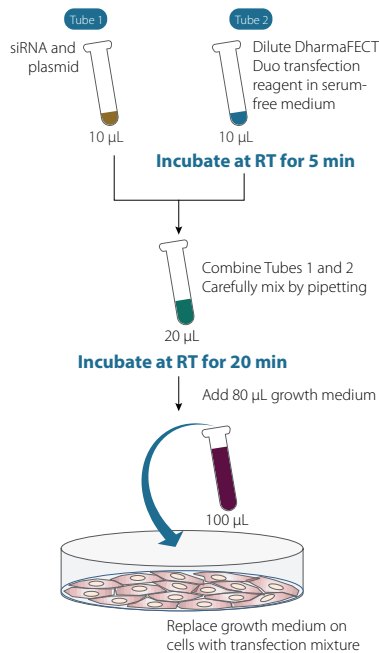


DharmaconTM DharmaFECTTM 1–4 transfection protocol

The following is a protocol for transfecting DharmaconTM synthetic siRNA or microRNA reagents into cultured mammalian cells using DharmaFECTTM transfection reagents: DharmaFECT 1 (Cat #T-2001-xx), DharmaFECT 2 (Cat #T-2002-xx), DharmaFECT 3 (Cat #T-2003-xx), or DharmaFECT 4 (Cat #T-2004-xx). For more details please see full [protocol](#).

This protocol is written for transfection of cells into 96-well tissue culture plates at 25 nM final concentration of siRNA (100 µL final volume).



If you have any questions

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

©2018 Horizon Discovery Group Company—All rights reserved. First published March 2018. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

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 5 µM in 1× siRNA buffer or another appropriate RNase-free solution	
		For one well	For multiple wells
Dilute siRNA for transfection	Tube 1	siRNA (5 µM)	0.5 µL
		Serum-free medium	To 10 µL
Prepare working solution of DharmaFECT for transfection	Tube 2	DharmaFECT transfection reagent	0.05–0.5 µL
		Serum-free medium	To 10 µL
Incubate at room temperature for 5 minutes before next step			
Combine Tube 1 and Tube 2 and carefully mix by pipetting			
Incubate at room temperature for 20 minutes before next step			
Combine transfection mixture	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		