



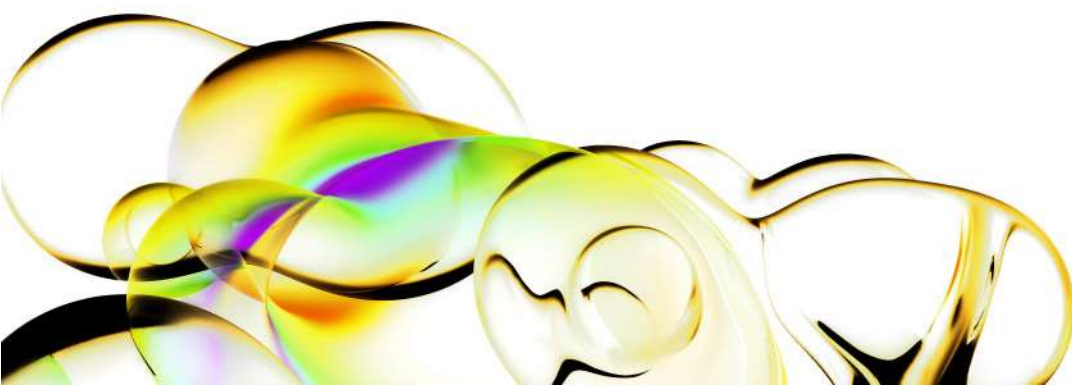
Protocol

Dharmacon™ E-dit-R™ guide RNA control kits via T7EI assay

Version: 1

May 2024

For research only. Not for use in diagnostic procedures.



Product Description

The Edit-R control kits provide validated guide RNAs and primer pairs for detection of cleavage for the designated housekeeping gene or safe harbor region.

These gene-specific positive control kits and region-specific cutting control kits are validated for mismatch detection assays to verify gene editing experiments. The following protocol is an example of thermal cycling and cleavage conditions for a DNA mismatch detection assay using T7 Endonuclease I (T7EI) and the Edit-R control primer set utilizing genomic DNA (gDNA) from direct cell lysis. The same primer pairs can also be used for “Tracking of Indels by Decomposition” (TIDE) analysis. The volumes in this protocol are for direct cell lysis in a 96-well plate format.

T7EI Protocol

1. Lyse cells in 100 µL of 1x Phusion High-Fidelity buffer with additives.

Reagent	Volume	Final concentration
5x Phusion HF Buffer	20 µL	1x
Proteinase K (~ 20 mg/mL)	5 µL	~ 1 mg/mL
RNase A (10 mg/mL)	5 µL	0.5 mg/mL
Water, nuclease free	70 µL	N/A

2. Seal the 96-well tissue culture plate with a plate seal to minimize evaporation and cross contamination. Incubate for 15-30 minutes at 56 °C.
3. Transfer cell lysate into a 96-well PCR-plate, seal plate with a plate seal, and place into a thermocycler for 5 minutes at 96 °C. Briefly centrifuge the plate to collect liquid at the bottom of each well.

Note: Due to variable depth of thermocycler heat blocks, deactivation time may need to be increased to 15 or 30 minutes.

4. Set up 50 µL PCR reaction for each sample to be analyzed.

Note: Other PCR reagents can be used provided that PCR optimization is performed for mismatch detection conditions and testing compatibility with direct cell lysis or purified gDNA.

Reagent	Volume	Final Concentration
5x Phusion HF buffer	10 µL	1x
Forward primer (50 µM)	0.5 µL	500 nM
Reverse primer (50 µM)	0.5 µL	500 nM
dNTPs (10 mM)	1 µL	200 µM each
Phusion Hot Start II High-Fidelity DNA Polymerase (2 U/µL)	1 µL	0.04 U/µL
Water, nuclease free	32 µL	N/A
Cell lysate	5 µL	N/A

5. Thermal cycling conditions:

Cycle step	Temperature	Time	Cycles
Initial denaturation	98 °C	3 min	1
Denature	98 °C	10 s	
Touchdown annealing	72 °C – 1 °C/cycle	15 s	10
Extension	72 °C	30 s	
Denature	98 °C	10 s	
Annealing	62 °C	15 s	25
Extension	72 °C	30 s	
Final extension	72 °C	10 min	1

6. Heat PCR samples to 95 °C for 10 minutes and then slowly (>15 min) cool to room temperature (~25 °C).

7. PCR product (~5 µl) can be run on a 2% agarose gel to confirm amplification.

8. Set up 15 µl reactions for mismatch detection assay using T7EI. Be sure to include at least one sample without the T7EI enzyme (No enzyme negative control and/or untreated (no gene editing negative control)).

Note: Other mismatch detection assays can be used following the manufacturer's recommended protocol.

9. Incubate for 25 minutes at 37 °C.

10. Immediately run entire reaction volume with appropriate gel loading buffer on 2% agarose gel, since the T7EI enzyme cannot be heat inactivated. Compare samples with expected gene editing to the no enzyme and/or untreated negative controls.

