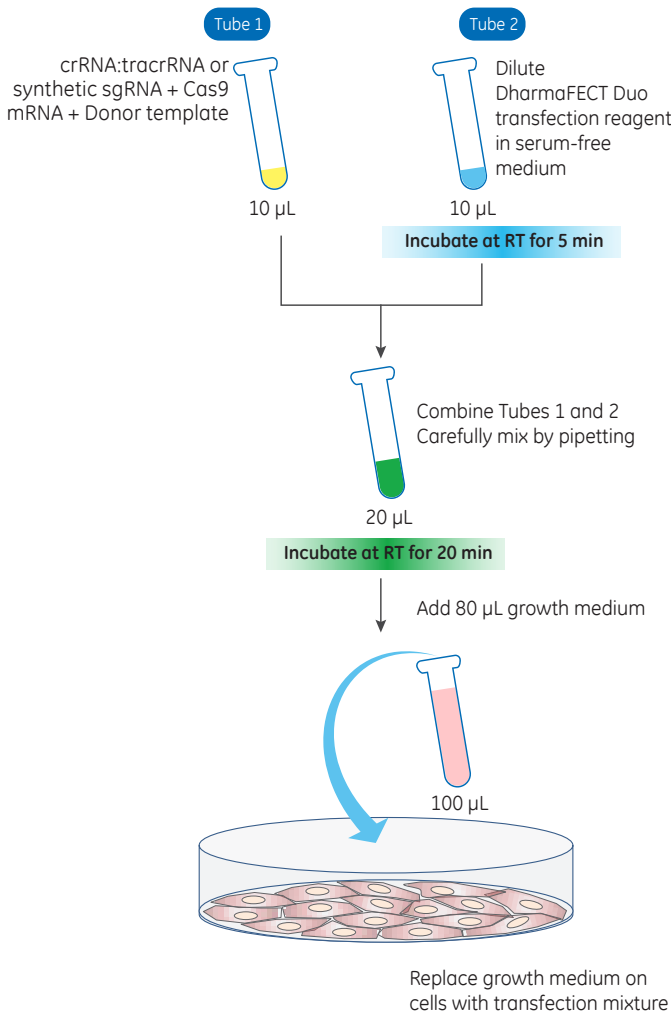


Dharmacon™ Edit-R™ Cas9 mRNA, synthetic guide RNA, and HDR donor template transfection protocol

The following is a protocol for transfecting Dharmacon™ Edit-R™ Cas9 mRNA (Cat #CAS11195) with synthetic guide RNA and homology-directed repair (HDR) donor template (ssDNA oligonucleotide or plasmid) 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). For a more detailed protocol please see this [protocol](#) and [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 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 7.4-7.5 or Dilute synthetic sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4-7.5		
	Donor template	Dilute donor oligo to a working concentration of 1.0 µM in 10 mM Tris-HCl pH 7.4-7.5; or Dilute donor plasmid to a working concentration of 100 ng/µL in 10 mM Tris-HCl pH 7.4-7.5		
	Cas9 mRNA	Dilute Cas9 mRNA to a working concentration of 100 ng/µL in serum-free medium		
Combine working solutions for transfection mixture		For one well	For multiple wells	
	Tube 1			
	synthetic guide RNA	1.25 µL	_ µL	
	Donor template	ssDNA donor oligo	1 µL	_ µL
		Donor plasmid	2 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 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	80 µL	_ µL	
	Total	100 µL	_ µL	
Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture			

GE, imagination at work and GE monogram are trademarks of General Electric Company. Dharmacon, Inc., a General Electric company doing business as GE Healthcare. All other trademarks are the property of General Electric Company or one of its subsidiaries. ©2017 General Electric Company—All rights reserved. First published November 2016. GE Healthcare UK Limited, Amersham Place, Little Chalfont, Buckinghamshire, HP7 9NA, UK.



Orders can be placed at:
gelifsciences.com/dharmacon

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