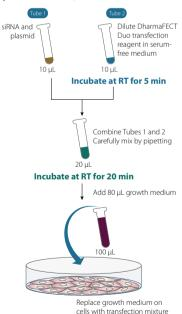


## DharmaFECT™ Duo co-transfection protocol

The following is a protocol for co-transfecting siRNA reagents with a plasmid into cultured mammalian cells using DharmaFECT™ Duo transfection reagent (Cat #T-2010-xx). For more details please see full protocol.

This protocol is written for transfection into 96-well tissue culture plates at 100 nM final concentration of siRNA with 100 ng of plasmid (100 µL final volume).



96-well protocol			
Day 1			
Cell plating	Seed cells at a density that is appropriate for your experiment		
Day 2			
Prepare working solutions of reagents for transfection	siRNA	Dilute siRNA to a working concentration of 2 μM in 1× siRNA buffer or another appropriate RNase-free solution	
	Plasmid	Dilute plasmid to a working concentration of 20 µg/mL in 10 mM Tris-HCl pH 7.4-buffered solution or another appropriate RNase-free solution	
Combine working solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	siRNA (2 μM)	5 μL	_ μL
	Plasmid (20 μg/mL)	5 μL	_ μL
Prepare working solution of DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	0.05-0.5 μL	_ µL
	Serum-free medium	Το 10 μL	_ μL
	Incubate at room temperature for 5 minutes before next step		
Combine transfection mixture	Combine Tube 1 and Tube 2 and carefully mix by pipeting		
	Incubate at room temperature for 20 minutes before next step		
	Add antibiotic-free full growth medium	80 μL	_ μL
	Total	100 μL	_ μL
Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture		

## If you have any questions, contact

- t +44 (0) 1223 976 000 (UK) or +1 800 235 9880 (USA); +1 303 604 9499 (USA)
- **f** + 44 (0) 1223 655 581
- w horizondiscovery.com/contact-us or dharmacon.horizondiscovery.com/service-and-support

Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom