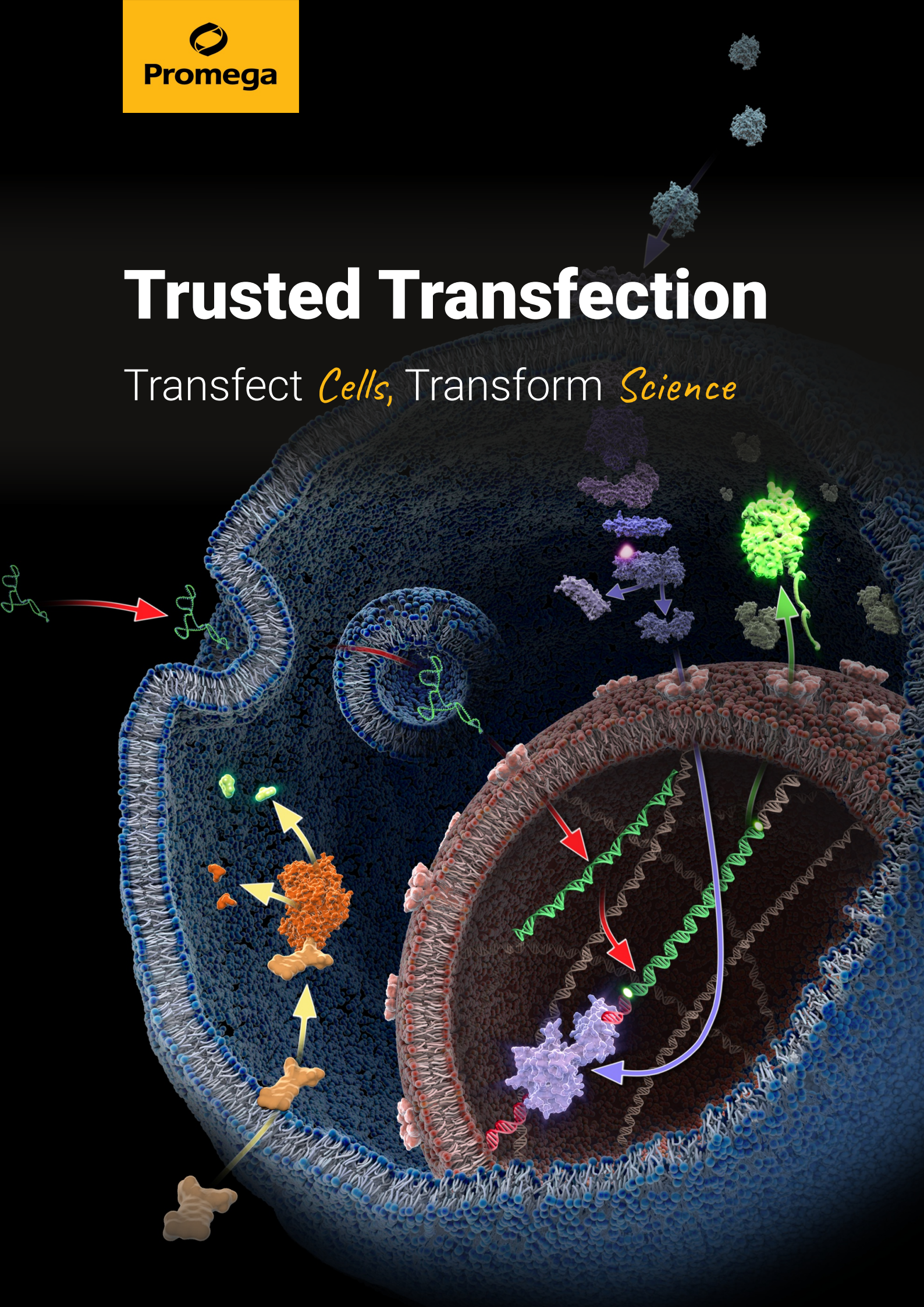
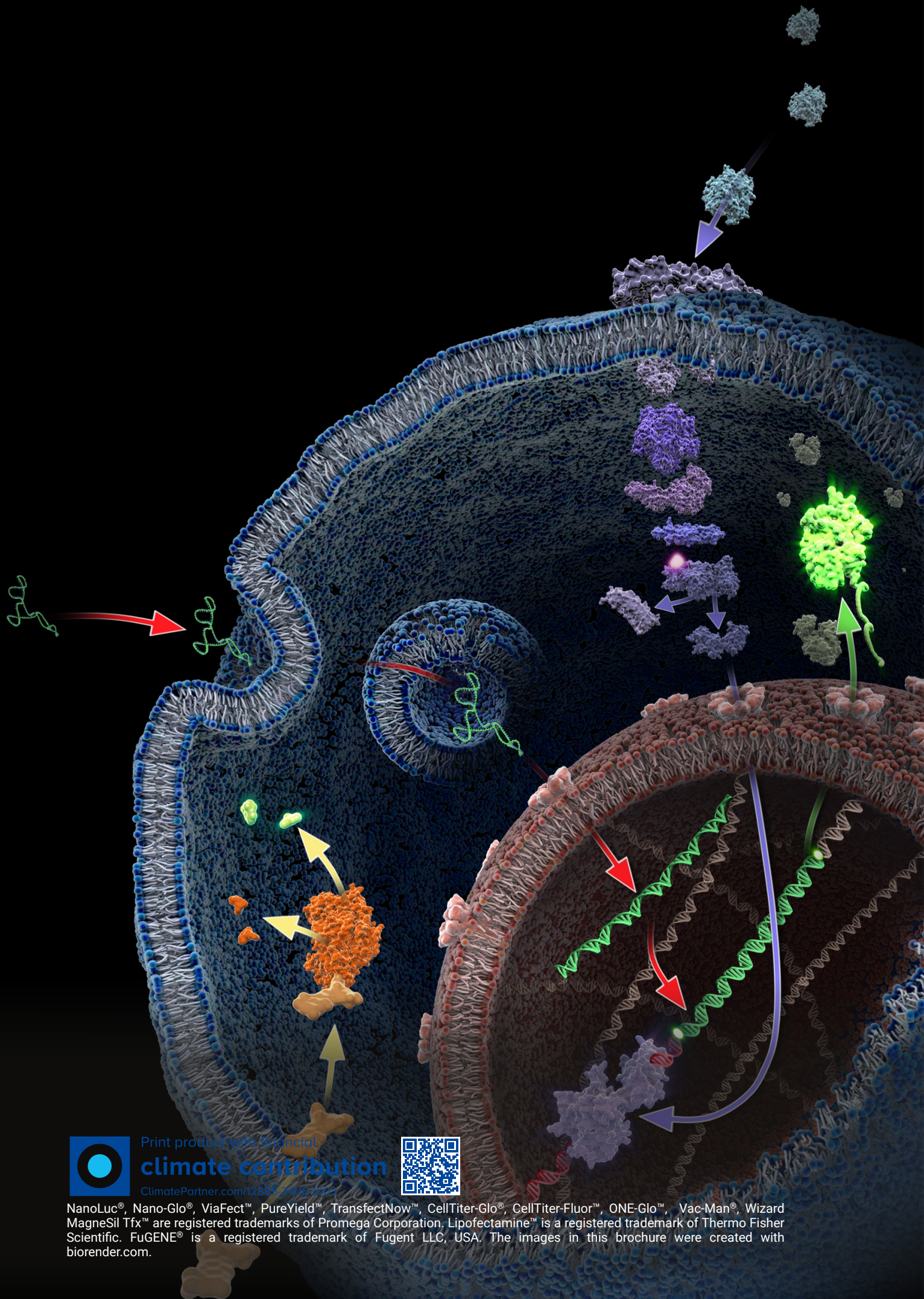


Trusted Transfection

Transfect *Cells*, Transform *Science*





Print product with financial
climate contribution

ClimatePartner.com/12885-2405-1023



NanoLuc®, Nano-Glo®, ViaFect™, PureYield™, TransfectNow™, CellTiter-Glo®, CellTiter-Fluor™, ONE-Glo™, Vac-Man®, Wizard MagneSil Tfx™ are registered trademarks of Promega Corporation. Lipofectamine™ is a registered trademark of Thermo Fisher Scientific. FuGENE® is a registered trademark of Fugent LLC, USA. The images in this brochure were created with biorender.com.

Table of Contents

Transfection 101	page 2
GloMax® Detection Systems	page 5
Best Choice Transfection Reagents	page 6
Best DNA Prep for your Transfection	page 8
Start Using HEK293 Cells as Reagents	page 9
Gentle and Efficient Transfection of dsRNA	page 10

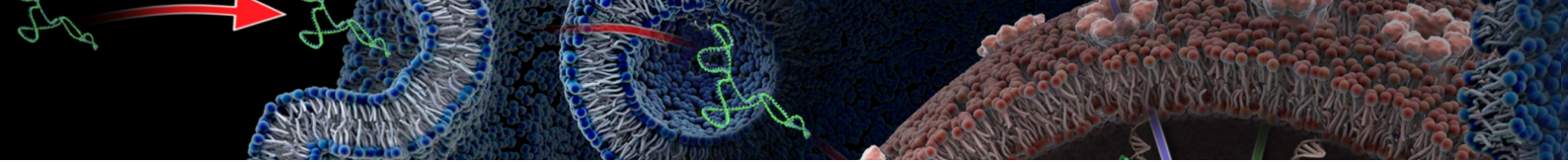


Are you interested in our latest research?

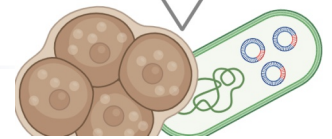
Check out our many on-demand webinars!

www.promega.com/webinars





Don't mix it up with transformation, that is transfer of DNA into bacteria or yeast.

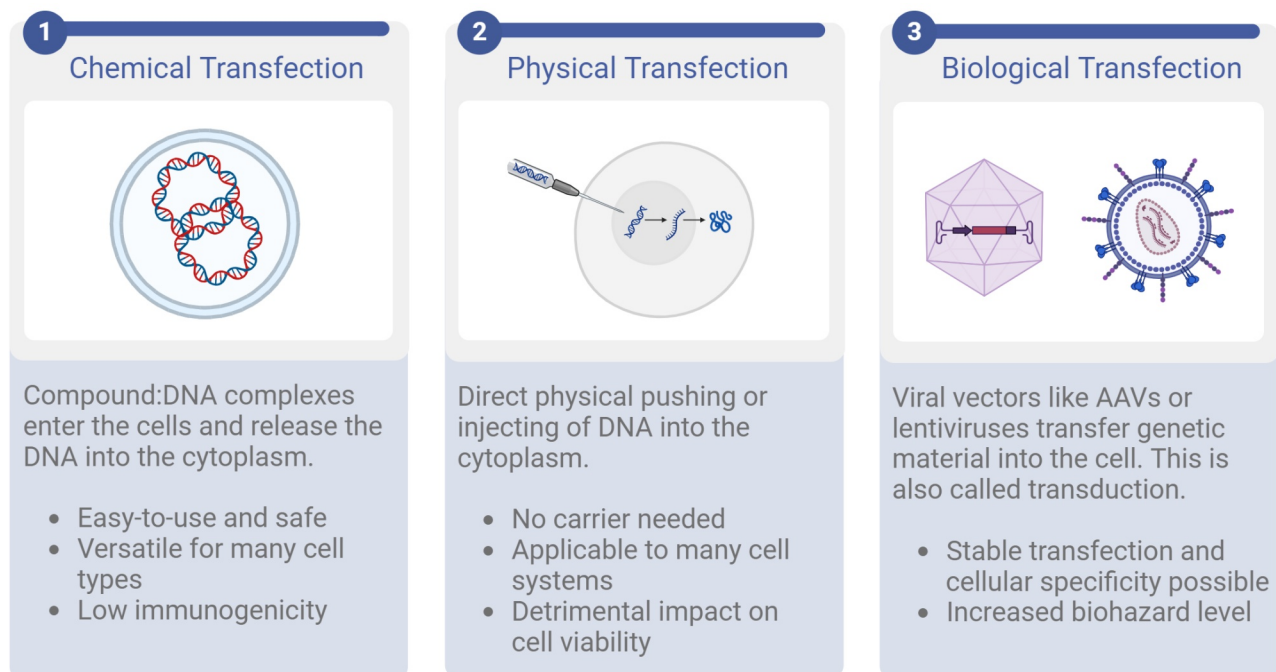


Transfection 101

Transfection Basics

Transfection allows for the direct manipulation of the genetic makeup of cells, providing a powerful means to explore genetic functions, develop new treatments, and advance scientific understanding in various fields of biology and medicine.

A large variety of transfection techniques for vertebrate and insect cells have been established.



Cell Type Specificity: Different reagents may perform better with specific cell types.

Cell Viability: Selecting a method that balances transfection efficiency with minimal cytotoxicity is key.

Stable vs Transient Transfection: Stable transfection is required for long-term studies. Transient transfection is suitable for short-term experiments and can be achieved with a broader range of reagents.

Scale of Transfection: Select a transfection technique that supports the scale of the experimental setup.

Safety and Regulatory Considerations: The safety profile of the transfection method and any regulatory compliance issues should be considered.

Key points to consider



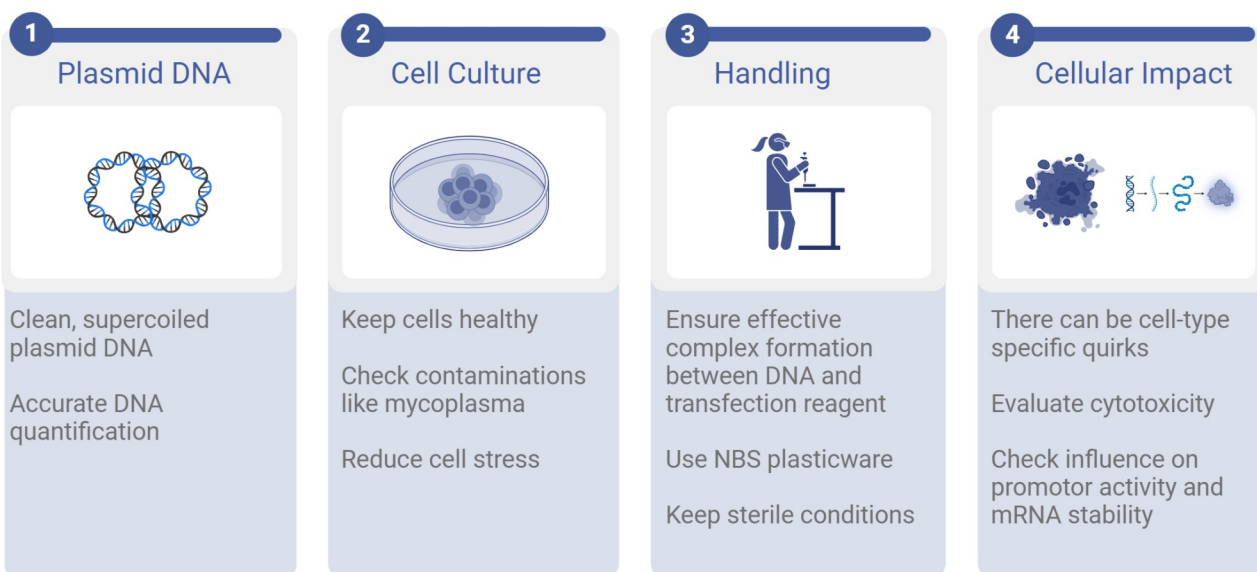
The right way to transfect ...

... can make or break your research

Transfection is crucial and significantly impacts experimental outcomes. Efficient transfection ensures effective DNA delivery, maintains cell health and supports gene expression.

Inconsistent or poor transfection can reduce assay sensitivity and may yield misleading data. Effective transfection reagents coupled with optimized protocols are essential for achieving reproducible and meaningful scientific results.

Mastering Transfection: Every Step Matters

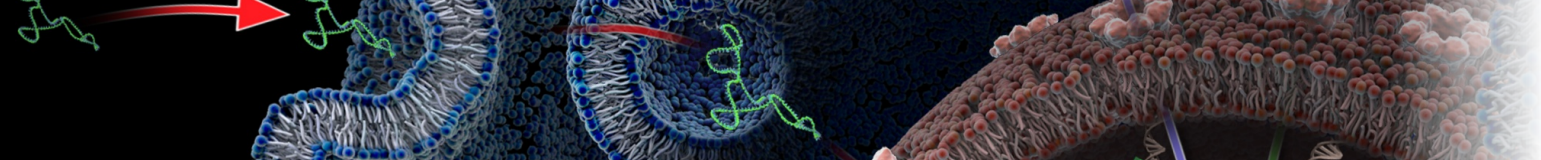


Do you want to learn more?

You'll find more insights in our transfection guide

www.promega.com/transfection-guide/





Trusted Transfection Reagents

Let biology, not transfectability dictate the cell line you work with

A diverse range of transfection reagents is necessary to meet the unique demands of each research project. Compatible with a wide array of downstream assays, these reagents are indispensable tools for cutting-edge research.

