



Custom Assay Services from Promega

Outline



Promega & Custom Assay
Services Team

What We Do

How We Do It

Case Studies

Support

Future Directions



World-Wide Connection



- Founded 1978 • Revenues ~\$250M USD • Over 50 distributors in Europe, Asia, Africa and the Americas
- ~1,000 employees worldwide • ISO 13485 certified • Manufacturing in US, Korea and Shanghai
- Over 2,000 products for life science research and applied science including drug discovery

40891A

Solid Infrastructure and Intellectual Foundation



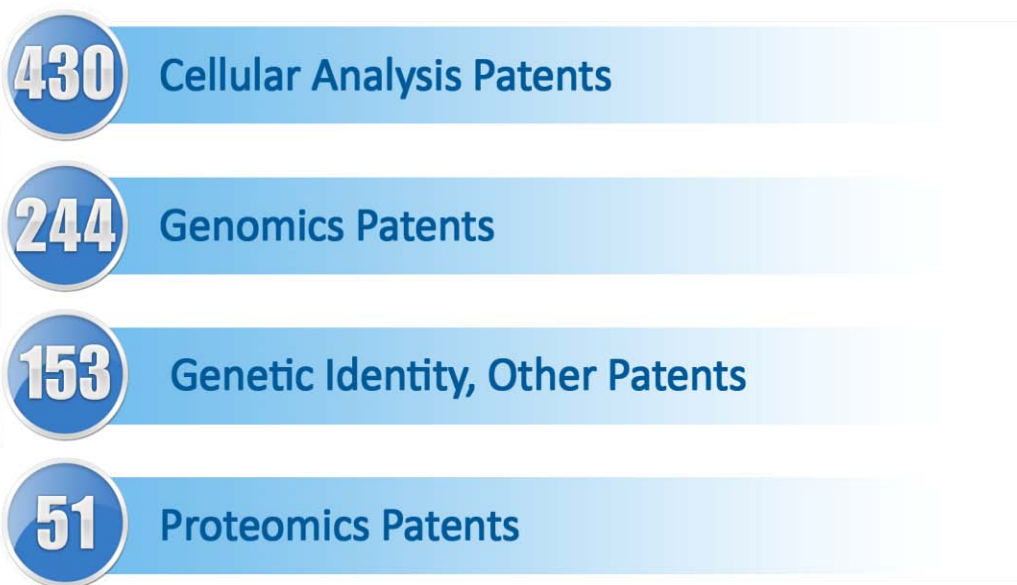
Quality System Overview

Promega Madison achieved ISO 9001 Certification in 1998, ISO 13485 in 2006

Quality system approach – complements and supports business objectives

Process driven with heavy emphasis on measures to drive performance

Intellectual Property



World-Class Capabilities in Custom Assay Development



Promega has the tools, expert staff and state-of-the-art facilities to support complete custom solutions for biochemical & cell-based assays

CELL ENGINEERING

Target expression in both normal & disease states

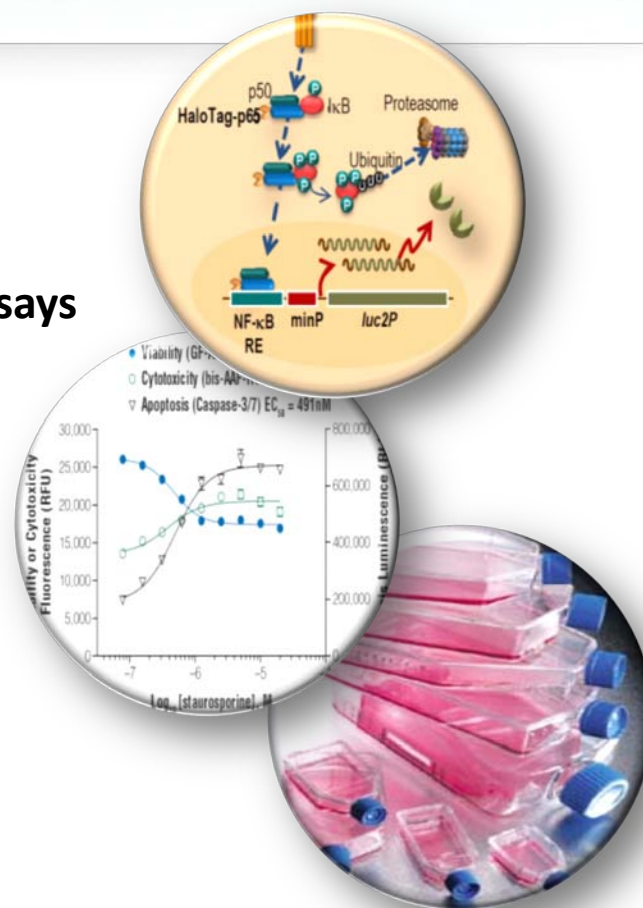
ASSAY DEVELOPMENT & QUALIFICATION

Target class expertise & multiple assay formats

ASSAY READY CELLS IN-SCALE

Pharmacological validation & unparalleled client support

And we provide all the post-delivery support to ensure your assay works in your hands



Value Proposition to Drug Developers



Let us partner with you to develop a complete biology-driven, Promega technology-enabled custom solution.

Faster time to data

Fewer in-house resources consumed & increased productivity

Flexible approach & offerings –
we can adapt our process to meet your needs

Flexible by Design



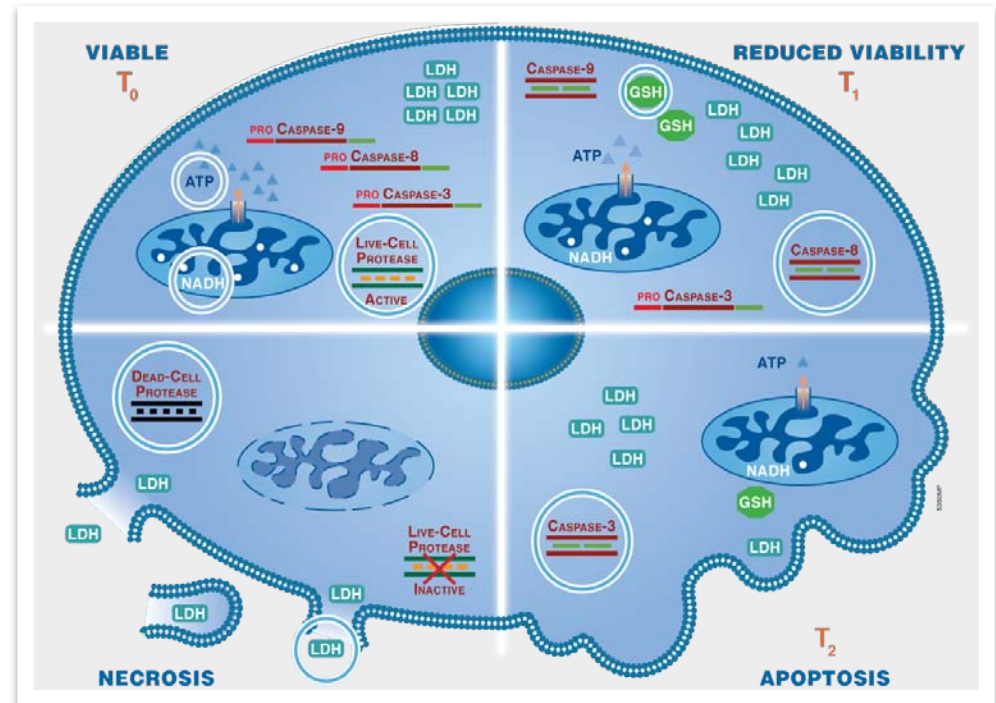
Whether your need is in small molecule drug discovery or developing biologics, we can create an assay that suits your requirements:

Cell-based assays

Biochemical assays

Multiple detection technologies

Multiplexed readouts for increased biological understanding



Forward Thinking...



