

jetOPTIMUS® Transfection Reagent

Short Protocol – DNA Transfection

Day 0: Cell Seeding

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

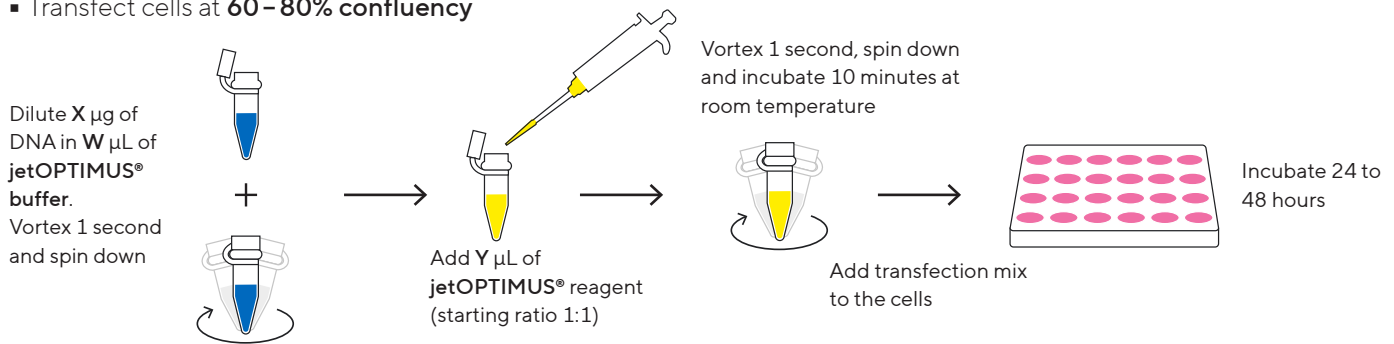
Culture vessel	Number of cells*	V = volume of medium during transfection
96-well	7,500 – 25,000	0.125 mL
24-well	40,000 – 100,000	0.5 mL
12-well	80,000 – 200,000	1 mL
6-well/35 mm	150,000 – 400,000	2 mL
60 mm/flask 25 cm ²	200,000 – 850,000	5 mL
100 mm/flask 75 cm ²	1 x 10 ⁶ – 4 x 10 ⁶	10 mL

Quantities per well, dish or flask.

*For specific cell type or suspension cells, please refer to the complete protocol.

Day 1: Transfection Using jetOPTIMUS® Reagent

- Use **jetOPTIMUS® buffer only**
- Transfect cells at **60 – 80% confluency**



Culture vessel	W = volume of jetOPTIMUS® buffer	X = amount of DNA added	Y = volume of jetOPTIMUS® reagent
96-well	12.5 µL	0.13 µg	0.13 – 0.19 µL
24-well	50 µL	0.5 µg	0.5 – 0.75 µL
12-well	100 µL	1 µg	1 – 1.5 µL
6-well/35 mm	200 µL	2 µg	2 – 3 µL
60 mm/flask 25 cm ²	500 µL	4 µg	4 – 6 µL
100 mm/flask 75 cm ²	1,000 µL	10 µg	10 – 15 µL

Quantities per well, dish or flask.

Day 2 – 3: Measure Gene Expression

See back page for optimization tips.

Download complete protocol on sartorius.com

Short Protocol – Optimization Tips

Protocol Optimization

- Test different DNA amounts: X, 0.5X and 1.5X
- Test different DNA/jetOPTIMUS® ratios, 1:1 to 1:1.5.
- For cell specific protocols, visit our website at www.sartorius.com.

Culture vessel	W=volume of jetOPTIMUS® buffer	X=amount of DNA added	Y=volume of jetOPTIMUS® reagent
96-well	12.5 µL	0.10-0.20 µg	0.10-0.30 µL
24-well	50 µL	0.25-0.75 µg	0.25-1 µL
12-well	100 µL	0.5-1.5 µg	0.5-2.25 µL
6-well/35 mm	200 µL	1-3 µg	1-4.5 µL
60 mm/flask 25 cm ²	500 µL	2-6 µg	2-9 µL
100 mm/flask 75 cm ²	1,000 µL	5-15 µg	5-22 µL

Quantities per well, dish or flask.

Tips to Increase Cell Viability of Sensitive Cells

- Replace medium 4 hours after transfection.
- Decrease DNA amount to 0.5X while maintaining the DNA/jetOPTIMUS® ratio previously used.
- Analyze transfection at an earlier time point (24 hours after transfection instead of 48 hours for instance).
- Perform transfection in reduced serum medium for sensitive cells.
- Check that the target gene does not affect cell viability.
- Ensure that cells have been passaged more than twice and less than 20 times prior to transfection.
- Discard overconfluent cells.

Good mRNA Transfection Practices

- Store appropriately jetOPTIMUS® (5 ± 3 °C).
- Regularly check for mycoplasma contamination.
- 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:

Please refer to the complete protocol available on 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