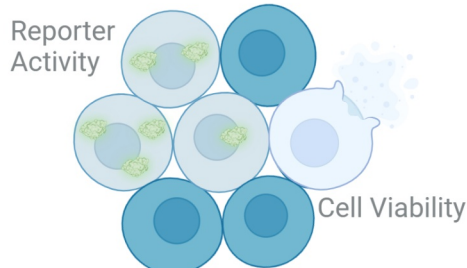
62% of test inquiries come from recommendations by fellow scientists, underscoring the trust placed in these reagents. Moreover, a significant proportion of those who test our reagents become regular buyers.

We're committed to ensuring that researchers find the perfect fit for their experiments. Thus, we provide both service support and sample requests to enable rigorous and informed testing.

How to Best Evaluate Transfection Reagents

1

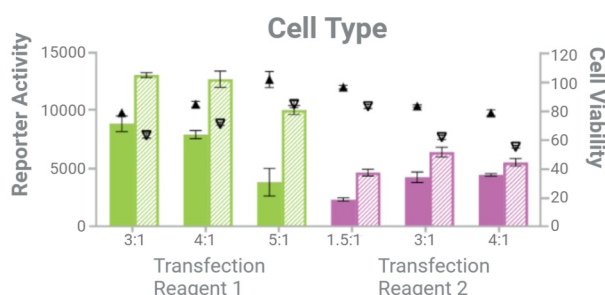
3 outcomes after transfection



1. Transfected Cells
2. Untransfected Cells
3. Dying/Damaged Cells

2

How to compare transfection reagents



1. Measure the **successful transfection rate** with a reporter
2. Measure **cell viability** with an accompanying assay

Get the highest degree of flexibility with the cell viability assays CellTiter-Glo® (when using a fluorescent reporter) and CellTiter-Fluor™ (when using a luminescent reporter) or profit from the ready-to-use solution for the firefly luciferase reporter:

ONE-Glo™ + Tox Luciferase Reporter and Cell Viability Assay



GloMax® Detection Systems

A Versatile, Reliable, and Intuitive Lab Companion

GloMax® Discover is an advanced multimode plate reader designed to provide optimal performance for Promega reagents with high-performance luminescence, fluorescence, UV-visible absorbance, BRET and FRET, two-color filtered luminescence, and kinetic measurement capabilities. GloMax® Discover can be used as a standalone plate reading instrument or integrated into high-throughput automated workflows. Results are easy to interpret using integrated data analysis software.

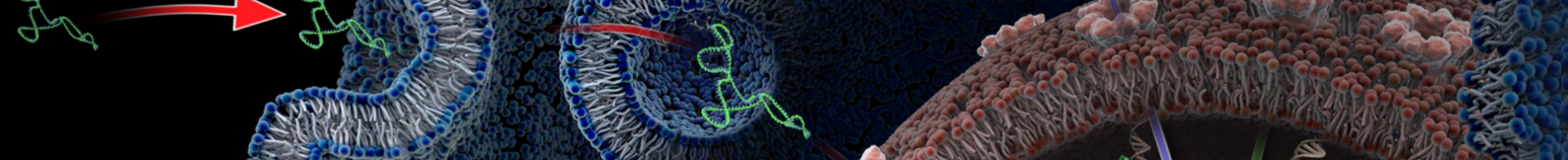
One Instrument. Numerous Applications.

- Reporter Gene Assays
- Cell Viability, Cytotoxicity & Apoptosis Assays
- Kinetic Measurements
- Multiplexing
- Oxidative Stress & Cell Metabolism
- ELISA
- BRET/FRET Analysis
- Lumit® Immunoassays

GloMax DISCOVER

A high-performance, easy-to-use multimode plate reader for luminescence, fluorescence, absorbance, BRET and FRET applications.





Best Choice Transfection Reagents

FuGENE® 6 – Reliable Routine Transfection

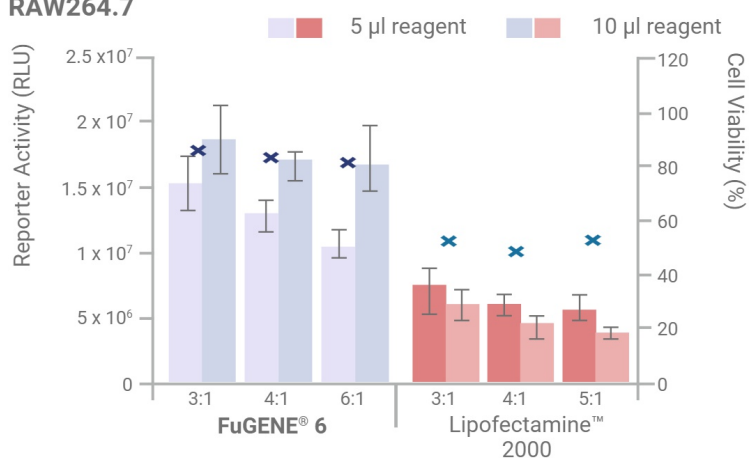
Simplify your daily transfections with FuGENE® 6, a versatile and gentle reagent designed to deliver consistent, reliable results across a broad range of routine laboratory cell lines.

