



[Return to Helix](#)



[Check Out](#)

[Quick Order](#)

[My Orders](#)

[Find My Gene™](#)

[My Helix](#)

[Register](#)

[My Account](#)



[Log In](#)

[Check Out](#)

[Quick Order](#)

Products + Services

My Account

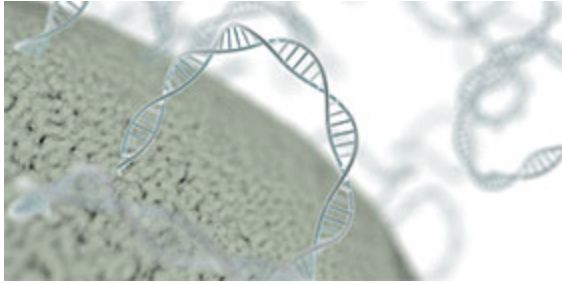
Support

Search



[Home »](#)

Products + Services



ViaFect™ Transfection Reagent

Generate more relevant data with efficient transfection of your optimal cell line without compromising cell viability.

[Product Details and Ordering](#)

[Learn More + Request a Sample](#)

ALL CATEGORIES

[Biobanking](#) >

[Biochemicals and Labware](#) >

[Biologics](#) >

[Cell Health and Metabolism](#) >

[Cell Signaling](#) >

[Cloning and DNA Markers](#) >

[DNA and RNA Purification](#) >

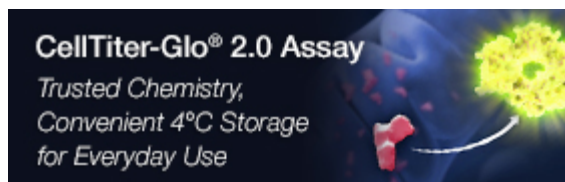
[Drug Discovery](#) >

[Epigenetics](#) >

[Genetic Identity](#) >

[Imaging and Immunological Detection](#) >

[Industrial and Environmental Monitoring](#) >



[View all Products »](#)

[What's New »](#)

FEATURED PRODUCTS + SERVICES

CellTiter-Glo® 2.0: A Luminescent Cell Viability Assay for Fast, Easy, Everyday Use

Promega GoTaq® G2 Polymerase at a Glance

Metabolism Assays

\$name

Quantus™ Fluorometer: Simple, Sensitive DNA and RNA Quantitation

NEED ASSISTANCE?



[Chat with Customer Service](#)

[Chat with a Scientist](#)

[Contact us](#)

About Us

[Careers](#) [Company Info](#) [Contact Us](#) [ISO Certification](#) [Press Releases](#) [Corporate Responsibility](#)

Connect With Us



Get Connected with Us

[Not your country?](#)

© 2015 Promega Corporation. All rights reserved.

[Legal and Trademarks](#) [Privacy Policy](#) [Terms and Conditions](#)