We have solutions for your needs today and innovative thinking for your needs tomorrow.

Our scientists and business teams continue to enter into research collaborations with leading laboratories world-wide to investigate novel solutions, such as stem cells & whole animal imaging

Scientific & Assay Development Expertise



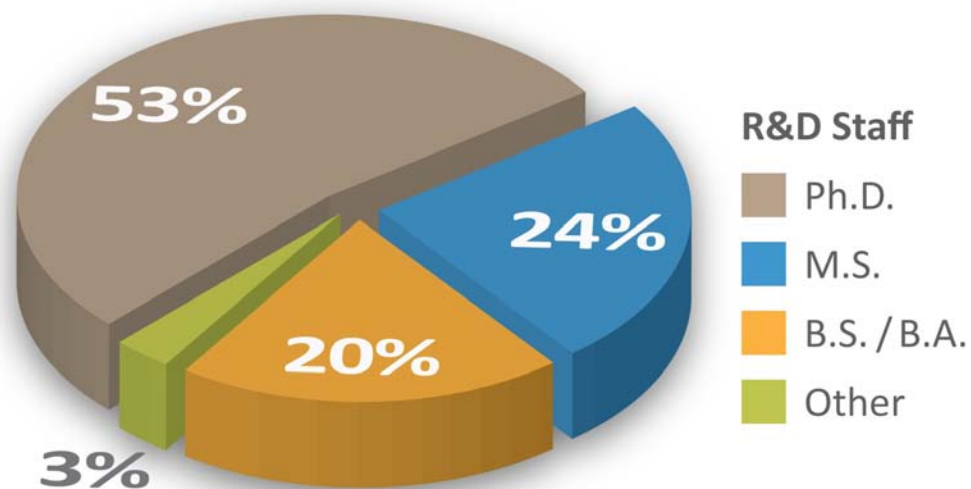
Global Leader in Bioluminescence Chemistry

R&D facilities located in:

Madison, WI
San Luis Obispo, CA
Seoul, Korea
Sunnyvale, CA

Extensive experience in developing assays for:

Genetic reporters
Cell viability & apoptosis
ADME/Tox
GPCRs
Kinases
In vivo imaging
Detection instrumentation



Transparent Collaboration Process



The Client is the Partner throughout all steps



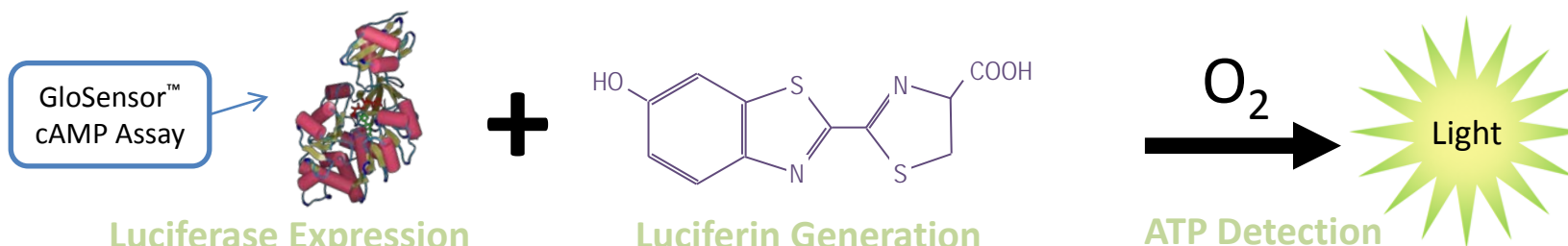
Comprehensive Portfolio of Assays for Cell Analysis

- CellTiter-Glo® Assay for monitoring cell viability
- Caspase-Glo® Assay for monitoring apoptosis
- Multiplexed viability, toxicity and apoptosis assays for more comprehensive monitoring of cellular health
- Luciferase reporter assays for cellular pathway analysis
- HaloTag® portfolio of live-cell protein trafficking, protein labeling, and purification tools
- GloSensor™ technology for the live-cell monitoring of cAMP, an important second messenger in GPCR signaling



And a comprehensive portfolio of reagents for genomics, proteomics and applied markets

Applying the Power of Bioluminescence to Biology



Cell-Based Assays

ONE-Glo™ Luciferase Assay System
Bright-Glo™ Luciferase Assay System
Steady-Glo® Luciferase Assay System
Dual-Glo™ Luciferase Assay System

Fluorescent Multiplex-Friendly Cell Health Assays

CellTiter-Fluor™ Cell Viability Assay
CytoTox-Fluor™ Cytotoxicity Assay
MultiTox-Fluor™ Multiplex Cytotoxicity Assay

Cell-Based Assays

ApoTox-Glo™ Triplex Assay
ApoLive-Glo™ Multiplex Assay
CytoTox-Glo™ Cytotoxicity Assay
Caspase-Glo® 3/7 Assay System
Caspase-Glo 8 & 9 Assay System
GSH-Glo™ Glutathione Assay
MultiTox-Glo Multiple Cytotoxicity Assay
P450-Glo™ Assay Screening System
Proteasome-Glo™ Cell Based Assays

Biochemical Assays

Calpain-Glo™ Assay
Caspase-Glo 2 & 6 Assay
DPPIV-Glo™ Assay
DUB-Glo™
MAO-Glo™ Assay
P450-Glo™ Assay Screening Systems
Proteasome-Glo™ Assays
UGT-Glo™ Assay