FuGENE® 6

Best Choice For

- ✓ COS7
- ✓ U2OS
- ✓ Raw264.2
- ✓ K562

RAW264.7



FuGENE® HD – For Challenging Cell Lines and Advanced Applications

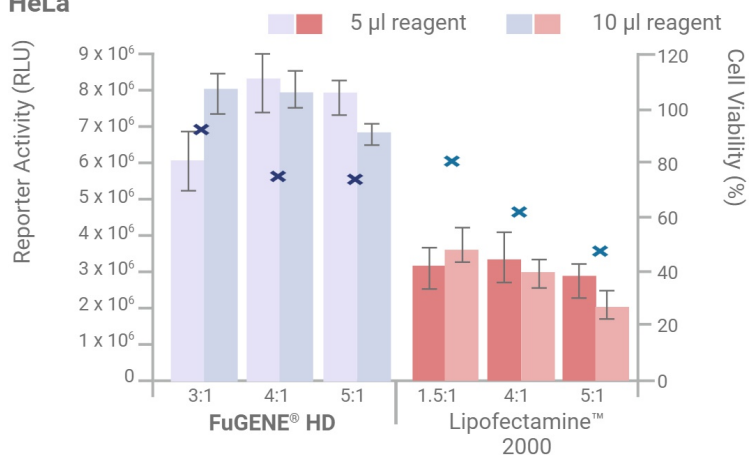
Elevate your research and tackle challenging cell lines and primary cells with FuGENE® HD, empowering you to explore new frontiers in stem cell studies as well as protein and virus production. FuGENE® HD's non-disruptive approach ensures high transfection efficiency without altering your cells' natural physiology.

FuGENE® HD

Best Choice For

- ✓ A549
- ✓ HeLa
- ✓ PC-3
- ✓ SH-SY5Y
- ✓ HepG2
- ✓ Sf9

HeLa





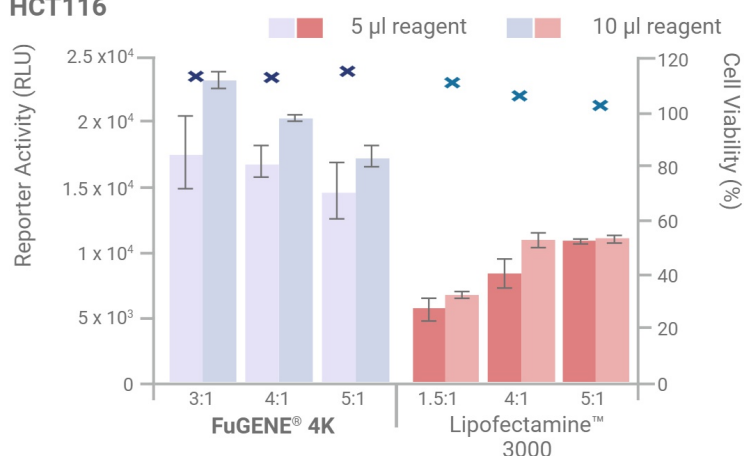
Are you working with other cell lines?
Contact Promega experts for a detailed recommendation
at techserv@promega.com



FuGENE® 4K – High-Efficiency Transfections Across Diverse Cells

Streamline your transfection processes with FuGENE® 4K's advanced, easy-to-use formula, perfect for achieving high-efficiency DNA delivery across a wide variety of cell lines and primary cells for advanced applications.

HCT116



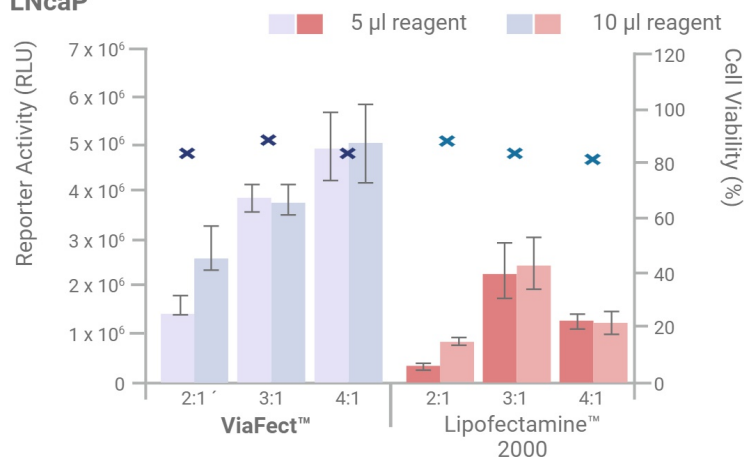
FuGENE® 4K Best Choice For

- ✓ HCT116
- ✓ H1299
- ✓ CHO-K1
- ✓ A375
- ✓ NIH3T3
- ✓ HEK293

ViaFect™ – Superior Performance in a Wide Range of Cells

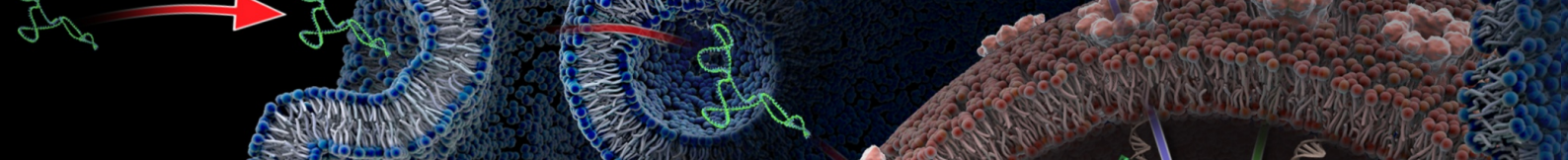
Enhance the efficiency and scope of your transfections with ViaFect™. It's engineered to deliver superior performance across a diverse array of cell lines, including those traditionally challenging to transfect. Optimize your experiments with the versatility and low toxicity of ViaFect™, ensuring high-level gene expression while maintaining the health and viability of your cells.

LNcaP



Viafect™ Best Choice For

- ✓ C2C12
- ✓ LNCaP
- ✓ Jurkat
- ✓ HT-29
- ✓ MCF7
- ✓ PC-12
- ✓ H9C2
- ✓ C2C12



Best DNA Prep For Your Transfection

PureYield™ Plasmid Systems

Get clean plasmid DNA for every day applications using the PureYield™ Plasmid System, available as Mini-, Midi-, and Maxipreps.

Benefit from the fast and streamlined workflow using the Vac-Man® Laboratory Vacuum Manifold together with a vacuum pump.



