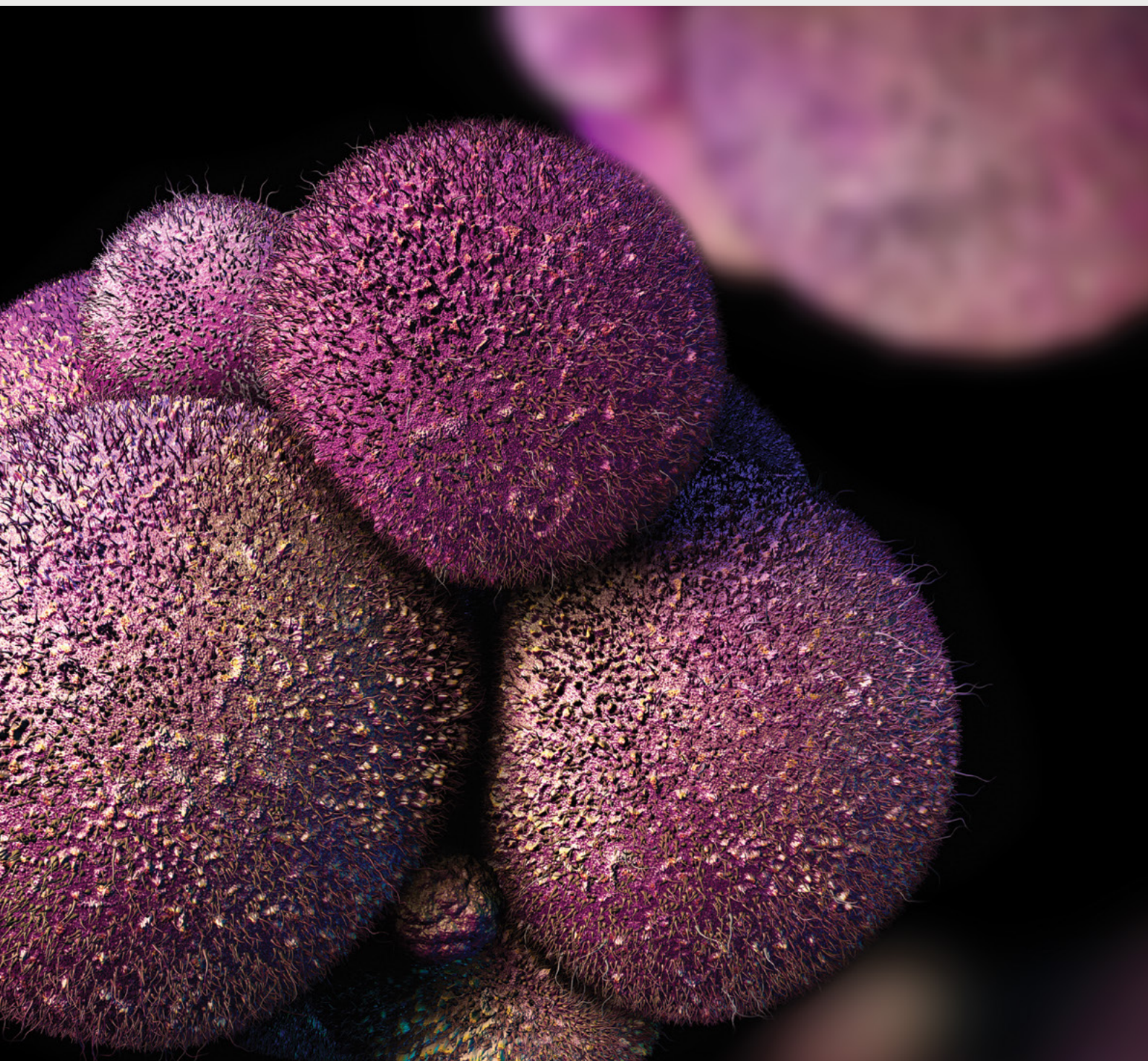




# Stem Cells

Integrated Solutions for  
Stem Cell Research



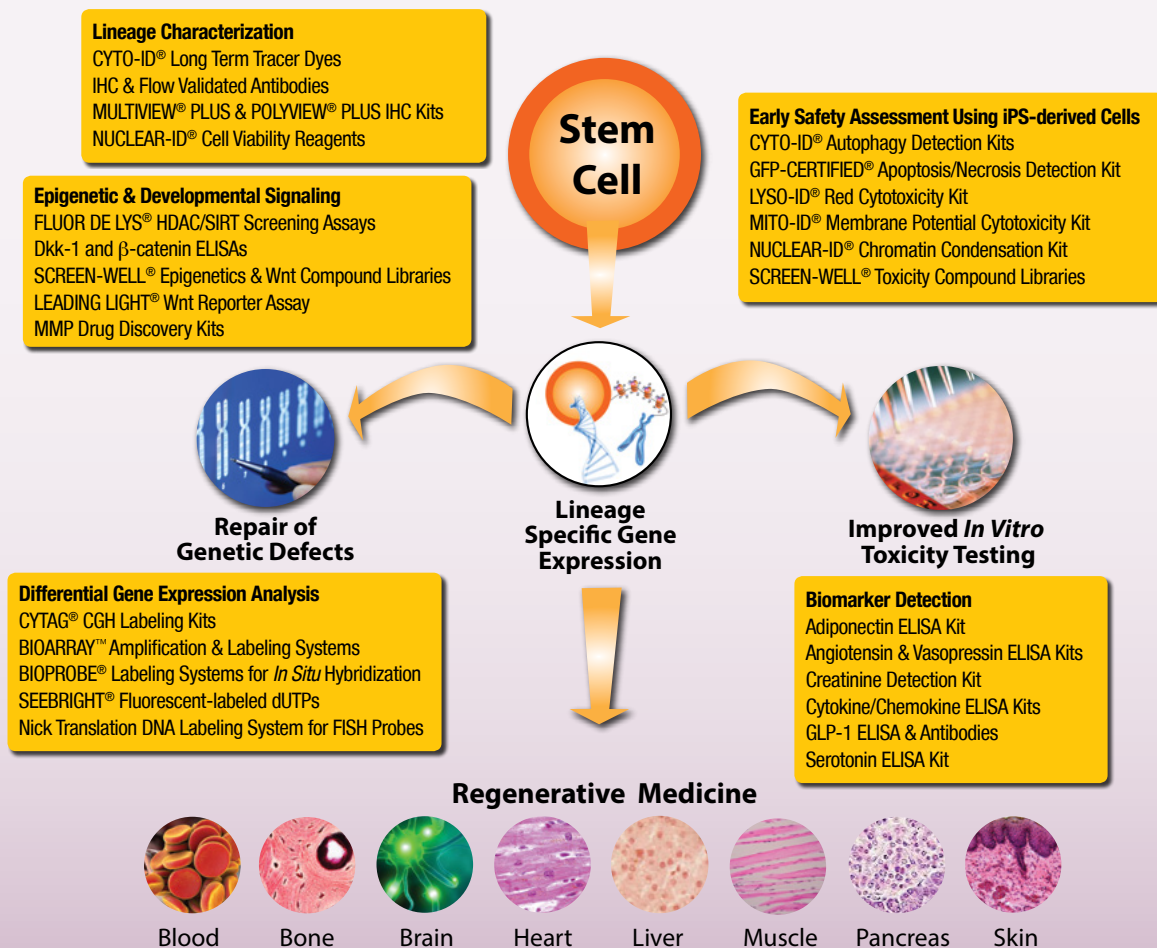


# DIFFERENTIATE YOUR RESEARCH

## ON THE PATH TO STEM CELL DISCOVERY

### Therapeutic Utility & Promise of Stem cells

Regardless of the specific path your stem cell research takes, we provide innovative tools to help you realize its full potential. Pluripotent stem cells are a renewable source of replacement cells which have the potential to repair damaged tissues or treat degenerative diseases without invoking immune rejection. With this, they hold the potential for treating a myriad of pathologies including cancer, diabetes, hematological disorders, cardiovascular, inflammatory and neurodegenerative diseases. Unlocking the cellular mechanisms that regulate self-renewal, reprogramming and differentiation holds promise for new cellular therapies, regenerative medicine, and improving model systems to assess drug efficacy and toxicity.