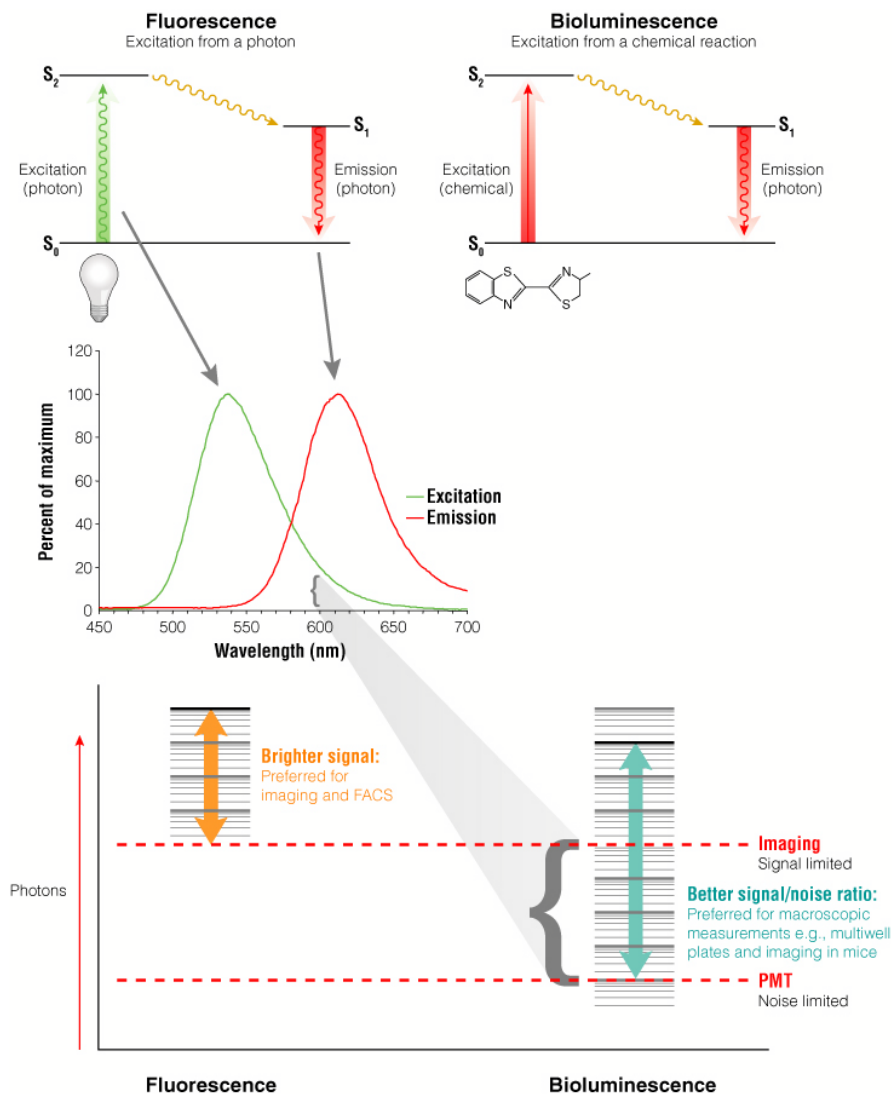
Cell-Based Assays

CellTiter-Glo® Luminescent Cell Viability
BacTiter-Glo® Microbial Cell Viability Assay
cAMP-Glo™ Assay

Biochemical Assays

ADP-Glo™ Kinase Assay
Kinase-Glo® Assay
Kinase-Glo® Plus Assay
Kinase-Glo® Max Kinase Assay
PDE-Glo™ Assay
Pgp-Glo™ Assay System

The Physics of Bioluminescence is Key to Better Biology



Luminescent assays avoid issues of fluorescence interference and offer a greater dynamic range when needed

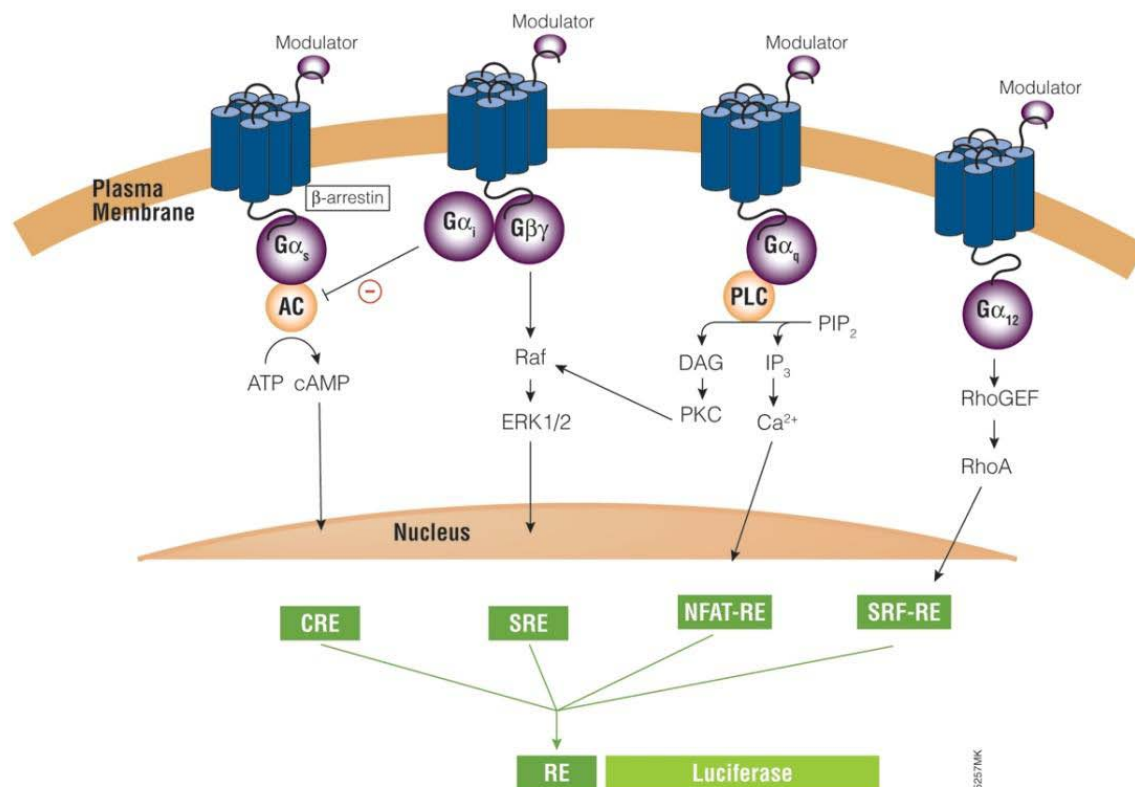
Better signal-to-noise ratio for multi-well plate assays & in vivo imaging in mice

Ultra-Glo is the best luciferase for cell-based luminescence protocols

Easier multiplexing allowed when luminescence & fluorescence combined

Luminescence detection transfers from biochemical to cell-based to animal studies seamlessly