Wizard MagneSil Tfx™ System

Do you need higher throughput but still want clean, transfection-ready plasmid DNA?

Prepare your plasmids in 96-well plates using the Wizard MagneSil Tfx™ System. DNA capture on paramagnetic particles enables complete automation of the workflow.

1

Compatibility



The Wizard MagneSil Tfx™ System is compatible with automation on all major platforms, including Hamilton Star/Nimbus, Tecan Freedom Evo/Fluent, Biomek i5/i7, Eppendorf epMotion, and particle movers such as KingFisher.

2

Field Support Access



When you purchase a Promega extraction solution, you gain free access to our elite field support scientists, method development services, and technical support. Contact us if you need help setting up the Wizard MagneSil Tfx™ System on your robot.

3

Customization



With decades of experience in building personalized automation solutions, we can help you customize the scale, application, and chemistry, regardless of platform, exactly to your needs—the only limit is your imagination.

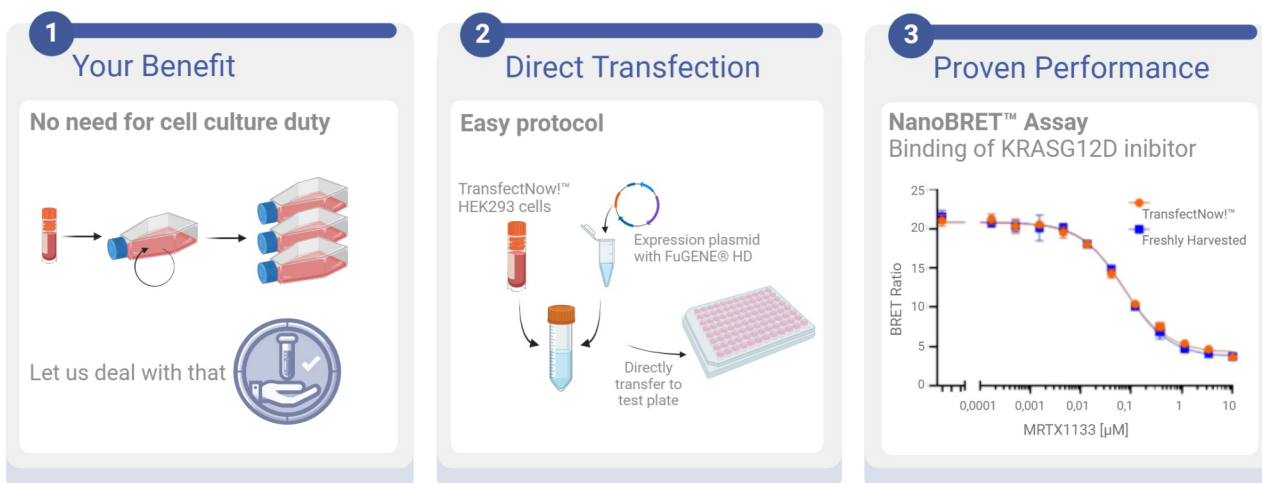
Start using HEK293 cells as reagents

TransfectNow™ HEK293 Cells

Did you know that the quality of HEK293 cells and their culture conditions have profound effects on transfection. Not only is prior cell culture work and time intensive, but it is also a major source of variability between experiments.

Transfection almost as easy as with *E. coli*. Obtain your results faster with TransfectNow™ HEK293 Cells. Transfect the cells right after thawing and without cell culture duty. Robust quality controls ensure that you can trust your transfection and focus on more important elements in your research. TransfectNow™ cells are best used with FuGENE® HD transfection reagent.

Inside View - How It Works



The TransfectNow™ HEK293 cells enable direct transfection and implementation of the downstream assays. Compare the exemplary results obtained using either TransfectNow™ cells or freshly harvested HEK293 cells. The binding of MRTX1133 to activated KRAS is tested using our NanoBRET™ Target Engagement Assay and virtually no difference is observed in the results between the two options (IC50 values: 0.071 with TransfectNow™ cells and 0.076 with freshly harvested cells)

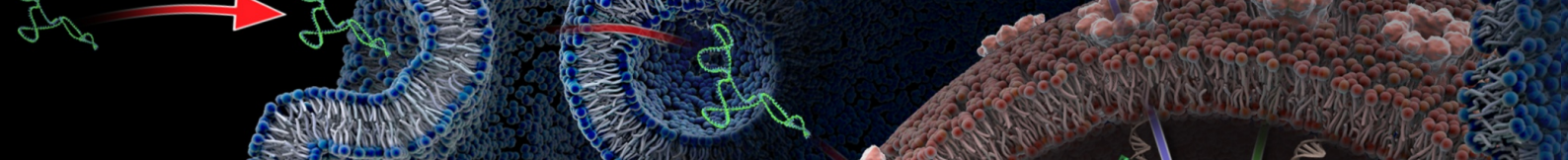


Are you interested in advancing drug discovery?

Then read our eBook **Assays for Drug Discovery & Development**

www.promega.com/drug-discovery-brochure



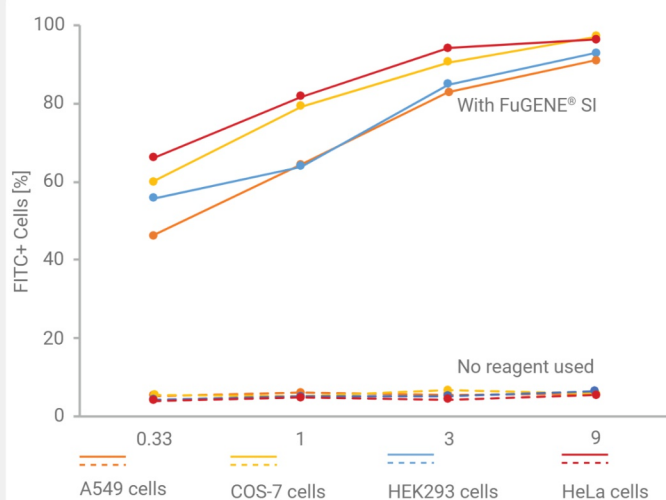


Gentle and Efficient Transfection of dsRNA

FuGENE® SI Transfection Reagent

The 100% synthetic FuGENE® SI Transfection Reagent is a high-efficiency, low-toxicity solution for delivering siRNA, miRNA, and other small RNA molecules into a wide range of eukaryotic cell lines.

