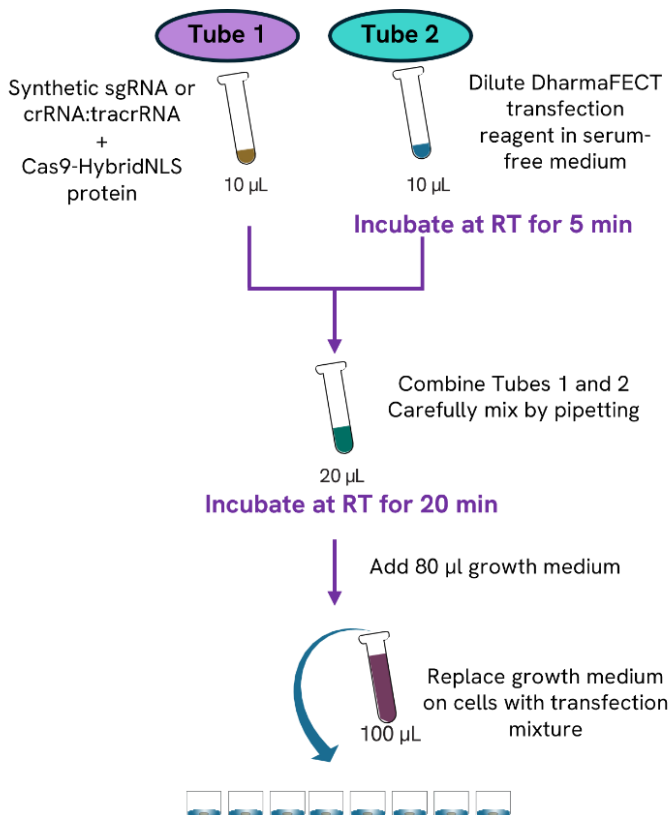


The following is a protocol for transfecting Dharmacon™ Edit-R™ synthetic guide RNA with Cas9 Protein (Cat# CAS135XX or CAS122XX) into cultured mammalian cells using [DharmaFECT™ transfection reagents](#). Synthetic guide RNA can be either [Edit-R predesigned](#) or [custom](#) synthetic single guide RNA (sgRNA), or Edit-R synthetic tracrRNA ([Cat #U-002005-xx](#)) complexed with [Edit-R predesigned](#) or [custom](#) crRNA. This protocol is intended for use after transfection conditions have been optimized for your cell line of interest. For full details, as well as optimization guidelines please see the full [technical manual](#).



96-well protocol				
Day 1	Plate cells	Seed cells at a density that generates 70-90% confluency on the next day		
	Prepare working solutions of reagents for transfection	Synthetic guide RNA	Dilute synthetic sgRNA to a working concentration of 2.5 µM in 10 mM Tris-HCl pH 7.4 or dilute and mix crRNA and tracrRNA to a working concentration of 2.5 µM in 10 mM Tris-HCl pH 7.4	
Day 2	Combine working solutions for transfection mixture	Cas protein	Dilute Cas9 protein to a working concentration of 2.5 µM in serum-free medium	
			For one well	For multiple wells
		Tube 1		
		Synthetic guide RNA	1 µl	— µl
	Prepare working solution of DharmaFECT for transfection	Cas9 protein	1 µl	— µl
		Serum-free medium	To 10 µl	— µl
		Incubate at room temperature for 5 minutes before next step		
	Combine transfection mixture	Tube 2		
		DharmaFECT transfection reagent	0.1-0.8 µl	— µl
		Serum-free medium	To 10 µl	— µl
		Incubate at room temperature for 20 minutes before next step		
	Transfect cells	Add full growth medium	80 µl	— µl
		Total volume	100 µl	— µl
	Replace growth medium on cells with 100 µl of transfection mixture			

If you have any questions, please visit us at horizondiscovery.com/contact-us
Horizon Discovery, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

Horizon Discovery Biosciences LTD is a Revvity company
Dharmacon Inc is a Revvity company