

# Dharmacon™ Edit-R™ synthetic gRNA positive control kits

## Product description

The Edit-R Synthetic Positive Control kits provide validated crRNAs and primer pairs for detection of cleavage for the designated human, mouse, or rat positive control gene.

These gene-specific positive controls and 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 crRNA Control primer set utilizing genomic DNA (gDNA) from direct cell lysis. For direct cell lysis, cells were treated for gene editing in a 96-well format.

### 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

- Seal the 96-well plate with a plate seal to minimize evaporation and cross contamination. Incubate for 15-30 minutes at 56 °C, followed by deactivation for 5 minutes at 96 °C. Briefly centrifuge plate to collect liquid at bottom of wells.

- Set up 50 µL PCR 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

### 4. Thermal cycling condition

Cycle step	Temperature	Time	Cycle(s)
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

- Heat PCR samples to 95 °C for 10 minutes and then slowly (> 15 minutes) cool to room temperature (~ 25 °C).
- PCR product (~ 5 µL) can be run on 2% agarose gel to confirm amplification.

7. Set up 15  $\mu$ L 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.**

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

8. Incubate for 25 minutes at 37 °C.
9. Immediately run entire reaction volume with appropriate gel loading buffer on 2% agarose gel, since 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/ $\mu$ L (Cat #F-549S)
- T7 Endonuclease I, 10 U/ $\mu$ L (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)

Species	Gene target	Species	Primer	Sequence	Tm* (°C)	MW	Extinction coefficient	Expected band sizes (bp)	
								No editing	With editing
UK-007050-01-xx	<i>PPIB</i>	Human	Forward	GAACCTAGGCTCCGCTCCTT	64.0	6044.0	176184	505	~ 330,
			Reverse	CTCTGCAGGTCAGTTTGCTG	64.3	6115.0	178151		~ 174
UK-007150-01-xx	<i>Ppib</i>	Mouse	Forward	CCAGCCTGTTTGTGGATT	64.2	6105.1	176905	572	~ 409,
			Reverse	ACTGTGAGGCTGCCAGATT	64.2	6148.1	191297		~ 163
UK-007250-01-xx	<i>Ppib</i>	Rat	Forward	TGCCATTTCCTCACTCCCTA	60.4	5938.9	165727	524	~ 342,
			Reverse	GGCTACCGCTACAAATGGAC	62.4	6111.0	197820		~ 170
UK-007060-01-xx	<i>DNMT3B</i>	Human	Forward	TGAGAAGGAGCCACTTGCTT	64.4	6157.1	197197	544	~ 335,
			Reverse	GACCAAGAACGGGAAAGTCA	64.4	6193.1	220800		~ 209
UK-007160-01-xx	<i>Dnmt3b</i>	Mouse	Forward	ACTTGGTGATTGGTGGAAAGC	64.3	6228.1	199164	561	~ 372,
			Reverse	TTAAGCCACCCCTCCTCTG	64.5	5972.9	174218		~ 189
UK-007260-01-xx	<i>Dnmt3b</i>	Rat	Forward	GGCATTGATCCCATGTAACC	60.4	6077.0	189330	532	~ 319
			Reverse	CTCCGAGCTGCTAAGAGGAA	62.4	6151.0	201754		~ 212

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

### If you have any questions, contact

**t** +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

**f** + 44 (0)1223 655 581

**w** [horizondiscovery.com/contact-us](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://dharmacon.horizondiscovery.com/service-and-support)

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

Phusion Hot Start II High-Fidelity, Proteinase K, RNase A, and 5x Phusion HF Buffer is a registered trademark of Thermo Scientific, Inc. T7 Endonuclease I and NEBuffer 2 is a registered trademark of NEB. All trademarks are the property of Horizon Discovery Company unless otherwise specified. ©2018 Horizon Discovery Group Company—All rights reserved. First published March 2015. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

**horizon**<sup>TM</sup>  
INSPIRED CELL SOLUTIONS

V30818