

RTF optimization plates

Introduction

Optimizing conditions is critical to the success of siRNA screens. In support of screens using RTF siRNA Libraries, we offer RTF Optimization Plates with either siGENOME™ or ON-TARGET*plus*™ Control siRNA Reagents.

These products are available in standard clear cell culture plates, as well as white or black-walled plates for assays with luminescent or fluorescent readouts.

Catalog numbers and descriptions of available products are listed in Table 1. RTF Optimization plates are provided in triplicate with human or mouse product line-specific controls (see Table 2). All controls are plated at 6.25 pmol per well, ready for resuspension and use at 50 nM concentration (see [RTF Optimization Plate Protocol](#) for details).

Table 1. RTF optimization plates: available products

Cat. #	Description	Plate color
H-102200	RTF Optimization Plates— ON-TARGET <i>plus</i> Human	Clear
H-102200B	RTF Optimization Plates (black)— ON-TARGET <i>plus</i> Human	Black
H-102200W	RTF Optimization Plates (white)— ON-TARGET <i>plus</i> Human	White
H-112200	RTF Optimization Plates— ON-TARGET <i>plus</i> Mouse	Clear
H-112200B	RTF Optimization Plates (black)— ON-TARGET <i>plus</i> Mouse	Black
H-112200W	RTF Optimization Plates (white)— ON-TARGET <i>plus</i> Mouse	White
H-002200	RTF Optimization Plates— siGENOME Human/Mouse	Clear
H-002200B	RTF Optimization Plates (black)— siGENOME Human/Mouse	Black
H-002200W	RTF Optimization Plates (white)— siGENOME Human/Mouse	White

Table 2. RTF optimization plate contents

ON-TARGET ^{plus} RTF optimization plates—human		
Row A	ON-TARGET ^{plus} Non-targeting siRNA #1	D-001810-01
Row B	ON-TARGET ^{plus} Non-targeting siRNA #2	D-001810-02
Row C	ON-TARGET ^{plus} Non-targeting siRNA #3	D-001810-03
Row D	ON-TARGET ^{plus} Non-targeting siRNA #4	D-001810-04
Row E	ON-TARGET ^{plus} Non-targeting pool	D-001810-10
Row F	ON-TARGET ^{plus} Human Cyclophilin B control pool	D-001820-10
ON-TARGET ^{plus} RTF optimization plates—mouse		
Row A	ON-TARGET ^{plus} Non-targeting siRNA #1	D-001810-01
Row B	ON-TARGET ^{plus} Non-targeting siRNA #2	D-001810-02
Row C	ON-TARGET ^{plus} Non-targeting siRNA #3	D-001810-03
Row D	ON-TARGET ^{plus} Non-targeting siRNA #4	D-001810-04
Row E	ON-TARGET ^{plus} Non-targeting pool	D-001810-10
Row F	ON-TARGET ^{plus} Mouse Cyclophilin B control pool	D-001820-20
siGENOME RTF optimization plates—human/mouse		
Row A	siGENOME Non-targeting siRNA #2	D-001210-02
Row B	siGENOME Non-targeting siRNA #3	D-001210-03
Row C	siGENOME Non-targeting siRNA #4	D-001210-04
Row D	siGENOME Non-targeting siRNA #5	D-001210-05
Row E	siGENOME Non-targeting pool #2	D-001206-14
Row F	siGENOME Cyclophilin B control pool (H, M, R) Control siRNA	D-001136-01

Shipping and storage

- RTF Optimization Plate reagents are shipped dried-down at ambient temperature in sealed 96-well plates. Under these conditions, the reagents are stable for at least four weeks.
 - » Clear plates (standard): Nunc™ (Thermo Scientific Cat. #152039); Fisher Scientific (Cat. #12-565-901)
 - » White plates: Nunc™ (Thermo Scientific Cat. #152028); Fisher Scientific (Cat. #12-565-903)
 - » Black plates: Nunc™ (Thermo Scientific Cat. #152037); Fisher Scientific (Cat. #2-565-902)
 - » Peelable Aluminum Seal (Agilent Cat #24210-001, white) **OR** Easy Pierce™ Heat Seal (Thermo Scientific Cat. #AB-3738, silver)
- Upon receipt, RTF library plates should be stored at +4 °C. Freezing is not recommended.

Related products

Product	Description
siGENOME and ON-TARGET ^{plus} RTF siRNA Libraries	<ul style="list-style-type: none"> Identify specific genes of interest from entire pathways and gene families Convenient high throughput and rapid screening technique by the effective delivery of siRNA to mammalian cells by reverse transfection SMARTpool siRNAs in a 96-well format Provided as six replicates Custom gene collections available upon request
DharmaFECT siRNA Transfection Reagents	<ul style="list-style-type: none"> siRNA transfection reagents available in four unique lipid-based formulations Optimized for efficient delivery of siRNA into mammalian cells with minimal impact on cell viability Catalog numbers: T-2001-0x, T-2002-0x, T-2003-0x, T-2004-0x

If you have any questions, contact

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 The Horizon logo and other trademarks are the property of Horizon Discovery Limited, unless otherwise stated. DHARMACON is a trademark of Dharmacon Inc.

horizon
a PerkinElmer company