Particularly suitable for researchers seeking maximum knockdown efficiency with minimal cell disruption.

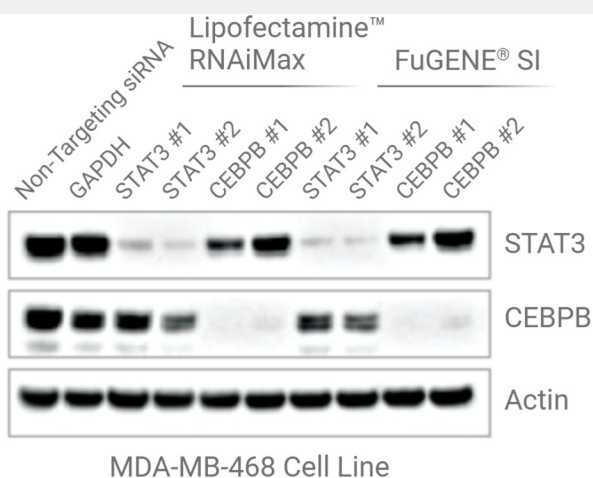


FuGENE® SI provides efficient transfection of labeled small RNAs

A549, COS-7, HEK293 and HeLa cells plated in 96-well microplates were transfected with 0.4 μ l of FuGENE® SI or a no-reagent negative control and various amounts of FITC-labeled control siRNA.

24 h after transfection, the cells were analyzed via flow cytometry for FITC-positive signal.

As expected, the FITC signal is near zero when no reagent is used and is high when using FuGENE® SI in diverse cell lines.



Effective knockdown of endogenous target genes with FuGENE® SI

Human MDA-MB-468 cells were reverse-transfected in 6-well plates with 40 pmol of the indicated targeting siRNAs or negative/positive controls with 7.5 μ l of either FuGENE® SI Transfection Reagent or Lipofectamine™ RNAiMAX per well. The cells were analyzed 72 h after transfection via Western blot for expression of target genes STAT3 and CEBPB.

Pay attention to the similar or better knockdown performance by FuGENE® SI.



Products in Each Section at a Glance

1 Transfection 101

Product	Quantity	Cat. #
Firefly Luciferase Control Vectors	20 µg	E5011
ONE-Glo™ EX Luciferase Assay System	10 ml	E8110
ONE-Glo™ + Tox Luciferase Reporter and Cell Viability Assay	1 plate	E7110
NanoLuc® Luciferase Control Vectors	20 µg	N1441
Nano-Glo® Luciferase Assay System	10 ml	N1110
CellTiter-Glo® 2.0 Cell Viability Assay	10 ml	G9241
CellTiter-Fluor™ Cell Viability Assay	10 ml	G6080
Transfection Carrier DNA	5 x 20 µg	E4881

2 Best Choice Transfection Reagents

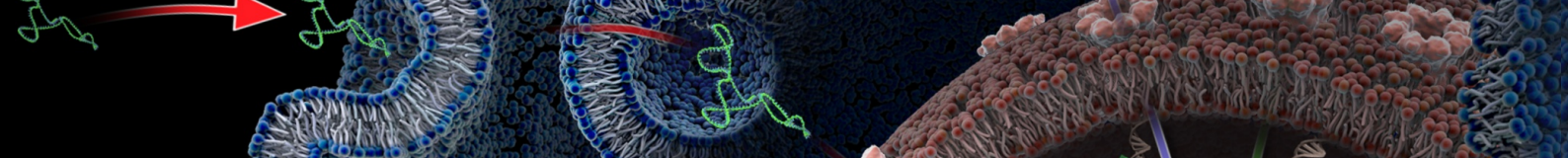
Product	Quantity	Cat. #
FuGENE® 6 Transfection Reagent	1 ml	E2691
	5 x 1 ml	E2692
FuGENE® HD Transfection Reagent	1 ml	E2311
	5 x 1 ml	E2312
FuGENE® 4K Transfection Reagent	1 ml	E5911
	5 x 1 ml	E5912
ViaFect™ Transfection Reagent	0.75 ml	E4981
	2 x 0.75 ml	E4982
FuGENE® SI Transfection Reagent	1 ml	E9311
	5 x 1 ml	E9312

3 Best DNA Prep For Your Transfection





Product	Quantity	Cat. #
PureYield™ Plasmid Miniprep System	100 / 250 preps	A1223 / A1222
PureYield™ Plasmid Midiprep System	25 / 100 preps	A2492 / A2495
PureYield™ Plasmid Maxiprep System	10 / 25 preps	A2392 / A2393
Vac-Man® Laboratory Vacuum Manifold	1 each	A7231
Wizard MagneSil Tfx™ System	4 x 96 preps	A2380

4 TransfectNow™ HEK293 Cells

Product	Quantity	Cat. #
TransfectNow™ HEK293 Cells	1 x 0.5 ml	NC1001
	2 x 1 ml	NC1002



