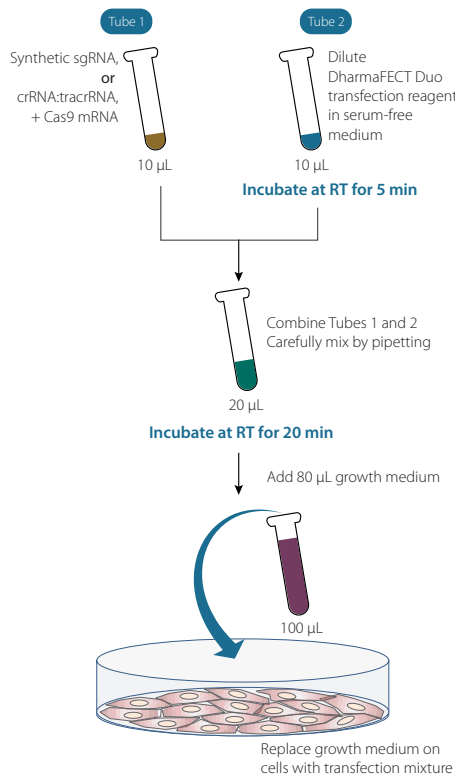


# Edit-R™ Cas9 mRNA and synthetic guide RNA transfection protocol

The following is an abbreviated protocol for transfecting 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 synthetic single guide RNA, or synthetic crRNA complexed with tracrRNA. 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 sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl, pH 7.4 <b>or</b> Dilute and mix crRNA and tracrRNA 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	<b>Tube 1</b>		
	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	<b>Tube 2</b>		
	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	<b>Incubate at room temperature for 20 minutes before next step</b>		
	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 sgRNA to a working concentration of 2 $\mu$ M in 10 mM Tris-HCl, pH 7.4 <b>or</b> Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 $\mu$ M in 10 mM Tris-HCl, pH7.4	
	Cas9 mRNA	Dilute Edit-R CRISPR-Cas9 mRNA to a working concentration of 100 ng/ $\mu$ L in serum-free medium	
Combine working concentration solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	Synthetic guide RNA	5 $\mu$ L	_ $\mu$ L
Cas9 mRNA	10 $\mu$ L	_ $\mu$ L	
Serum-free medium	To 50 $\mu$ L	_ $\mu$ L	
Prepare working concentration DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	0.5–0.8 $\mu$ L	_ $\mu$ L
	Serum-free medium	To 50 $\mu$ L	_ $\mu$ L
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 $\mu$ L	_ $\mu$ L
	Total	500 $\mu$ L	_ $\mu$ L
Transfect cells	Replace growth medium on cells with 100 $\mu$ L 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 sgRNA to a working concentration of 2 $\mu$ M in 10 mM Tris-HCl, pH 7.4 <b>or</b> Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 $\mu$ M in 10 mM Tris-HCl, pH7.4	
	Cas9 mRNA	Dilute Edit-R CRISPR-Cas9 mRNA to a working concentration of 100 ng/ $\mu$ L in serum-free medium	
Combine working concentration solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	Synthetic guide RNA	25 $\mu$ L	_ $\mu$ L
Cas9 mRNA	50 $\mu$ L	_ $\mu$ L	
Serum-free medium	To 250 $\mu$ L	_ $\mu$ L	
Prepare working concentration DharmaFECT Duo for transfection	Tube 2		
	DharmaFECT Duo transfection reagent	2.5–20 $\mu$ L	_ $\mu$ L
	Serum-free medium	To 250 $\mu$ L	_ $\mu$ L
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 $\mu$ L	_ $\mu$ L
	Total	2,500 $\mu$ L	_ $\mu$ L
Transfect cells	Replace growth medium on cells with 100 $\mu$ L 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](http://horizondiscovery.com/contact-us)

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

All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2019 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**<sup>™</sup>  
INSPIRED CELL SOLUTIONS