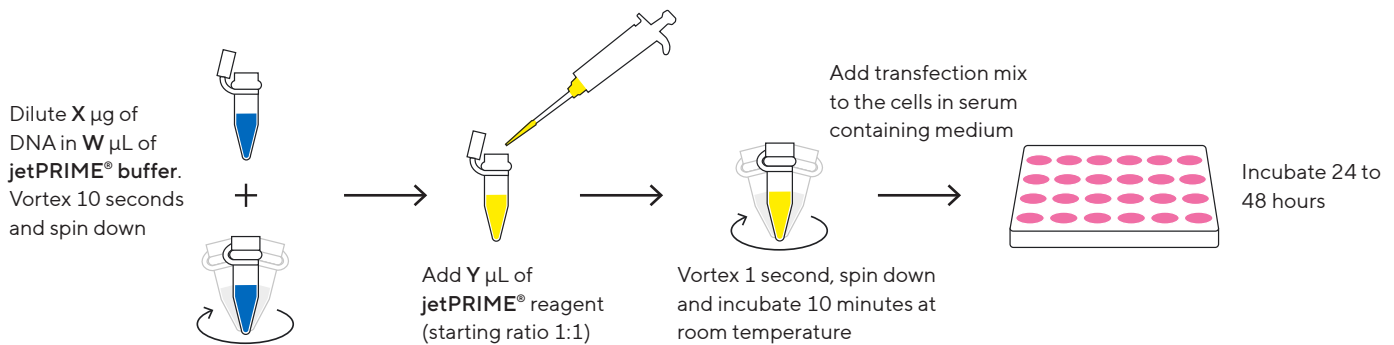
- Seed cells in **V** mL of cell growth medium according to the table below

Culture vessel	Number of cells*	V = volume of medium during transfection
96-well	7,500 - 10,000	0.1 mL
24-well	50,000 - 80,000	0.5 mL
12-well	80,000 - 150,000	1 mL
6-well/35 mm	150,000 - 250,000	2 mL
100 mm/flask 75 cm <sup>2</sup>	1 x 10 <sup>6</sup> - 2 x 10 <sup>6</sup>	10 mL

Quantities per well, dish or flask.

#### Day 1: Transfection

- Perform transfection **in the presence of serum**
- Use **jetPRIME<sup>®</sup> buffer only**
- Transfect cells at **60 - 80% confluency**



Culture vessel	1. W = volume of jetPRIME <sup>®</sup> buffer [µL]	2. X = amount of DNA added [µg]	3. Y = volume of jetPRIME <sup>®</sup> reagent [µL]
96-well	10	0.1	0.2
24-well	50	0.5	1
12-well	75	0.8	1.6
6-well/35 mm	200	2	4
100 mm/flask 75 cm <sup>2</sup>	500	10	20

Quantities per well, dish or flask.

#### Day 2 - 3: Measure Gene Expression

See back page for optimization tips.

Download complete protocol on [sartorius.com](http://sartorius.com)

# Short Protocol – Optimization Tips (DNA)

## Protocol Optimization

- Test different DNA amounts: X, 0.5X and 1.5X.
- Test different DNA/jetPRIME® ratios, 1:2 to 1:3.
- For cell specific protocols, visit [www.sartorius.com](http://www.sartorius.com).

Culture vessel	W = volume of jetPRIME® buffer [µL]	X = amount of DNA added [µg]	Y = volume of jetPRIME® reagent [µL]
96-well	10	0.05 - 0.20	0.10 - 0.60
24-well	50	0.25 - 0.75	0.50 - 2.25
12-well	75	0.4 - 1.2	0.8 - 3.6
6-well/35 mm	200	1 - 3	2 - 9
100 mm/flask 75 cm <sup>2</sup>	500	5 - 15	10 - 45

Quantities per well, dish or flask.

For HEK-293 and HeLa cells, you may decrease the DNA amount to 0.5X and use the 1:2 DNA/jetPRIME® ratio.

## Tips to Increase Cell Viability of Sensitive Cells

- Replace medium 4 hours after transfection.
- Decrease DNA amount to 0.5X while maintaining the DNA/jetPRIME® ratio previously used.
- Analyze transfection at an earlier time point (24 hours after transfection instead of 48 hours for instance).
- Check that the target gene does not affect cell viability.

## Good DNA Transfection Practices

- Store appropriately jetPRIME® (5 ± 3 °C).
- Ensure that cells have been passaged more than twice and less than 20 times prior to transfection.
- Discard overconfluent cells.
- Regularly check for mycoplasma contaminations.
- Use a reporter gene to set up and optimize transfection conditions.
- Serum quality may drastically affect transfection efficiency. When purchasing a new batch of serum or trypsin, check cell viability as well as transfection efficiency.

## NOTE:

jetPRIME® is also recommended for DNA | siRNA co-transfection. Please refer to the complete protocol on [www.sartorius.com](http://www.sartorius.com).

## 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](http://sartorius.com)