

Dharmacon™ Edit-R™ HDR donor plasmid kits

Product description

The Dharmacon™ Edit-R™ HDR donor plasmid kit contains all the DNA components to build and quality check a custom DNA donor plasmid to introduce a fluorescent tag or custom sequence to any genomic location when coupled with Edit-R HDR DNA donor plasmid homology arm (HA) custom primer pairs. Specifically, the EGFP and mKate2 kits come with the Edit-R HDR donor backbone, the appropriate donor plasmid insert, and the Edit-R HDR donor plasmid colony PCR primers specific to the backbone and insert. The Universal kit for custom inserts differs in that it lacks the insert specific colony PCR primers and the insert fragment. Reagents are supplied as dried pelleted oligos with enough material for ten (10) reactions, when used as directed.

The HA PCR primers are designed using the free, online [HDR Donor Designer](#) to amplify HAs with lengths ranging from 500 to 1000 bp; additionally, the primers will have an adapter sequence appended, which is necessary for compatibility with the HDR donor plasmid backbone and mKate2. The HDR Donor Designer additionally designs these primers to introduce silent mutations to disrupt the intact CRISPR sequence to avoid unintentional cutting of gDNA previously edited to contain the fluorescent reporter.

Kit contents

Cat. #	Description	Item #	Amount
UK-008100-01-10	Edit-R HDR Plasmid Donor Kit - EGFP		
	Edit-R HDR Plasmid Donor Backbone	U-008300-01-1200	1200 ng
	Edit-R HDR insert - EGFP	U-008100-01-250	250 ng
	Edit-R Colony PCR Primer EGFP Forward	U-008100-FB-05	5 nmol
	Edit-R Colony PCR Primer EGFP Reverse	U-008100-RA-05	5 nmol
	Edit-R Colony PCR Primer Backbone Forward	U-008300-FW-05	5 nmol
UK-008200-03-10	Edit-R HDR Plasmid Donor Kit -mKate2		
	Edit-R HDR Plasmid Donor Backbone	U-008300-01-1200	1200 ng
	Edit-R HDR insert - mKate2	U-008200-01-250	250 ng
	Edit-R Colony PCR Primer Backbone Reverse	U-008300-RV-05	5 nmol
	Edit-R Colony PCR Primer Backbone Forward	U-008300-FW-05	5 nmol
	Edit-R Colony PCR Primer mKate2 Reverse	U-008200-RA-05	5 nmol

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 **or** dharmacon.horizondiscovery.com/service-and-support

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

©2018 Horizon Discovery Group Company—All rights reserved. All trademarks are the property of Horizon Discovery Company unless otherwise specified. First published June 2014. UK Registered Head Office: Building 8100, Cambridge Research Park, Cambridge, CB25 9TL, United Kingdom.

Cat. #	Description	Item #	Amount
	Edit-R Colony PCR Primer mKate2 Forward	U-008200-FB-05	5 nmol
	Edit-R Colony PCR Primer mKate2 Reverse	U-008200-RA-05	5 nmol
	Edit-R Colony PCR Primer Backbone Forward	U-008300-FW-05	5 nmol
	Edit-R Colony PCR Primer Backbone Reverse	U-008300-RV-05	5 nmol
UK-008300-01-10	Edit-R HDR Plasmid Donor Kit - Universal		
	Edit-R HDR Plasmid Donor Backbone	U-008300-01-1200	1200 ng
	Edit-R Colony PCR Primer Backbone Forward	U-008300-FW-05	5 nmol
	Edit-R Colony PCR Primer Backbone Reverse	U-008300-RV-05	5 nmol

Shipping and storage

Store at -20 °C. The Edit-R HDR Donor Plasmid Kit is shipped at ambient temperature and should be stored in a -20° C freezer.

Handling precautions

Oligonucleotides and purified DNA fragments are susceptible to enzymatic degradation by nucleases and to chemical degradation by extreme pH and temperature. We recommend wearing gloves and maintaining nuclease-free conditions when handling.

Related products

- Dharmacon™ DharmaFECT™ Duo Transfection Reagent is optimized for transfecting plasmid DNA and small RNA into a wide variety of cell lines. For more information, click [here](#).

Related documents

Go to dharmacon.horizondiscovery.com to find:

- [Edit-R HDR Plasmid Donor Kit Technical Manual](#)
- [Application Notes](#)
- [Posters](#)