Solutions for GPCR Biology



GloSensor™ cAMP Assay:
Live-cell HTS assay

cAMP-Glo™ Assay:
Lytic HTS assay

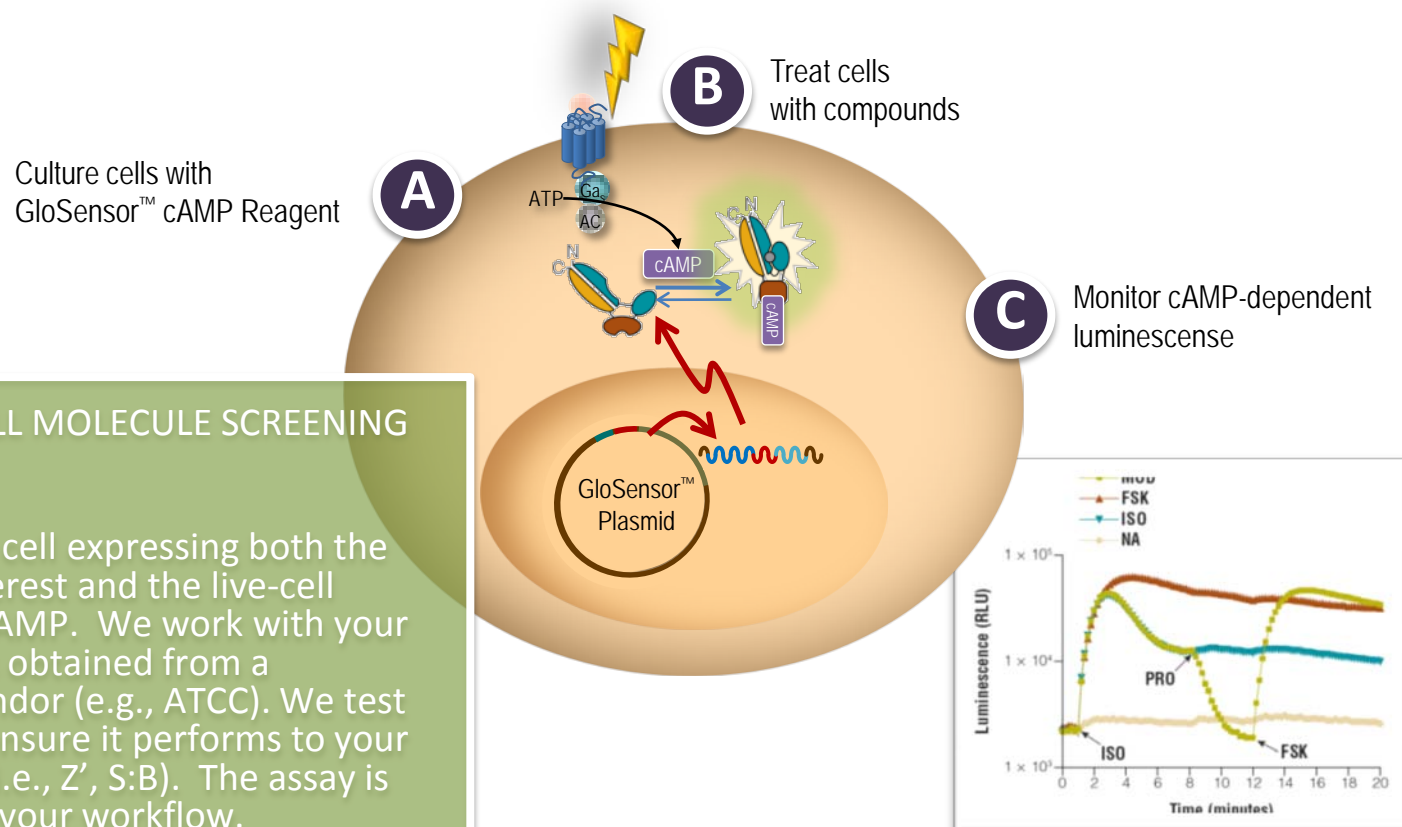
Luciferase reporter assays for:
G12/13
cAMP-RE
Ca $^{2+}$ -RE
ERK: SRE-RE

HaloTag® Assay:
Receptor internalization via protein trafficking

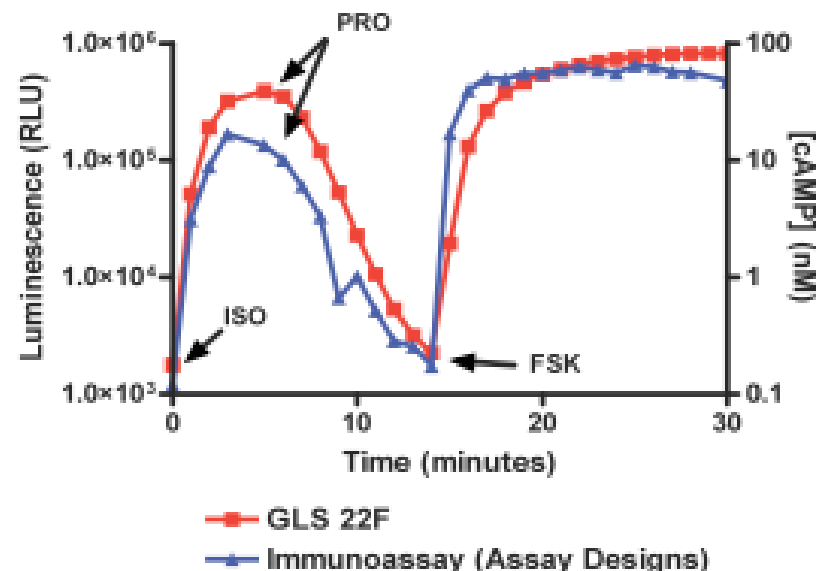
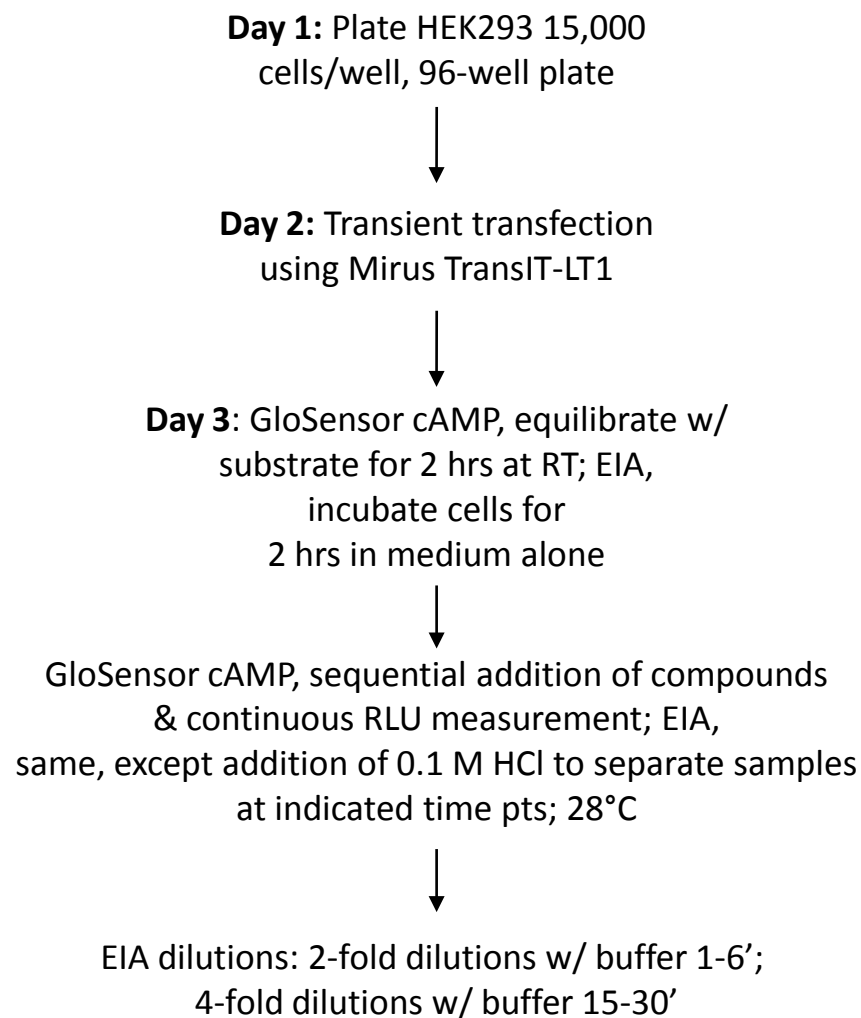
Monitor cAMP Levels in Real-time with a Live-cell, Non-lytic Bioluminescent Assay



