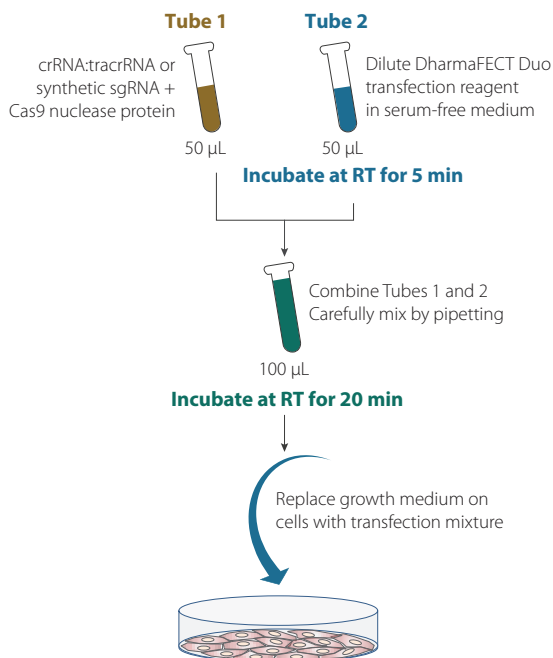


Dharmacon™ Edit-R™ Cas9 nuclease protein and synthetic guide RNA transfection protocol

The following is a protocol for transfecting [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).



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		