

DharmaFECT™ siRNA transfection reagents

Product description

- Panel of four transfection reagents optimized for efficient delivery of siRNA into mammalian cells.
- Provided as a liquid at a concentration of 1 mg/mL.
- Functionally verified to produce at least 75% transfection efficiency by assessing mRNA knockdown of positive control siRNAs.
- Confirmed to be free of microbial contaminants and ribonucleases.

Shipping and storage

- DharmaFECT siRNA Transfection Reagents are 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 <u>Stability Testing Product Bulletin</u> 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

Basic siRNA Transfection Protocol.

Publication reference guide

When referencing the use of DharmaFECT siRNA transfection reagents, please include the following information: DharmaFECT (1, 2, 3, or 4) siRNA Transfection Reagent, Horizon Discovery Group Company, Lafayette, CO.

Product	Size	Cat.#
DharmaFECT 1	0.2 mL 0.75 mL 1.5 mL 1.5 mL × 5 tubes	T-2001-01 T-2001-02 T-2001-03 T-2001-04
DharmaFECT 2	0.2 mL 0.75 mL 1.5 mL 1.5 mL × 5 tubes	T-2002-01 T-2002-02 T-2002-03 T-2002-04
DharmaFECT 3	0.2 mL 0.75 mL 1.5 mL 1.5 mL × 5 tubes	T-2003-01 T-2003-02 T-2003-03 T-2003-04
DharmaFECT 4	0.2 mL 0.75 mL 1.5 mL 1.5 mL × 5 tubes	T-2004-01 T-2004-02 T-2004-03 T-2004-04
DharmaFECT 1, 2, 3, and 4	0.2 mL each 0.75 mL each 1.5 mL each	T-2005-01 T-2005-02 T-2005-03

If you have any questions

- 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

