## **Product Manual**

# StemTAG™ Alkaline Phosphatase Staining Kit (Purple)

**Catalog Number** 

**CBA-306** 

100 assays

FOR RESEARCH USE ONLY Not for use in diagnostic procedures



### **Introduction**

Embryonic stem (ES) cells are continuous proliferating stem cell lines of embryonic origin first isolated from the inner cell mass (ICM). Two distinguishing features of ES cells are their ability to be maintained indefinitely in an undifferentiated state and their potential to develop into any cell within the body. Based on previous methods developed for mouse ES cells, human ES cell lines were first established by Dr. James Thomson and colleagues. Like mouse ES cells, human ES cells express high levels of membrane alkaline phosphatase (AP) and Oct-4, a transcriptional factor critical to ICM and germline formation. However, unlike mouse ES cells, hES cells do not express stage-specific embryonic antigen (SSEA-1). In addition, prolonged propagation of hES cells is typically achieved by coculture with primary mouse embryonic fibroblasts (MEFs) serving as feeder cells. Human ES cell lines are not able to maintain their undifferentiated state in the absence of supporting feeder layer cells, even when exogenous cytokines such as leukemia inhibitory factor (LIF) and gelatin-coated plates are used.

Marker	Mouse ES	Mouse EG	Human ES	Human EG	Human EC
Name	Cells	Cells	Cells	Cells	Cells
AP	$\sqrt{}$	$\sqrt{}$	$\sqrt{}$		$\sqrt{}$
SSEA-1		$\sqrt{}$	_		1
SSEA-4	_	_	$\sqrt{}$		$\checkmark$
TRA-1-					
60	_	1	$\sqrt{}$		$\sqrt{}$
TRA-1-					
81	_	1	$\sqrt{}$		$\sqrt{}$
Oct-4		$\sqrt{}$	$\sqrt{}$	unknown	$\sqrt{}$
ES Cell = Embryonic stem cell					
EG Cell = Embryonic germ cell					
EC Cell = Embryonic carcinoma cell					

Table 1. Comparison of Mouse and Human Pluripotent Stem Cells.

Although stem cells from different origins require different growth conditions for self-renewal and display different cell surface markers (see Table 1), AP is the most widely used stem cell marker. The StemTAG<sup>TM</sup> Alkaline Phosphatase Staining Kit provides an efficient system for monitoring ES cell undifferentiation/differentiation through AP activity by immunocytochemistry staining.

# **Related Products**

- 1. CBA-300: StemTAG<sup>TM</sup> Alkaline Phosphatase Staining Kit (Red)
- 2. CBA-301: StemTAG<sup>TM</sup> Alkaline Phosphatase Activity Assay Kit (Colorimetric)
- 3. CBA-312: MEF Feeder Cells (Puromycin-resistant)
- 4. CBA-316: SNL Feeder Cells
- 5. CBA-320: CytoSelect<sup>TM</sup> 96-Well Hematopoietic Colony Forming Cell Assay

# **Kit Components**

- 1. Fixing Solution (Part No. 30001): One bottle 50 mL
- 2. StemTAG<sup>TM</sup> AP Staining Solution (Part No. 130601): One amber bottle 40 mL



# **Materials Not Supplied**

- 1. Human or Mouse Embryonic Stem Cells and Culture Medium
- 2. 1X PBS
- 3. 1X PBST (1X PBS containing 0.05% Tween-20)

#### Storage

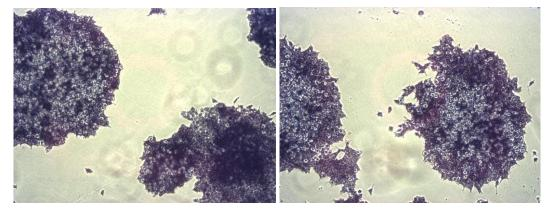
Store all components at 4°C.