### Biochemical Assays

Detection of protein activity related to transcriptional regulation and cell signaling is critical to understanding the mechanisms associated with stem cell pluripotency and differentiation. Our innovative assays in the area of epigenetics and Wnt signaling provide stem cell research with robust screening assays, featuring unrivaled sensitivity and specificity for investigating these regulatory pathways in stem cell biology.

### Nucleic Acid Labeling and Amplification

We are a recognized pioneer and innovator of life sciences tools, backed by patented DNA and RNA labeling chemistries for genomics research and development. Our unique offering of kits for genomic analysis enables monitoring the genomic integrity of long-term stem cell cultures or to assess expression of genes regulating stem cell differentiation and fate.

### Cell Analysis & Detection Reagents

Our expertise in fluorescent probe chemistry and cellular analysis combine to provide high-quality, flexible detection platforms for stem cell profiling and functional analysis. Our unique CELLESTIAL® assays and widely cited antibodies enable multiplex detection of cellular markers and key parameters of cell viability, cell signaling, death pathways, and toxicity.

# MODULATE STEM CELL FATE WITH SMALL MOLECULES & PROTEINS

Stem cell self-renewal and differentiation involves complex events which lead to the generation of various progeny cells and tissues. We offer an extensive collection of small molecules and proteins that can be used to modulate differentiation, biochemical activity, and cell growth.

## Small Molecules

Differentiation	Product #
Compound E	ALX-270-415
Cyclopamine	BML-GR334
DAPT	ALX-270-416
Dexamethasone	BML-EI126
DIF-1	BML-GR324
Forskolin	BML-CN100
IDE-1	BML-SC100
IDE-2	BML-SC101
Purmorphamine	ALX-420-045
Retinoic acid, all trans	BML-GR100
Sodium butyrate	ALX-270-301
Valproic acid	ALX-550-304

Proliferation & Self-Renewal	Product #
BI-D1870	BML-EI407
BML-275 (Dorsomorphin)	BML-EI369
ID-8	BML-SC105
XAV939	BML-WN100
Y-27632 . Dihydrochloride	ALX-270-333

Re-programming	Product #
ALK5 inhibitor I	ALX-270-448
AM-580	BML-GR104
BIX01294	ALX-270-517
Doxycycline hyclate	ALX-380-273
Pifithrin- $\alpha$ (cyclic)	ALX-270-287
Pifithrin- $\alpha$	BML-GR325
Pifithrin- $\mu$	BML-AP503
Tranylcypromine	BML-EI217

SCREEN-WELL® Compound Libraries	Product #
Cardiotoxicity Library	BML-2850
Epigenetics Library	BML-2836
FDA Approved Drugs Library	BML-2843
Hematopoietic Toxicity Library	BML-2852
Hepatotoxicity Library	BML-2851
Myotoxicity Library	ENZ-LIB101
Nephrotoxicity Library	ENZ-LIB100
Stem Cell Library	ENZ-LIB103
Wnt Pathway Library	BML-2838

## Proteins

Growth Factors	Product #
Activin-A (recombinant)	ALX-201-809
Axin 1/2 (recombinant)	ENZ-PRT130/-131
BMP-4 (recombinant)	ENZ-PRT146
Dkk-1 (recombinant) (His)	ADI-KPR-CP100
EGF (recombinant)	ALX-201-812
Flt-3 Ligand (recombinant)	ALX-201-800
Follistatin (recombinant)	ALX-201-813
GDF-3 (recombinant)	ENZ-PRT176
GDF-5 (recombinant)	ALX-201-802
KGF (recombinant)	ENZ-PRT182
m-CSF (recombinant)	ENZ-PRT143
Noggin (recombinant)	ENZ-PRT127
Stem Cell Factor (recombinant)	ALX-201-804
Stem Cell Factor (recombinant)	ALX-201-805
Stem Cell Factor (recombinant)	ALX-201-806
Wnt-3a (recombinant)	ENZ-60001

