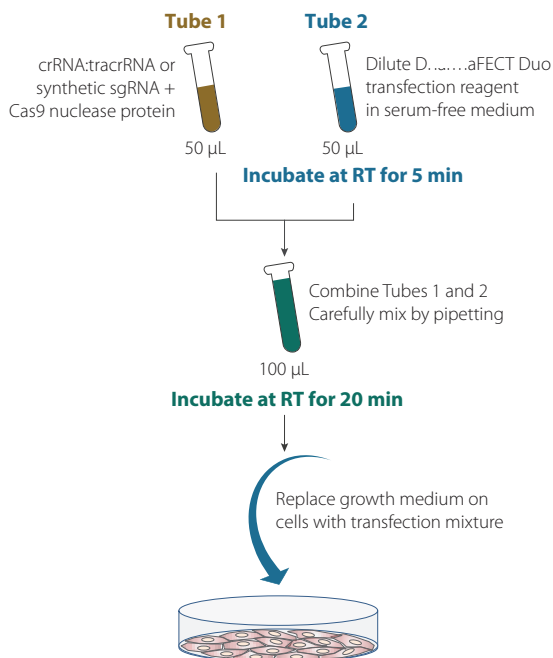


Edit-RTM Cas9 nuclease protein and synthetic guide RNA transfection protocol

The following is a protocol for transfecting Edit-RTM Cas9 Nuclease protein NLS, (Cat #CAS11XXX) with synthetic guide RNA into cultured mammalian cells using DharmaFECTTM transfection reagents (Cat #T-20XX-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 sgRNA (custom). For a more detailed protocol please see the [Technical Manual](#).

The protocol is written for transfection into 96-well tissue culture plates (100 µL final volume).



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 September 2016. 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 gives 70-90% confluency on the next day

Day 2

Prepare working solutions for 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 nuclease protein	Dilute Cas9 nuclease protein to a working concentration of 2.5 µM in serum-free medium	
Combine working solutions for transfection mixture		For one well	For multiple wells
	Tube 1		
	Synthetic guide RNA	2.5 µL	_ µL
Cas9 nuclease protein	1 µL	_ µL	
Serum-free medium	To 50 µL	_ µL	
Prepare working solution of DharmaFECT for transfection	Tube 2		
	DharmaFECT transfection reagent	0.1–0.8 µL	_ µL
Serum-free medium	To 50 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 pipetting		
	Incubate at room temperature for 20 minutes before next step		
Total	100 µL	_ µL	
Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture		
Return to full media	After 14-18 hours; replace transfection mixture on the cells with typical cellular growth medium		

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	crRNA:tracrRNA	Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 μM in 10 mM Tris-HCl (pH7.5)	
	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		
	crRNA:tracrRNA	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			
Prepare transfection mixture	Combine Tube 1 and Tube 2 and carefully mix by pipeting		
	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	crRNA:tracrRNA	Dilute and mix crRNA and tracrRNA to a working concentration of 2.5 μM in 10 mM Tris-HCl (pH7.5)	
	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		
	crRNA:tracrRNA	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			
Prepare transfection mixture	Combine Tube 1 and Tube 2 and carefully mix by pipeting		
	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		

©2017 Horizon Discovery Group Company—All rights reserved. First published September 2016. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom

Orders can be placed at dharmacon.com

Customer Support: cs.dharmacon@horizondiscovery.com

Technical Support: ts.dharmacon@horizondiscovery.com or

If you have any questions: +1 800 235 9880 (option 1); +1 303 604 9499 (option 1)

t +44 (0) 1223 976 000 (UK) or +1 855 772 4252 (USA)

f +44 (0)1223 655 581

e info@horizondiscovery.com

w www.horizondiscovery.com

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

ABgene, ALPS 3000, Easy Pierce, Displace, Fisherbrand and Thermo Scientific are trademarks of Thermo Fisher Scientific, Inc.
©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.

horizonTM
INSPIRED CELL SOLUTIONS

V1-0816