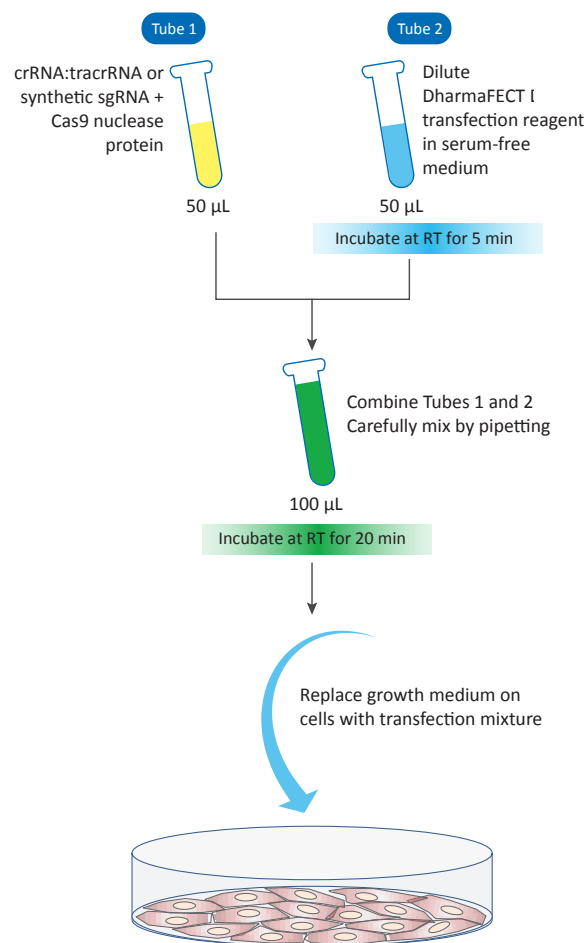


# Dharmacon™ Edit-R™ Cas9 nuclease protein and synthetic guide RNA transfection protocol

The following is a protocol for transfecting Dharmacon™ Edit-R™ Cas9 Nuclease protein NLS, (Cat #CAS11XXX) with synthetic guide RNA into cultured mammalian cells using DharmaFECT™ 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).



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
	<b>Tube 1</b>		
	Synthetic guide RNA	2.5 µL	_ µL
	Cas9 nuclease protein	1 µL	_ µL
Prepare working solution of DharmaFECT for transfection	<b>Tube 2</b>		
	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		

©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

**Dharmacon™**  
A Horizon Discovery Group Company

Orders can be placed at:  
dharmacon.com

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