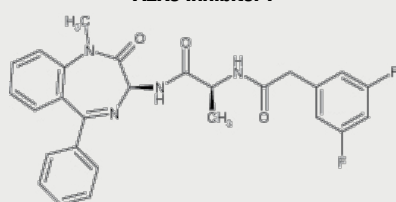
Cytokines/Chemokines	Product #
FGF-4 (recombinant)	ENZ-PRT173
FGF-8 (recombinant)	ENZ-PRT174
FGF basic (recombinant)	ALX-201-824
FGF-basic 147 (recombinant)	ALX-201-822
FGF-basic 154 (recombinant)	ALX-201-823
IL-3 (recombinant)	ALX-201-820
IL-3 (recombinant)	ALX-201-821
IL-6 (recombinant)	ALX-201-818
IL-6 (recombinant)	ALX-201-819
IL-12 (recombinant)	ALX-201-815
TGF- $\beta$ 3 (recombinant)	ALX-201-817

Stem Cell Markers	Product #
Thy-1 (Fc) (recombinant)	ALX-522-091
Thy-1 (Fc) (recombinant)	ALX-522-096

Cell Culture Tools	Product #
Collagen 1, rat tail (thin plate coating)	ALX-522-440
Fibrinogen (plasma)	ENZ-PRT157
Laminin (fibroblasts)	ENZ-PRT159
Vitronectin (recombinant)	ENZ-PRT179

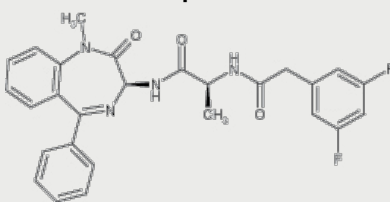
Human (H) Rat (R) Mouse (M)

ALK5 Inhibitor I



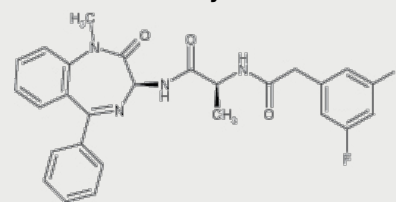
TGF- $\beta$  RI KINASE INHIBITOR (ALX-270-448)

Compound E



$\gamma$ -SECRETASE INHIBITOR (ALX-270-415)

Y-27632 . dihydrochloride



ROCK INHIBITOR (ALX-270-333)

# ACHIEVE YOUR STEM CELL POTENTIAL

## LIVE CELL ANALYSIS

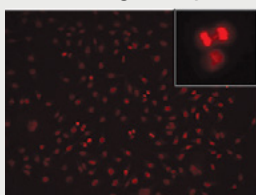
### MULTIPLEX ANALYSIS OF CELL FATE AND DIFFERENTIATION BY FLOW CYTOMETRY

Our fluorescent probe-based CELLESTIAL® assays and high-specificity antibodies enable multiplex characterization of stem cell populations, including real-time functional assays to assess cell viability, cell death, and toxicity by flow cytometry. Our GFP-CERTIFIED® dyes enable simultaneous functional analysis in GFP-expressing cells or with other green-emitting dyes/labels.

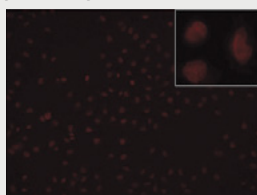
#### NUCLEAR-ID® Red Cell Cycle Kit (GFP-CERTIFIED®) (ENZ-51008)

- Brighter, more stable red fluorescence yields improved sensitivity
- Highly cell permeable dye minimizes cell density optimization
- Suitable for multiplexing with green fluorophores, including GFP and FITC
- Available as a complete kit with controls or as a stand-alone dye

##### Brightest Dye for Cell Cycle Analysis



NUCLEAR-ID® Dye (2.5 µM)

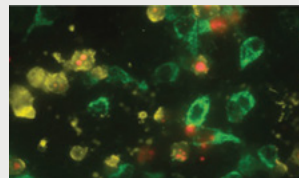


Competitor Dye (2.5 µM)

#### GFP-CERTIFIED® Apoptosis/Necrosis Detection Kit (ENZ-51002)

- True multiplexing capabilities with GFP and other green fluorescent probes
- Optimized for both fluorescence microscopy and flow cytometry applications
- Assess effects of stem cell isolation and preservation techniques, drug treatment effects, reprogramming modulators

##### Detect Cell Death in GFP-Expressing Cell Lines



GFP-compatible multiplex assay for distinguishing between healthy, early apoptotic, late apoptotic and necrotic cells.