Monitor Gs- and Gi-coupled GPCRs Using GloSensor™ cAMP Assay



GloSensor™ cAMP Assay is as Sensitive as an Immunoassay

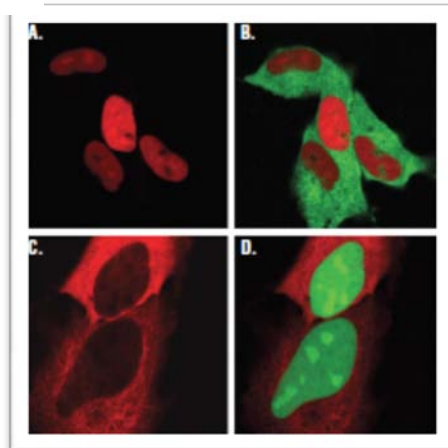
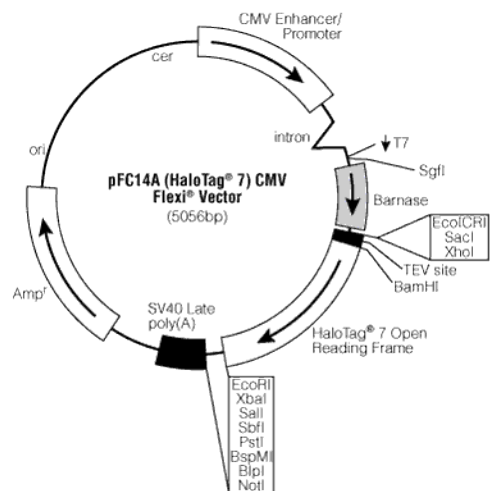
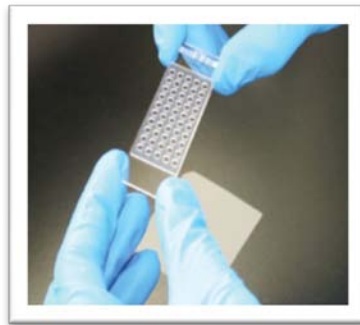
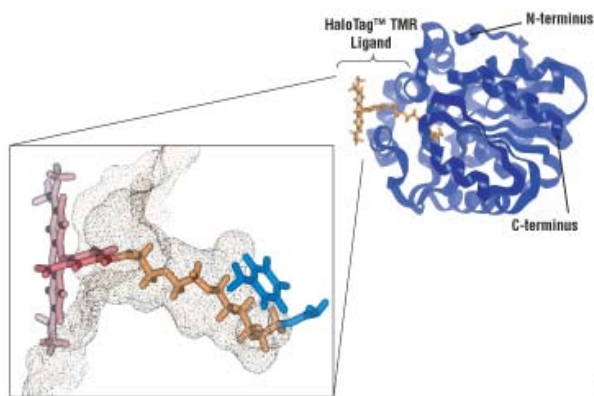


Assay Designs, Cat.# 901-066, acetylated

Monitor Protein Trafficking, Label Live Cells and So Much More



Using HaloTag® Protein Labeling Technology



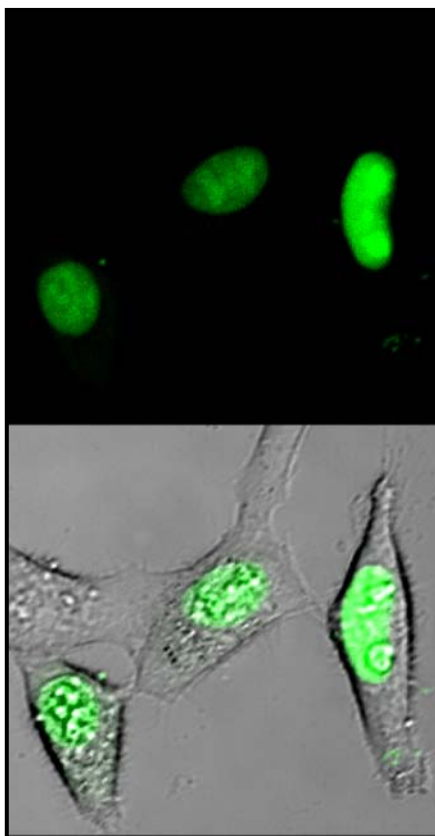
**A wide variety of Flexi Vectors,
Ligands and other label tools
allow:**

- Live cell imaging
- Protein localization
- Protein trafficking
- Protein capture
- Cell-to-gel analysis
- Protein:protein interaction analysis
- Protein:DNA interaction analysis
- Protein purification

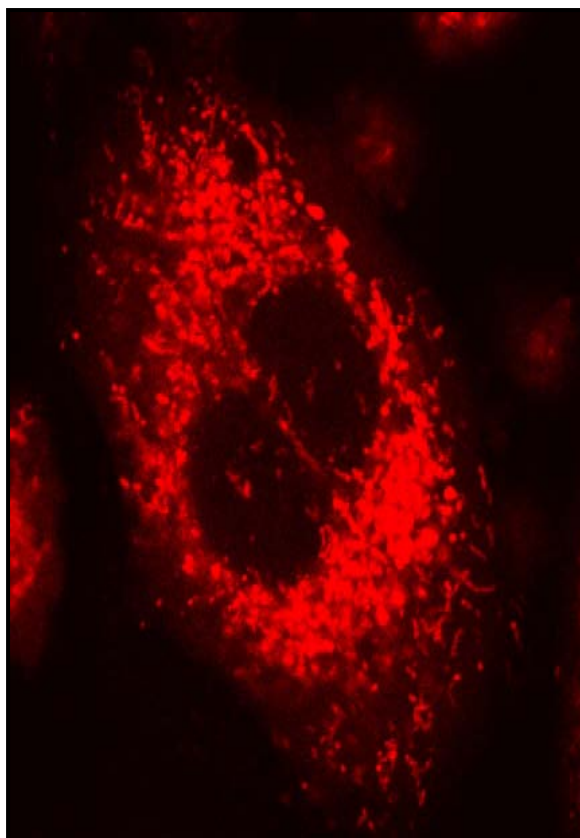
Sub-cellular targeting of HaloTag fusions



