

TOX transfection control

Product information

Molecular weight	Sequence	Cat. #
Proprietary	Proprietary	D-001500-01-XX

XX=05,20, or 50 for 5, 20, and 50 nmol amounts

The Dharmacon TOX transfection control is recommended as a first indicator of the success of transfection. It is also an ideal control for optimizing transfection conditions, as it provides a quantifiable measure of cellular uptake by the easily visualized characteristics of cell death or apoptosis.

Shipping and storage

This reagent is shipped and stored under the same conditions as our siRNA controls. siRNA reagents are shipped as dry pellets at room temperature (23 °C). Under these conditions, they are stable for at least four weeks.

- Upon receipt, siRNA reagents should be stored at –20 °C to –80 °C. Under these conditions, they are stable for at least one year.
- Upon resuspension in 1x siRNA Buffer, aliquot the siRNA into small volumes and store at –20 °C to –80 °C. For best results, limit freeze-thawing of each tube to five events. Under these conditions, the siRNA is stable for at least six months.

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.

Handling precautions

Oligonucleotides 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 the oligonucleotides.

Related products

It is recommended to include a positive and negative control, such as Dharmacon RNAi Controls Reagents, in every RNAi experiment. For more information, click [here](#) for RNAi controls.

Publication reference guide

When referencing the use of Dharmacon siRNA reagents, please include the following information: product name, catalog number, Dharmacon, Inc., Lafayette, CO.