## ANTIBODIES

### IDENTIFY STEM CELL POPULATIONS WITH MARKERS OF DIFFERENTIATION

We offer a broad range of antibodies for use in all areas of stem cell research. Many are validated for use in key applications such as flow cytometry, immunohistochemistry, immunoprecipitation and Western blotting. Every antibody is backed by our **Worry-Free Antibody Trial Program**, allowing you to evaluate any of our antibodies of interest for your specific application or species without risk.

Stem Cell Markers	Embryonic	Mesenchymal	Hematopoietic	Endothelial Progenitors	Neural Progenitors	Cancer
β1-Integrin	•					
βIII-Tubulin					•	
BCRP	•		•			
CD9	•					
CD14		•		•		
CD31				•		•
CD40		•				
CD44		•				
CD45		•		•		
CD49b/CD29	•					
c-Kit			•			
c-Myc	•					
CXCR4			•			
Dnmt3b	•					
E-cadherin	•					
eNOS				•		
EpCAM						•
Galanin	•					
GFAP					•	
NCAM					•	
Neutrophil Elastase			•			
NGFR		•				
Oct-3/4	•					
PSCA						•
Serpin			•			

## FOCUSED COMPOUND LIBRARIES FOR SCREENING

### ESSENTIAL STANDARDS FOR STEM CELL REGULATORY AND SIGNALING PATHWAY SCREENING

We have a long and successful history assembling focused libraries for chemical genomics, receptor de-orphaning, drug re-purposing, pathway analysis, and more.

#### SCREEN-WELL® Stem Cell Library (ENZ-LIB103)

- Contains 130 compounds available in a 100 µL/well format
- Compounds can be used for diverse stem cell research to investigate cellular mechanisms that regulate self-renewal, reprogramming, and differentiation
- We provide detailed information in our comprehensive documentation for each compound
- Bulk compounds available for re-supply

## GENOME ANALYSIS

### MAXIMIZE GENOMIC DATA WITH MINIMAL SAMPLE INPUT

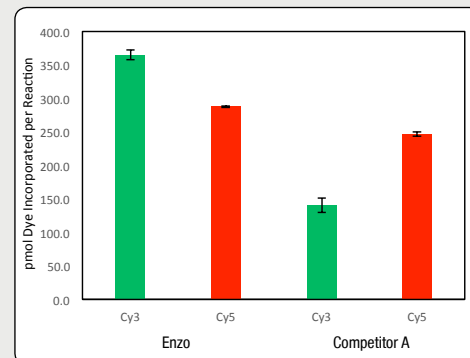
Detect DNA copy number gains and losses associated with chromosome abnormalities.

#### CYTAG® SuperCGH Labeling Kit (ENZ-GEN120)

- Perform total genomic DNA analysis without amplification or complexity reduction
- Increase resolution for comprehensive, unbiased analysis of DNA copy number changes
- Produce superior DLR scores below 0.12

DNA samples with 50ng input were labeled with Enzo's CYTAG® SuperCGH Labeling Kit for Oligo Arrays or a leading competitor's kits. Enzo's proprietary labeling technology generates a superior dye incorporation.

#### CYTAG® SuperCGH Provides Superior Dye Incorporation



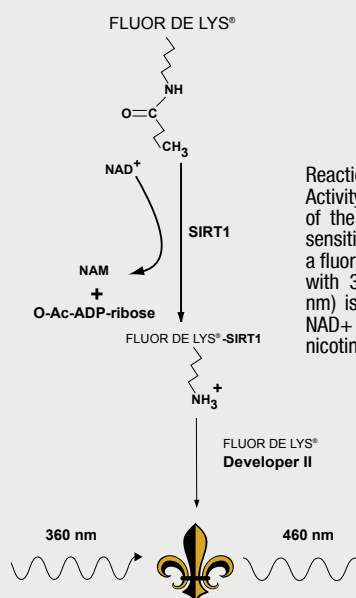
## BIOCHEMICAL & REPORTER ASSAYS

### DISCOVER NOVEL EPIGENETIC AND WNT PATHWAY MODULATORS

Our assays for screening modulators of HDAC/SIRT activity and the Wnt pathway offer the sensitivity, specificity, and reproducibility researchers need to reveal the epigenetic fingerprints and developmental signals regulating stem cell maintenance and self-renewal.

