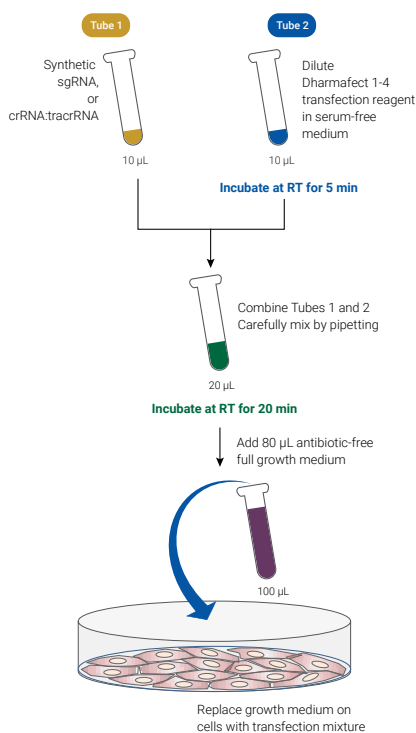


CRISPRmod synthetic guide RNA transfection protocol for dCas9-VPR or dCas9-SALL1-SDS3 expressing cells

The following is an abbreviated protocol for transfecting CRISPRmod CRISPRa/CRISPRi synthetic guide RNA into cultured mammalian cells expressing dCas9-VPR or dCas9-SALL1-SDS3 using DharmaFECT™ 1-4 transfection reagent (Cat. T-2001, T-2002, T-2003, T-2004). Synthetic guide RNA can be either synthetic single guide RNA, or synthetic crRNA complexed with tracrRNA. Intended for use after optimization for your cell line has been completed. For full details, as well as optimization guidelines please see the appropriate CRISPRmod technical manual [CRISPRa tech manual](#) and [CRISPRi technical manual](#)).

This protocol is written for transfection of dCAS9 expressing cells into 96, 24, or 6-well tissue culture plates at 25 nM final concentration of synthetic guide RNA.



96-well protocol

Day 1

Cell plating Seed cells at a density that is optimal for specific downstream phenotypic assay(s)

Day 2

Prepare working solutions of reagents for transfection Synthetic guide RNA Dilute sgRNA to a working concentration of 1 µM in 10 mM Tris-HCl, pH 7.4 **or** Dilute and mix crRNA and tracrRNA to a working concentration of 1 µM in 10 mM Tris-HCl, pH 7.4

	For one well		For multiple wells	
	Tube 1			
Combine working solutions for transfection mixture	Synthetic guide RNA	2.5 µL	_ µL	_ µL
	Serum-free medium	To 10 µL	_ µL	_ µL
Prepare working solution of Dharmafect 1-4 for transfection	DharmaFECT 1-4 transfection reagent	0.05-0.8 µL	_ µL	_ µL
	Serum-free medium	To 10 µL	_ µL	_ µL
Incubate at room temperature for 5 minutes before next step				
Combine Tube 1 and Tube 2 and carefully mix by pipeting				
Incubate at room temperature for 20 minutes before next step				
Combine transfection mixture	Add full growth medium	80 µL	_ µL	_ µL
	Total	100 µL	_ µL	_ µL
Transfect cells	Replace growth medium on cells with 100 µL of transfection mixture			
Incubate cells for 48-72 hours before performing downstream phenotypic assay(s) or gene expression analysis.				

24-well protocol

Day 1

Cell plating Seed cells at a density that is optimal for specific downstream phenotypic assay(s)

Day 2

Prepare working solutions of reagents for transfection Synthetic guide RNA Dilute sgRNA to a working concentration of 1 μ M in 10 mM Tris-HCl, pH 7.4 **or** Dilute and mix crRNA and tracrRNA to a working concentration of 1 μ M in 10 mM Tris-HCl, pH7.4

		For one well		For multiple wells	
		Tube 1			
Combine working solutions for transfection mixture	Synthetic guide RNA	12.5 μ L			– μ L
	Serum-free medium	To 50 μ L			– μ L

Tube 2					
Prepare working solution of Dharmafect 1-4 for transfection	DharmaFECT 1-4	0.24 - 4 μ L			– μ L
	Serum-free medium	To 50 μ L			– μ L

Incubate at room temperature for 5 minutes before next step

Combine Tube 1 and Tube 2 and carefully mix by pipeting

Incubate at room temperature for 20 minutes before next step					
Combine transfection mixture	Add full growth medium	400 μ L			– μ L
	Total	500 μ L			– μ L

Transfect cells Replace growth medium on cells with 500 μ L of transfection mixture

Incubate cells for 48-72 hours before performing downstream phenotypic assay(s) or gene expression analysis.

6-well protocol

Day 1

Cell plating Seed cells at a density that is optimal for specific downstream phenotypic assay(s)

Day 2

Prepare working solutions of reagents for transfection Synthetic guide RNA Dilute sgRNA to a working concentration of 1 μ M in 10 mM Tris-HCl, pH 7.4 **or** Dilute and mix crRNA and tracrRNA to a working concentration of 1 μ M in 10 mM Tris-HCl, pH7.4

		For one well		For multiple wells	
		Tube 1			
Combine working solutions for transfection mixture	Synthetic guide RNA	50 μ L			– μ L
	Serum-free medium	To 200 μ L			– μ L

Tube 2					
Prepare working solution of Dharmafect 1-4 for transfection	DharmaFECT 1-4	1-20 μ L			– μ L
	Serum-free medium	To 250 μ L			– μ L

Incubate at room temperature for 5 minutes before next step

Combine Tube 1 and Tube 2 and carefully mix by pipeting

Incubate at room temperature for 20 minutes before next step					
Combine transfection mixture	Add full growth medium	1600 μ L			– μ L
	Total	2000 μ L			– μ L

Transfect cells Replace growth medium on cells with 2000 μ L of transfection mixture

Incubate cells for 48-72 hours before performing downstream phenotypic assay(s) or gene expression analysis.

For more information

To find the contact information in your country for your technology of interest, please visit us at horizondiscovery.com/contact-us

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

©2022 Horizon Discovery Group Ltd and its affiliated companies. All rights reserved. PerkinElmer is a trademark of PerkinElmer, Inc., registered in the United States and other countries. Horizon Discovery is a trademark mark of Horizon Discovery Ltd. Dharmacon is a trademark of Dharmacon, Inc. All PerkinElmer, Inc., Horizon Discovery Ltd and Dharmacon, Inc. trademarks are used with permission. Other product names and brand names may be trademarks or registered trademarks of their respective owners.

horizon
a PerkinElmer company