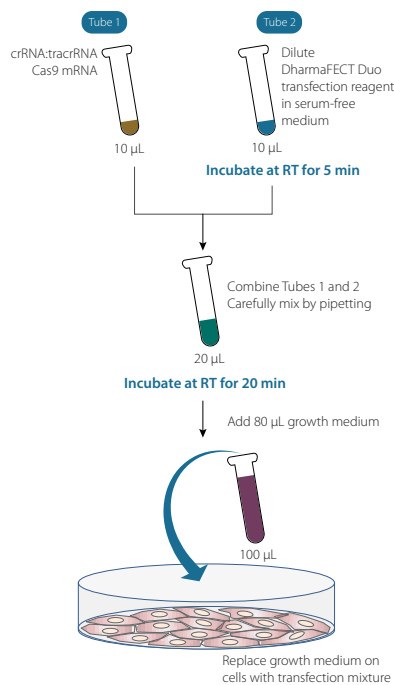


# Edit-R™ Cas9 mRNA, synthetic guide RNA, and HDR donor template transfection protocol

The following is a protocol for transfecting **Edit-R™ Cas9 mRNA** with synthetic guide RNA and homology-directed repair (HDR) donor template (ssDNA oligonucleotide or plasmid) into cultured mammalian cells using **DharmaFECT™ Duo transfection reagent**. Synthetic guide RNA can be either Edit-R synthetic tracrRNA 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](#).



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 <b>or</b> Dilute synthetic sgRNA to a working concentration of 2 µM in 10 mM Tris-HCl pH 7.4	
	Donor template	Dilute donor oligo to a working concentration of 1.0 µM in 10 mM Tris-HCl pH 7.4 <b>or</b> Dilute donor plasmid to a working concentration of 100 ng/µL 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	
Combine working solutions for transfection mixture		For one well	For multiple wells
	<b>Tube 1</b>		
	Synthetic guide RNA	1.25 µL	_ µL
	Donor template	ssDNA donor oligo 1 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	<b>Tube 2</b>		
	DharmaFECT Duo transfection reagent	0.1–0.8 µL	_ µL
	Serum-free medium	To 10 µL	_ µL
<b>Incubate at room temperature for 5 minutes before next step</b>			
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		

## For more information

To find the contact information in your country for your technology of interest, please visit us at [horizondiscovery.com/contact-us](https://horizondiscovery.com/contact-us)

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