

Scaling Up or Down Lipofectamine® 3000 Reagent Transfections

Use the following table to scale the volumes for your transfection experiment. The most common sizes factor are listed below.

Culture Vessel	Multiplication Factor ¹	Shared Reagents		DNA Transfection			siRNA Transfection	
		Vol. Growth Medium	Vol. Opti-MEM for complexing	DNA (μg)	P3000™ (μL)	Lipofectamine® 3000 Reagent ² (μL)	siRNA (pmol)	Lipofectamine® 3000 Reagent ² (μL)
96-well	0.17	100 μL	2 × 5 μL	0.1	0.2	0.15, 0.3	3	0.3
48-well	0.50	250 μL	2 × 12.5 μL	0.25	0.5	0.37, 0.75	7.5	0.75
24-well	1.00	500 μL	2 × 25 μL	0.5	1	0.75, 1.5	15	1.5
12-well	2.00	1 mL	2 × 50 μL	1	2	1.5, 3	30	3
6-well	5.00	2 mL	2 × 125 μL	2.5	5	3.75, 7.5	75	7.5
60 mm	11.05	5 mL	2 × 250 μL	5.5–11	11–22	8.25, 16.5	166	17
10 cm	28.95	10 mL	2 × 500 μL	14–28	28–56	21.7, 43.4	434	43
T75	39.47	15 mL	2 × 750 μL	20–40	40–80	29.6, 59.2	592	59
T175	92.11	35 mL	2 × 1.75 mL	46–90	92–180	69, 138	1382	138

¹ After determining the optimum reagent amount, use the multiplication factor to determine the reagent amount needed for your new plate format.

² Optimum amount needed is determined from the Lipofectamine® 3000 Transfection Reagent Protocol.