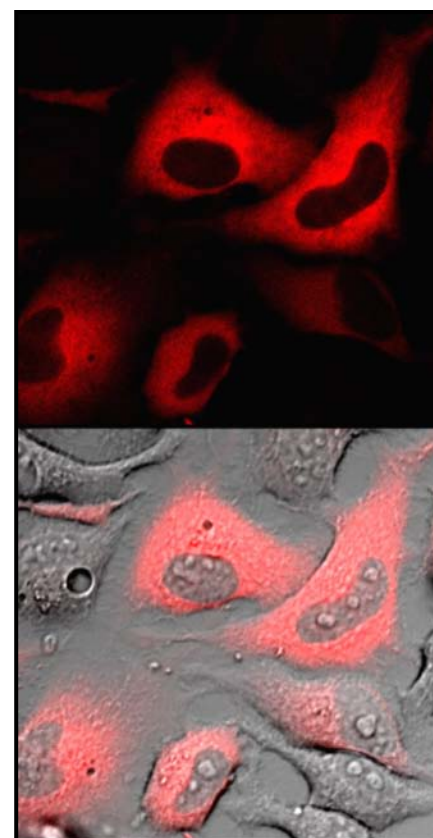
Nucleus
HaloTag - NLS₃



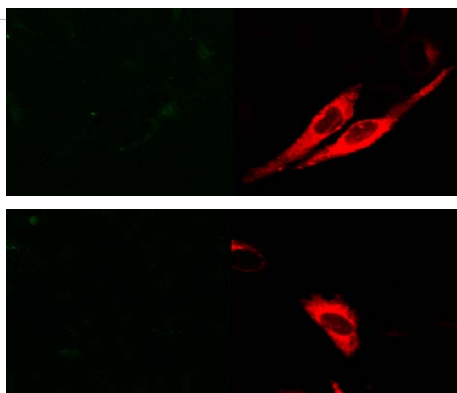
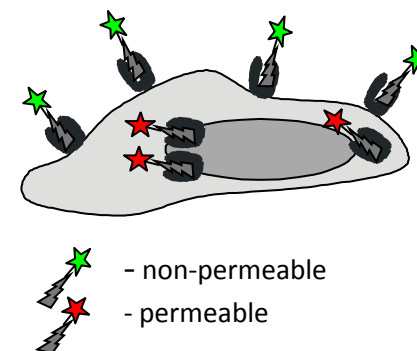
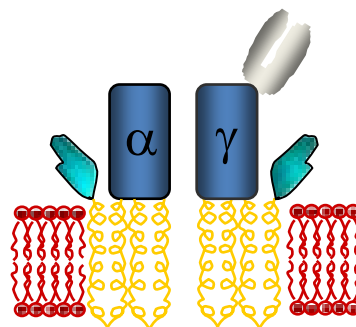
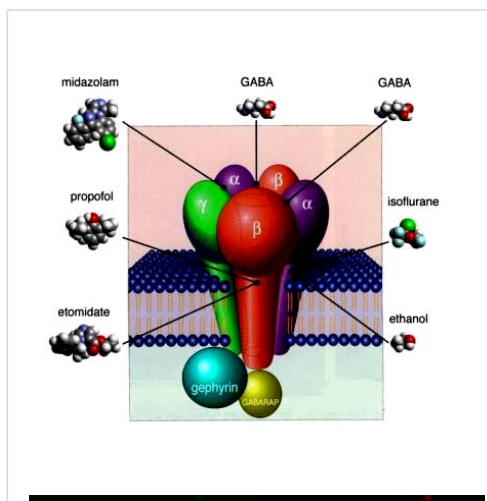
Mitochondria
Mito - HaloTag



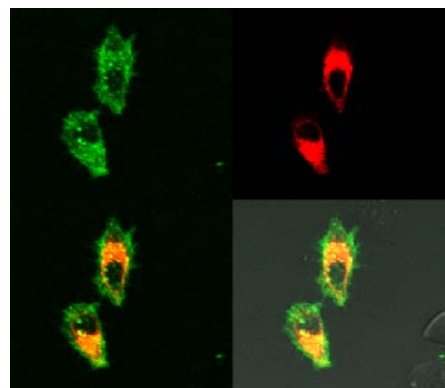
Cytosol
p65 - HaloTag



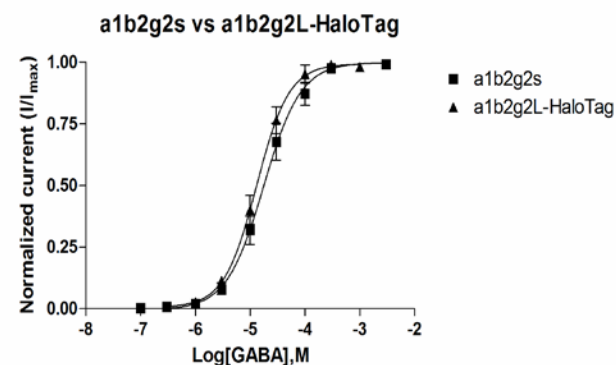
GABA_A receptor, a ligand-gated ion channel super family