Materials

- Thermo Scientific Phusion Hot Start II High-Fidelity DNA Polymerase, 2 U/µL (Cat #F-549S)
- T7 Endonuclease I, 10 U/µL (NEB, Cat #M0302S)
- Thermo Scientific Proteinase K, ~ 20 mg/mL (Cat #EO0492)
- Thermo Scientific RNase A, 10 mg/mL (Cat #EN0531)
- Thermo Scientific 5x Phusion HF Buffer (Cat #F-518L)
- NEBuffer 2 (NEB, Cat #B7002S)

Reagent	Volume	Final concentration
PCR from gDNA (300-500 ng/reaction)	10 µL	variable
Water, nuclease free	3 µL	N/A
NEBuffer 2 (10x)	1.5 µL	1x
T7 Endonuclease I (10 U/µL)	0.5 µL	0.33 U/µL

Target gene/ region	Species	Guide RNA target sequence	Primer	Primer sequence	Tm* (°C)	Expected band sizes (bp)	
						No editing	With editing
<i>PP1B</i>	Human	GUGUAUUUUGACCUACGAAU	Forward	GAAGCTTAGGCTCCGCTCCTT	64.0	505	~ 330
			Reverse	CTCTGCAGGTCAGTTTGCTG	64.3		~ 174
<i>DNMT3B</i>	Human	GCUGAAUUACUCACGCCCA	Forward	TGAGAAGGAGCCACTTGCTT	64.4	544	~ 335
			Reverse	GACCAAGAACGGGAAAGTCA	64.4		~ 209
<i>AASV1_sg1</i>	Human	GACGCAAGGAGACATCCGT	Forward	ACTAGCTGAGCTCTCGGACC	63.1	481	~272
			Reverse	GCTGCCCAAGGATGCTCTTT	68.0		~209
<i>AASV1_sg2</i>	Human	GCGGCTCCAATTCGGAAGTG	Forward	GGCCAGGCAGATAGACCAGA	66.9	514	~303
			Reverse	CCAACGCCGACGGTATCAG	69.1		~211
<i>AASV1_sg3</i>	Human	GAGTTCGGCGGCACAGCAA	Forward	GAACCCAGCGAGTGAAGACG	67.6	590	~427
			Reverse	CTCTCCGATGTTGAGCCCC	68.4		~163
<i>Ppib</i>	Mouse	GUUACUUUGAUUUACAAU	Forward	CCAGCCTGTTTGTGGATT	64.2	572	~409
			Reverse	ACTGTGAGGCTGCCAGATT	64.2		~163
<i>Dnmt3b</i>	Mouse	GCUGAAUUUACCCGCCCA	Forward	ACTTGGTGATTGGTGAAGC	64.3	561	~372
			Reverse	TTAAGCCACACCCTCTCTG	64.5		~189

*Tm calculated specifically for use with Phusion DNA polymerases. Tm should be calculated for specific DNA polymerases and PCR reagents using appropriate parameters.

For more information

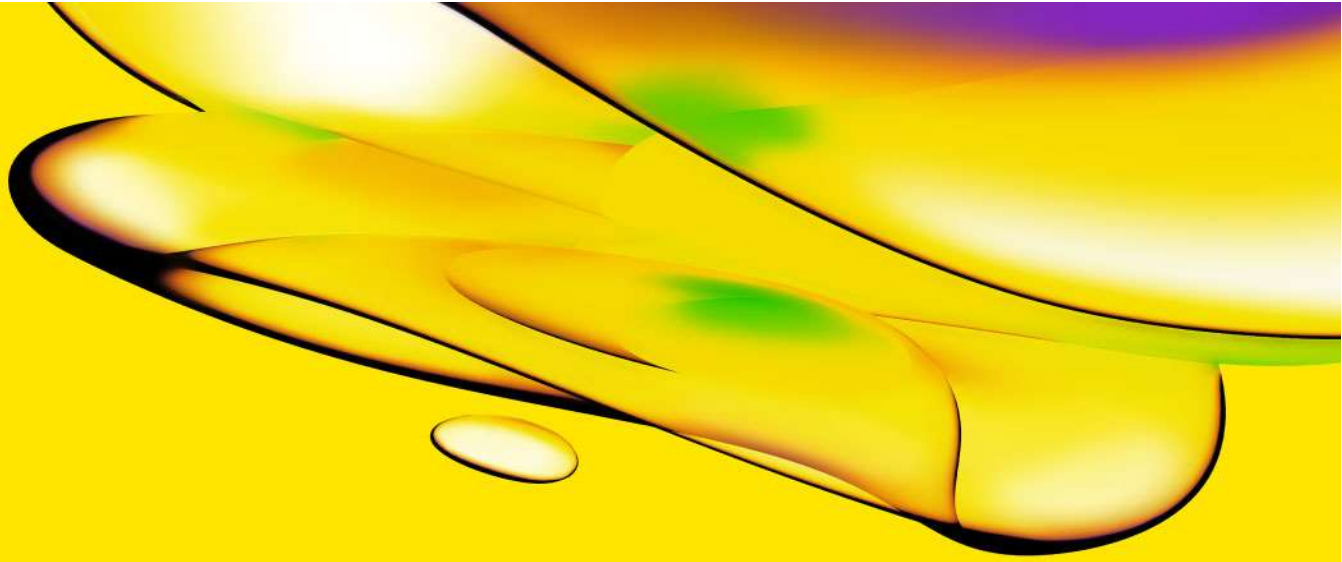
To find the contact information in your country for your technology of interest, please visit us at

horizondiscovery.com/contact-us

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

©2020 The Horizon logo and other trademarks are the property of Horizon Discovery Limited, unless otherwise stated.

DHARMACON is a trademark of Dharmacon Inc.



revvity

Revvity
940 Winter Street
Waltham, MA 02451 USA

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity. All rights reserved.