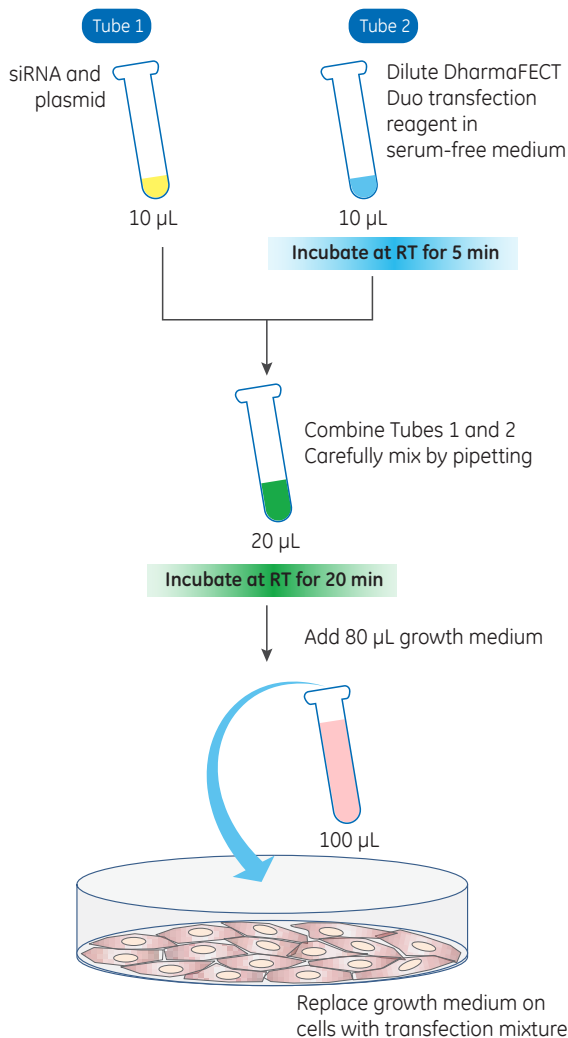


# Dharmacon™ DharmaFECT™ Duo Transfection Reagent—for siRNA and Plasmid Co-transfection protocol

The following is a protocol for co-transfecting Dharmacon™ 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 1x 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
	<b>Tube 1</b>		
	siRNA (2 µM)	5 µL	_ µL
	Plasmid (20 µg/mL)	5 µL	_ µL
Prepare working solution of DharmaFECT Duo for transfection	<b>Tube 2</b>		
	DharmaFECT Duo transfection reagent	0.05–0.5 µL	_ µL
	Serum-free medium	To 10 µL	_ µL
<b>Incubate at room temperature for 5 minutes before next step</b>			
Combine transfection mixture	Combine Tube 1 and Tube 2 and carefully mix by pipetting		
	<b>Incubate at room temperature for 20 minutes before next step</b>		
	Add antibiotic-free full growth medium	80 µL	_ µL
	Total	100 µL	_ µL
Transfect cells	<b>Replace growth medium on cells with 100 µL of transfection mixture</b>		

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon, Inc., a General Electric company doing business as GE Healthcare. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2017 General Electric Company—All rights reserved. First published June 2017. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK.



Orders can be placed at:  
[gelifesciences.com/dharmacon](http://gelifesciences.com/dharmacon)

Customer Support: [cs.dharmacon@ge.com](mailto:cs.dharmacon@ge.com)  
 Technical Support: [ts.dharmacon@ge.com](mailto:ts.dharmacon@ge.com) or  
 +1 800 235 9880; 303 604 9499 if you have any questions.