#### FLUOR DE LYS® SIRT1 Fluorometric Drug Discovery Assay Kit (BML-AK555)

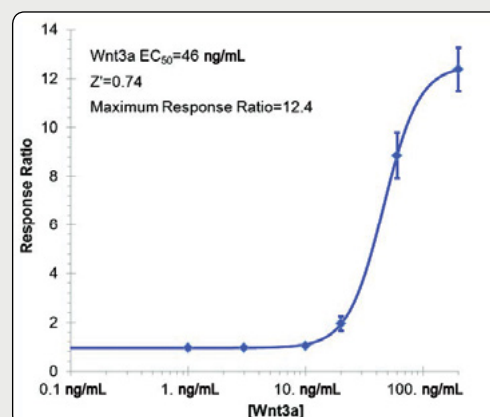
- Useful for inhibitor screening or characterizing enzyme kinetics
- Includes optimal substrate selected from a panel of acetylated sites in p53 and histones
- Supplied with enough recombinant enzyme for 96 assays



#### LEADING LIGHT® Wnt Reporter Assay (ENZ-61001)

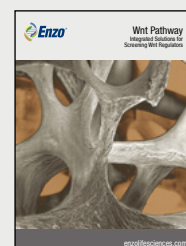
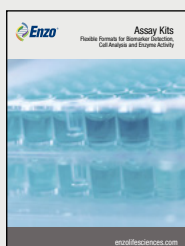
- No transfection required
- Luciferase-based assay accelerates screening analysis of Wnt pathway effectors
- Highest sensitivity in response to Wnt effectors without the use of lithium chloride
- Highly active Wnt-3a protein included to ensure robust signal 96- or 384-well format allows high-throughput analysis of Wnt pathway effectors

#### Dose-dependent Activation by Wnt-3a Protein



Luciferase activity in the LEADING LIGHT® Wnt reporter cell line increases in a dose-dependent manner in response to treatment with Wnt-3a protein without the need for LiCl.

***You may also be interested in:***



**Global Headquarters**  
**ENZO LIFE SCIENCES, INC.**  
 10 Executive Blvd.  
 Farmingdale, NY 11735  
 Ph: 800.942.0430  
 Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

**European Sales Office**  
**ENZO LIFE SCIENCES (ELS) AG**  
 Industriestrasse 17  
 CH-4415 Lausen, Switzerland  
 Ph: +41 61 926 8989  
 Fax: +41 61 926 8979  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

## LOCAL EUROPEAN OFFICES

**Belgium, The Netherlands & Luxembourg**  
 Enzo Life Sciences BVBA  
 Avenue Louise 65/Box 11  
 1050 Bruxelles  
 Belgium  
 Ph: +32 3 466 0420  
 Fax: +32 3 808 7033  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

**France**  
 Enzo Life Sciences (ELS) AG  
 Branch Office Lyon  
 13, avenue Albert Einstein,  
 F-69100 Villeurbanne, France  
 Ph: +33 472 440 655  
 Fax: +33 481 680 254  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

**Germany**  
 Enzo Life Sciences GmbH  
 Basler Strasse 57a  
 DE-79540 Lörrach  
 Germany  
 Ph: +49 7621 5500 526  
 Fax: +49 7621 5500 527  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

**UK & Ireland**  
 Enzo Life Sciences (UK) Ltd.  
 1 Colleton Crescent  
 Exeter EX2 4DG  
 Ph: 0845 601 1488 (UK customers)  
 Ph: +44 1392 825900  
 Fax: +44 1392 825910  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

For local distributors and detailed product information visit us online:  
**[www.enzolifesciences.com](http://www.enzolifesciences.com)**

**Put our experience to work for you!**

Our broad range of scientific expertise and manufacturing capabilities enables us to provide a comprehensive set of solutions for stem cell research.