γ2L-HaloTag co-expressed with α1 or β2 does not traffic to cell surface

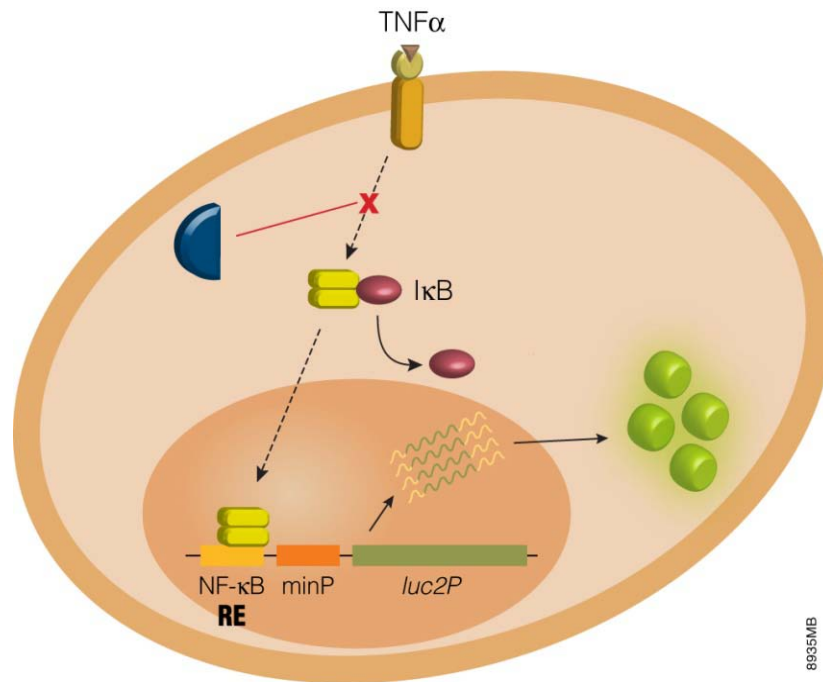


γ2L-HaloTag co-expressed with both α1 and β2 traffics to the cell surface



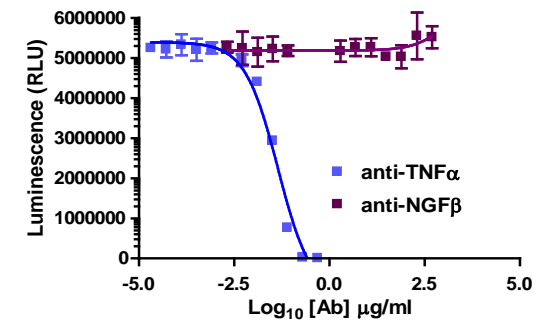
GABAAR pharmacology is not affected

NF- κ B-Response Element Reporter Gene Assay

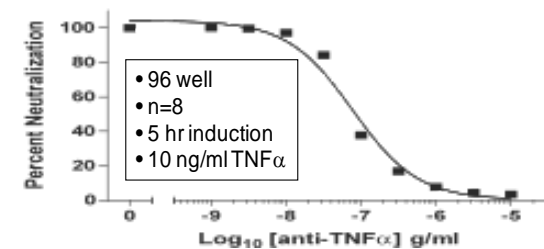
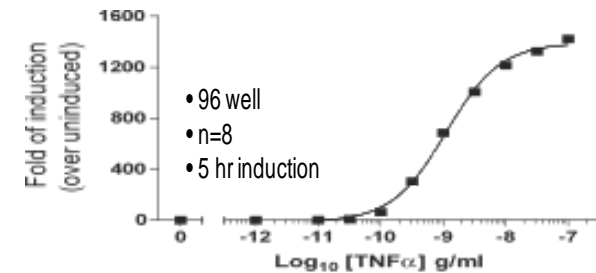


8935MB

Specificity



TNF α /anti-TNF α bioactivity

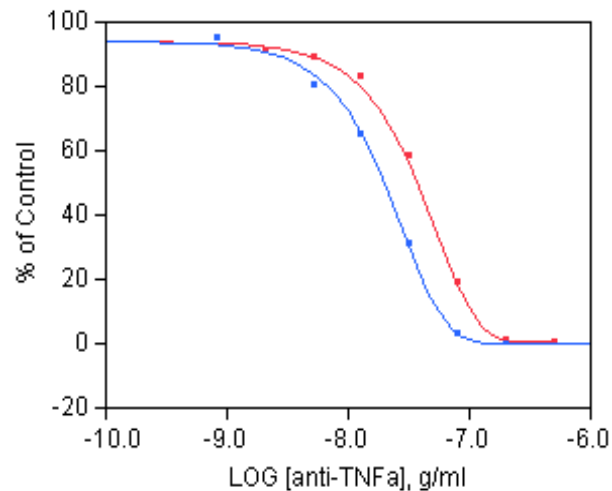


Parallelism & Linearity Across Potencies



Anti-TNF α Bioassay Using NF- κ B-RE Luciferase Assay

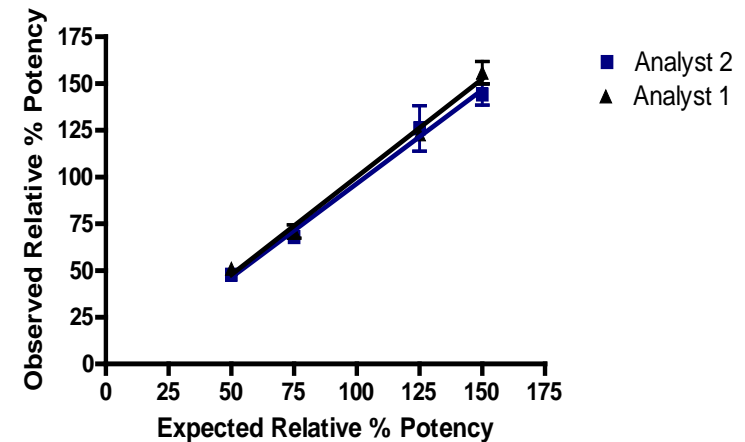
Example: Parallelism is observed for a 50% potency sample relative to a 100% reference sample



Relative Potency

Parameter	Estimate
Log Standard EC50	-7.6939
Log Unknown EC50	-7.4249
Relative Potency	0.53823
Parallel Check (p > F)	0.91758

Linearity is evident for the potency range



Parallelism analysis performed using JMP software (SAS Institute)

