

DharmaFECT™ Duo Transfection Reagent—for siRNA and plasmid co-transfection

Introduction

DharmaFECT™ Duo is a lipid-based reagent specially formulated for co-transfection of plasmid and siRNA. Under optimized conditions, efficient delivery of both plasmid and siRNA can be achieved with minimal cell toxicity. As is generally observed with plasmid transfection, cell viability is reduced compared to siRNA transfection alone. When evaluating delivery conditions, a threshold of 70% cell viability and 80% gene silencing is set for successful, non-toxic co-transfection of plasmid and siRNA.

Product description

- Provided as a liquid at a concentration of 1 mg/mL.
- Confirmed to be free of microbial contaminants and ribonucleases.

Product	Size	Cat. #
DharmaFECT Duo	0.2 mL	T-2010-01
	0.75 mL	T-2010-02
	1.5 mL	T-2010-03
	1.5 mL × 5 tubes	T-2010-04

Shipping and storage

- DharmaFECT Duo co-transfection reagent is shipped on ice packs. If the ice melts during transit there is no risk to the product. Stability testing demonstrates no loss in product functionality after prolonged storage at warm temperatures (see the Stability Testing Product Bulletin for additional details).
- Upon receipt, store at 4 °C. **DO NOT FREEZE.** DharmaFECT transfection reagents are labeled with a date of manufacture and are good for 24 months following that date when stored as recommended.

Related products

We recommend including a positive and negative control, such as Dharmacon™ RNAi Control Reagents, in every RNAi experiment. For more information, click [here](#).

Accompanying documents

siRNA/Plasmid Co-Transfection Protocol

Publication reference guide

When referencing the use of DharmaFECT Duo Co-Transfection Reagents, please include the following information: DharmaFECT Duo Transfection Reagent, Horizon Discovery, Lafayette, CO.

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.