Worldwide Promega Contact Information

-  Headquarters
-  R&D and Production
-  Branches
-  Distributors



Europe

Germany/Austria

Tel: +49 6227 6906 0
Fax: +49 6227 6906 222
de_custserv@promega.com

Switzerland

Tel: +44 878 90 00
Fax: +44 878 90 10
ch_custserv@promega.com

France

Tel: +33 (0)437 2250 00
Fax: +33 (0)437 2250 10
contactfr@promega.com

Italy

Tel: +39 0254 0501 94
Fax: +39 0254 0501 45
customer.service.italia@promega.com

Poland

Tel: +48 22 531 0667
Fax: +48 22 531 0669
pl_custserv@promega.com

Central/Eastern Europe

Tel: +49 6227 6906 0
Fax: +49 6227 6906 222
de_ceesales@promega.com

Spain

Tel: +34 902 538 200
Fax: +34 902 538 300
esp_custserv@promega.com

Belgium/Luxembourg/Netherlands

Tel: +31 71 532 42 44
Fax: +31 71 532 49 07
benelux@promega.com

Denmark, Estonia, Finland, Iceland, Norway, Sweden

Tel: +46 8 452 2450
Fax: +46 8 452 2455
sweorder@promega.com

United Kingdom

Tel: +44 23 8076 0225
Fax: +44 20 8076 7014
ukcustserve@promega.com

North America

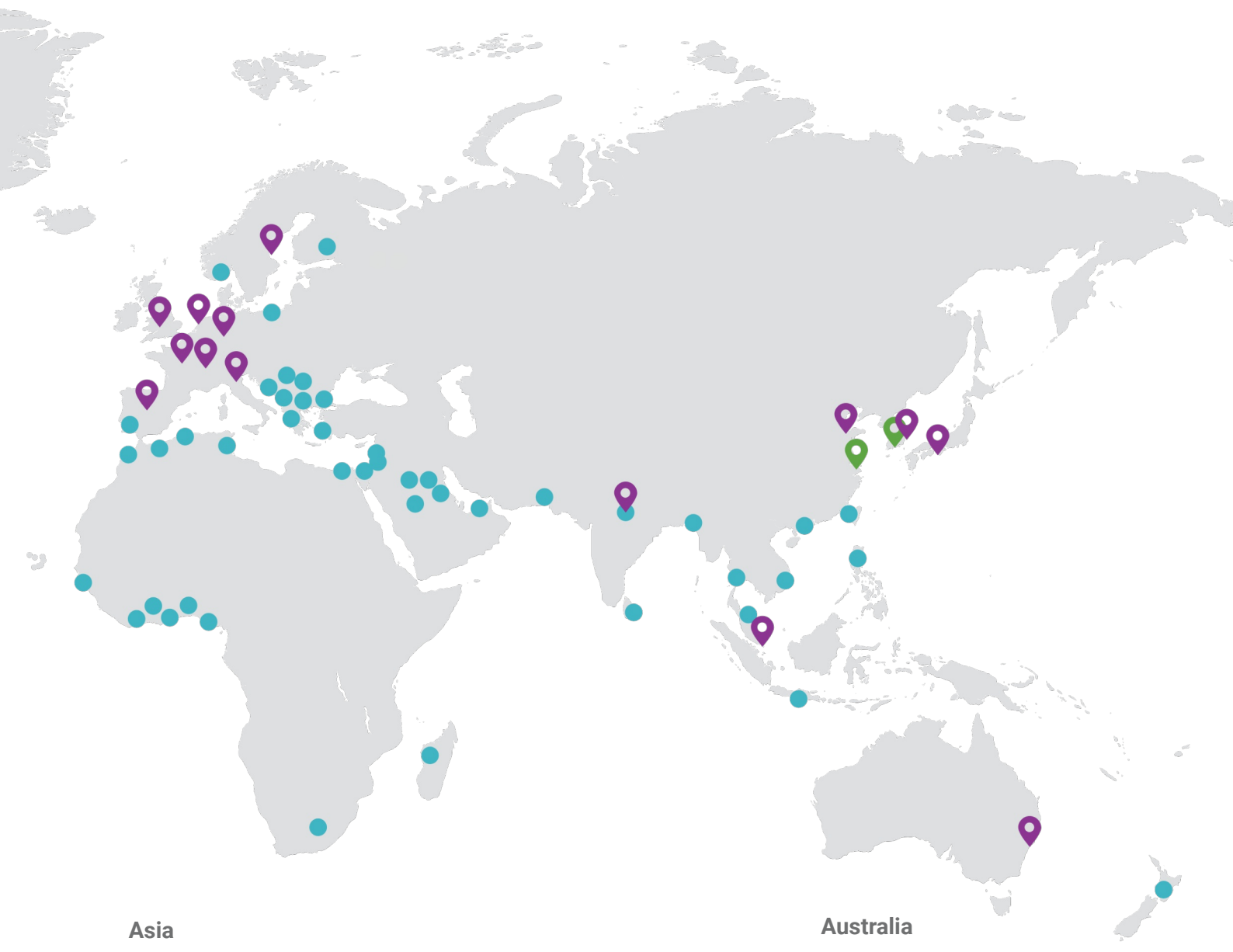
USA/Canada

Tel: +1 608 274 4330
Fax: +1 608 277 2516
custserv@promega.com

South America

Brazil

Tel: +55 11 5090 3780
Fax: +55 11 5096 3780
promega.brasil@promega.com



Asia

China

Tel: +86 10 5825 6268
Fax: +86 10 5825 6160
info@promega.com.cn

Japan

Tel: +81 3 3669 7981
Fax: +81 3 3669 7982
jpmarketing@jp.promega.com

Korea

Tel: +82 2158 83718
Fax: +82 2262 85418
CustServiceKR@promega.com

India

Tel: +91 11 43005814/15/16/17
Fax: +91 11 41035028
ind_custserv@promega.com

Asia-Pacific region, Singapore

Tel: +65 6513 3450
Fax: +65 6773 5210
sg_custserv@promega.com

Australia

Australia

Tel: +61 2 8838 3800
Fax: +61 2 8838 3855
austcustserv@promega.com

For more information about Promega branches
and distributors:
www.promega.com/support/contact-promega

