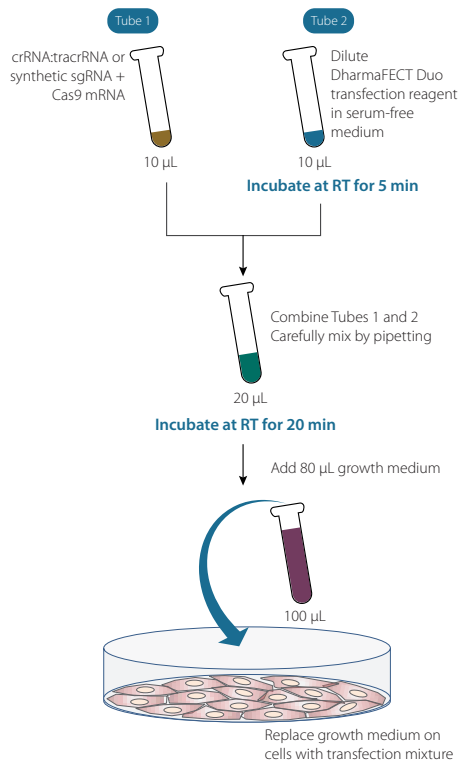


Dharmacon™ Edit-R™ Cas9 mRNA and synthetic guide RNA transfection protocol

The following is an abbreviated protocol for transfecting Dharmacon™ Edit-R™ Cas9 mRNA (Cat #CAS11195, #CAS11859, or #CAS11860) with synthetic guide RNA into cultured mammalian cells using DharmaFECT™ Duo transfection reagent (Cat #T-2010-xx). Synthetic guide RNA can be either Edit-R synthetic tracrRNA (Cat #U-002005-xx) complexed with crRNA (predesigned or custom) or Edit-R synthetic single guide RNA (sgRNA, custom). Intended for use after optimization for your cell line has been completed. For full details, as well as optimization guidelines please see the [Technical Manual](#).



96-well protocol			
Day 1			
Cell plating	Seed cells at a density that gives 70-90% confluency on the next day		
Day 2			
Prepare working solutions of reagents for transfection	synthetic guide RNA	Dilute and mix crRNA and tracrRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4 or Dilute synthetic sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4	
	Cas9 mRNA	Dilute Cas9 mRNA to a working concentration of 100 ng/µL in serum-free medium	
		For one well	For multiple wells
Combine working solutions for transfection mixture	Tube 1		
	synthetic guide RNA	1.25 µL	_ µL
	Cas9 mRNA	2 µL	_ µL
	Serum-free medium	To 10 µL	_ µL
Prepare working solution of DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	0.1–0.8 µL	_ µL
	Serum-free medium	To 10 µL	_ µL
Incubate at room temperature for 5 minutes before next step			
Combine Tube 1 and Tube 2 and carefully mix by pipetting			
Combine transfection mixture	Incubate at room temperature for 20 minutes before next step		
	Add full growth medium	80 µL	_ µL
	Total	100 µL	_ µL
Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture		

24-well protocol			
Day 1			
Cell plating	Seed cells at a density that gives 70-90% confluency on the next day		
Day 2			
Prepare working concentration solutions of materials for transfection	synthetic guide RNA	Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 µM in 10 mM Tris-HCl (pH7.4) or Dilute synthetic sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4	
	Cas9 mRNA	Dilute Edit-R CRISPR-Cas9 mRNA to a working concentration of 100 ng/µL in serum-free medium	
Combine working concentration solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	synthetic guide RNA	5 µL	_ µL
	Cas9 mRNA	10 µL	_ µL
	Serum-free medium	To 50 µL	_ µL
Prepare working concentration DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	0.5–0.8 µL	_ µL
	Serum-free medium	To 50 µL	_ µL
	Incubate at room temperature for 5 minutes before next step		
	Combine Tube 1 and Tube 2 and carefully mix by pipeting		
Prepare transfection mixture	Incubate at room temperature for 20 minutes before next step		
	Add full growth medium	400 µL	_ µL
	Total	500 µL	_ µL
	Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture	

6-well protocol			
Day 1			
Cell plating	Seed cells at a density that gives 70-90% confluency on the next day		
Day 2			
Prepare working concentration solutions of materials for transfection	synthetic guide RNA	Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 µM in 10 mM Tris-HCl (pH7.4) or Dilute synthetic sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4	
	Cas9 mRNA	Dilute Edit-R CRISPR-Cas9 mRNA to a working concentration of 100 ng/µL in serum-free medium	
Combine working concentration solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	synthetic guide RNA	25 µL	_ µL
	Cas9 mRNA	50 µL	_ µL
	Serum-free medium	To 250 µL	_ µL
Prepare working concentration DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	2.5–20 µL	_ µL
	Serum-free medium	To 250 µL	_ µL
	Incubate at room temperature for 5 minutes before next step		
	Combine Tube 1 and Tube 2 and carefully mix by pipeting		
Prepare transfection mixture	Incubate at room temperature for 20 minutes before next step		
	Add full growth medium	2,000 µL	_ µL
	Total	2,500 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

All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published October 2017. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

horizon[™]
INSPIRED CELL SOLUTIONS