Accuracy of Bioluminescent Anti-TNF α Bioassay



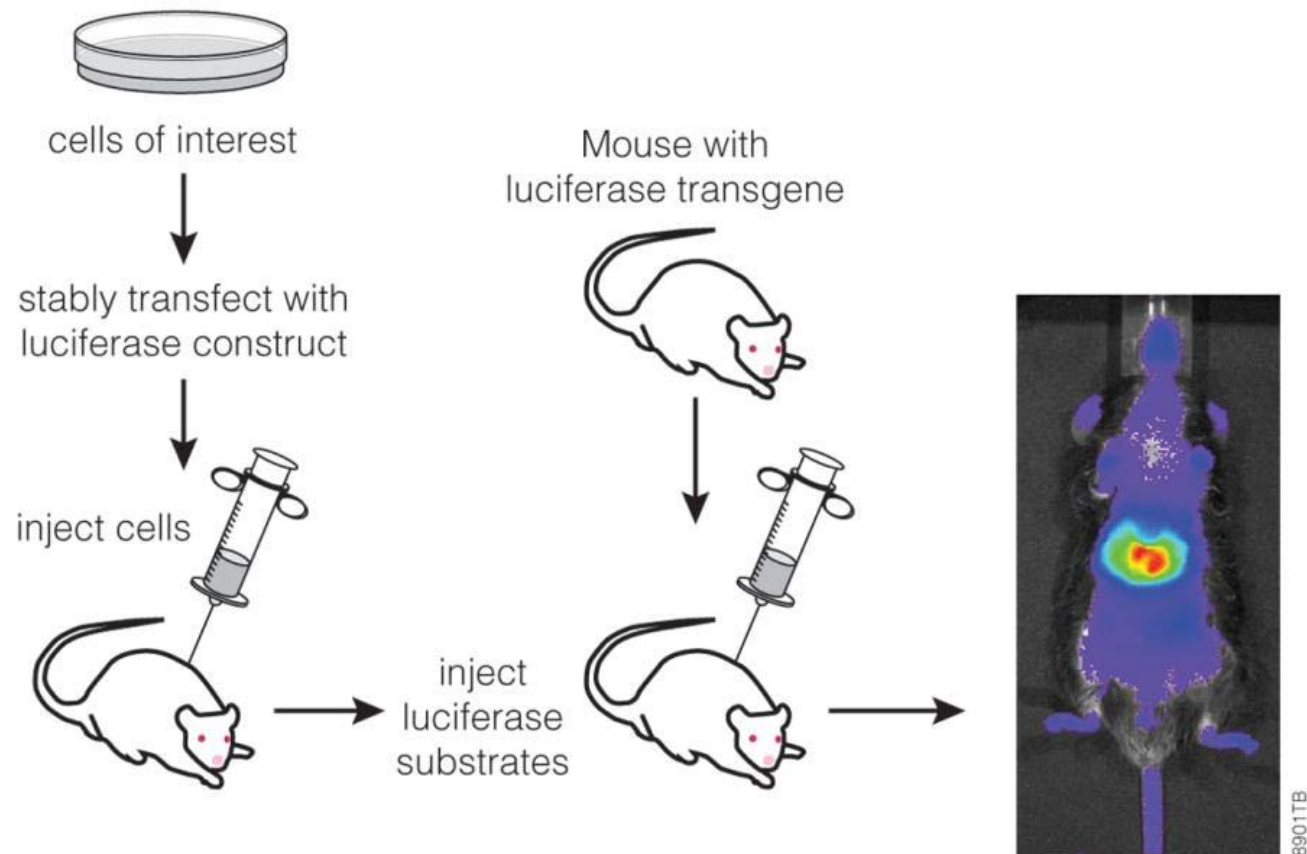
NF- κ B-RE reporter gene assay

Potency sample	Recovery Analyst 1	Recovery Analyst 2
50%	101.8%	95.4%
75%	94.6%	90.5%
125%	98.2%	100.8%
150%	103.9%	96.1%

Data are means from 3 separate assays performed on 3 different days

From Cells to In Vivo Imaging

Luciferase technology is ideal for carrying assays into live animals



CASE STUDIES: Solutions for Small Molecule & Biologics Drug Discovery



CASE STUDY: Custom Stable Cell Line Need for Small Molecule Drug Discovery Program



A large pharmaceutical company was having difficulty developing a stable cell line in a specific cellular background due to technical issues.

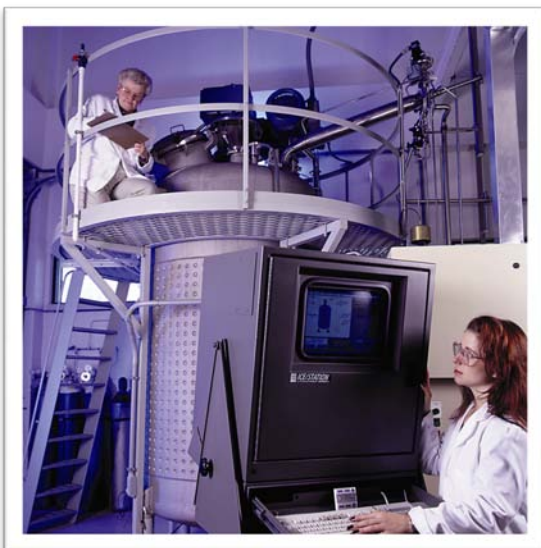
Problem: Difficulty developing a stable cell; needed a cell line with sensitive readout for specific cell signal response

Proposal: Create and validate HaloTag-containing stable cell line

Solution: Validated stable cell line in 12 weeks

Result: Fewer resources expended by Client; successful assay when previous efforts failed; Client spent less in saved resources and effort in not troubleshooting assay development.

CASE STUDY: Custom Bioassay Need for Drug Discovery in Biologics



A global pharmaceutical company needed a more sensitive and easier-to-perform bioassay for pre-clinical and clinical phases of new drug development.

Problem: Laborious, poorly sensitive assay to support potential blockbuster drug in clinical trials and post-launch

Proposal: Create a qualified luciferase reporter assay including training & transfer of assay to Client

Solution: Qualified potency bioassay developed in 21 weeks

Result: More sensitive & easier to perform lot release assay; Client not experienced in developing cell-based assays and concentrated on other aspects of the project while Promega developed the assay solution.

Strategic Collaborations Team



Eric Muhr (Europe/ROW)

Strategic Collaborations Director

Promega Corporation

Cell: 908-391-5774

eric.muhr@promega.com

Mike Bodnaruk (US East & South)

Strategic Collaborations Manager

Promega Corporation

Office: 617-276-2432

Cell: 617-276-2432

michael.bodnaruk@promega.com

Richard Somberg, PhD (US Central & West)

Strategic Collaborations Manager

Promega Corporation

Office: 608-298-4864

Cell: 608-358-1589

richard.somberg@promega.com

The Strategic Collaborations Team is supported by a network of professionals within Promega. From the experienced R&D staff, to Custom Order Specialists, to business and legal management, we're committed to meeting your needs and exceeding your expectations.

General Inquiries:

cas@promega.com

Worldwide Support



Headquarters:

Promega Corporation

2800 Woods Hollow Road
Madison, WI 53711-5399 USA
Tel: 608-274-4330
Fax: 608-277-2516
Toll-Free: 800-356-9526
Toll-Free Fax: 800-356-1970



www.promega.com

Manufacturing:

Promega BioSciences, Inc.

San Luis Obispo, CA, USA

Promega BioSystems, Inc.

Seoul, South Korea

Promega BioSystems

Sunnyvale, Inc.

Sunnyvale, CA, USA

Shanghai Promega Biological Products, Ltd.

Shanghai, China

Terso Solutions, Inc.

Madison, WI, USA

Branch Offices:

Promega Corporation

Sydney, Australia

Promega (Beijing)

Biotech Co., Ltd

Beijing, China

Promega France

Lyon, France

Promega GmbH

Mannheim, Germany
(Austria)

Promega Italia, S.r.l.

Milan, Italy

Promega KK

Tokyo, Japan

Promega Korea, Ltd

Seoul, South Korea

Promega Benelux b.v.

Leiden, The Netherlands
(Belgium/Luxembourg)

Promega Pte Ltd.

Pacific Asia Region, Singapore

Promega Biotech Ibérica, SL

Madrid, Spain

Promega Biotech AB

Stockholm, Sweden
(Denmark)

Promega AG

Dübendorf, Switzerland

Promega UK

Southampton, United Kingdom



PROMEGA CORPORATION • 2800 WOODS HOLLOW ROAD • MADISON, WI 53711-5399 USA • TELEPHONE 608-274-4330

www.promega.com • ©2010 PROMEGA CORPORATION • ALL RIGHTS RESERVED • PRICES AND SPECIFICATIONS SUBJECT TO CHANGE WITHOUT PRIOR NOTICE • PRINTED IN USA 6/10 • 18482-061110