## **Assay Protocol (24-Well Plate)**

- 1. Culture mouse ES cells in medium containing LIF; alternatively, culture human ES cells on a MEF feeder layer.
- 2. Gently aspirate the medium from the ES cells and wash the cells with 1 mL of 1X PBST. Aspirate the wash solution.
- 3. Add Fixing Solution to the cells, 0.4 mL per well for a 24-well plate. Incubate at room temperature for 2 minutes.
- 4. Remove the fixing solution and wash the fixed cells twice with 1 mL of 1X PBST.
- 5. Aspirate the final wash, and add 0.4 mL per well of StemTAG<sup>TM</sup> AP Staining Solution
- 6. Incubate the cells at room temperature for 15-30 minutes, protected from light.
- 7. Remove the AP Staining Solution, and then wash the stained cells twice with 1 mL of 1X PBS. Store cells in 1X PBS at 4°C. For long-term storage, overlay the cells with 1X PBS containing 20% Glycerol. Store at 4°C.
- 8. Count the purple stained cell colonies (undifferentiated ES cells) vs. colorless colonies (differentiated ES cells) using a light microscope.

# **Example of Results**

The following figures demonstrate typical results with the StemTAG<sup>™</sup> Alkaline Phosphatase Staining Kit (Purple). One should use the data below for reference only. This data should not be used to interpret actual results.



**Figure 1: AP staining of ES Cells.** Murine embryonic stem cells (ES-D3) are maintained in an undifferentiated stage on gelatin-coated dishes in the presence of LIF. AP staining of undifferentiated ES cells was performed as described in the Assay Protocol.



### References

- 1. Wobus AM, Holzhausen H, Jäkel P et al. (1984). Exp Cell Res 152:212–219.
- 2. Thomson JA, Itskovitz-Eldor J, Shapiro SS et al. (1998). Science 282:1145–1147.
- 3. Smith AG, Nichols J, Robertson M et al. (1992). Dev Biol 151:339–351.
- 4. Reubinoff BE, Pera MF, Fong CY et al. (2000). *Nat Biotechnol* **18**:399–404.

### **Recent Product Citations**

- 1. Kim, H. et al. (2022). Bioactive Surface of Zirconia Implant Prepared by Nano-Hydroxyapatite and Type I Collagen. *Coatings*. **12**(9):1335. doi: 10.3390/coatings12091335.
- 2. Mitchell, M.W. et al. (2022). An induced pluripotent stem cell line (CIMRi001-A) from a Vici syndrome donor with a homozygous recessive c.1007A>G (p.Q336R) mutation in the EPG5 gene. *Stem Cell Res.* doi: 10.1016/j.scr.2022.102833.
- 3. Chan, C.D. et al. (2021). Co-localisation of intra-nuclear membrane type-1 matrix metalloproteinase and hypoxia inducible factor-2α in osteosarcoma and prostate carcinoma cells. *Oncol Lett.* **21**(2):1-7. doi: 10.3892/ol.2020.12419.
- 4. Oh, J.E. et al. (2020). Isolation and characterization of dental follicle-derived Hertwig's epithelial root sheath cells. *Clin Oral Investig*. doi: 10.1007/s00784-020-03481-4.
- 5. Zhang, Y. et al. (2020). SENP3 Suppresses Osteoclastogenesis by De-conjugating SUMO2/3 from IRF8 in Bone Marrow-Derived Monocytes. *Cell Rep.* **30**(6):1951-1963.e4. doi: 10.1016/j.celrep.2020.01.036.
- 6. Li, N. et al. (2020). miR-144-3p suppresses osteogenic differentiation of bone marrow mesenchymal stem cells from patients with aplastic anemia through repression of TET2. *Mol Ther Nucleic Acids*. doi: 10.1016/j.omtn.2019.12.017.
- 7. Khan, M.I. et al. (2016). Comparative gene expression profiling of primary and metastatic renal cell carcinoma stem cell-like cancer cells. *PLoS One* **11**:e0165718.

### Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

# **Contact Information**

Cell Biolabs, Inc. 5628 Copley Drive San Diego, CA 92111

Worldwide: +1 858-271-6500 USA Toll-Free: 1-888-CBL-0505 E-mail: tech@cellbiolabs.com

www.cellbiolabs.com

©2012-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

