

SUMMARY

Mouse embryonic stem (ES) cells are permanent lines of undifferentiated pluripotent cells derived from early mouse embryos. They exhibit two important properties: the ability to self-renew and give rise to new pluripotent ES cells, and the ability to differentiate into all specialized cell types found in the adult mouse. Research in the 1980's and 1990's focused on combining techniques for manipulating the mouse genome with ES cell technology to produce chimeric and knock-out mice for evaluating the function of mammalian gene products *in vivo*. More recently, advances in controlling ES cell *in vitro* differentiation are providing insights into the mechanisms involved in embryonic development, tissue lineage specification and the generation of cells for potential therapeutic applications.

CHARACTERISTICS AND CULTURE OF MOUSE ES CELLS

ES cells are isolated from the inner cell mass of pre-implantation embryos or blastocysts at day 3.5 of mouse development. When mouse blastocysts are cultured on a feeder layer of mitotically inactivated embryonic fibroblasts, the outer layer of cells attach and undifferentiated cells from the inner cell mass (ICM) spontaneously form clumps that can be further passaged to yield ES cell lines.^{1,2} These cells are considered pluripotent as they can be maintained indefinitely in the undifferentiated state in culture and have the ability to contribute to all of the tissues, including the germ cells, when injected back into mouse blastocysts.³ Maintenance of ES cells *in vitro* is achieved by co-culture on irradiated mouse fibroblasts² or on gelatinized dishes with a differentiation inhibitory factor called Leukemia Inhibitory Factor (LIF).⁴

In addition to LIF, a number of other extrinsic and intrinsic regulators of self-renewal have been identified (for review, see Chambers and Smith).⁵ It has recently been reported that Bone Morphogenetic Protein 4 (BMP-4) can replace the requirement of serum during ES cell line derivation and maintenance through suppression of neural differentiation.⁶ The secreted Wnt proteins may also play a role in maintaining pluripotency, as Wnt signaling is down-regulated upon differentiation.^{7,8}

The transcription factor Oct-4 (Oct-3, Oct-3/4, Pou5f1) is also required for regulation of cell fate in the early embryo; it is expressed in the ICM and down-regulated upon differentiation into trophoblast.⁹ Induction of differentiation through withdrawal of LIF results in concomitant reduction in Oct-4 levels.⁵ A second homeodomain-containing protein, Nanog, is expressed in the inner cells of a compacted morula and blastocyst (as well as early germ cells) and is down-regulated just prior to implantation. Nanog overexpression can maintain self-renewal of ES cells independently of the LIF/gp130/Stat3 pathway.^{10,11} Other factors proposed to have a role in maintaining "stemness" include Sox2 (SRY-related HMG box 2) and FoxD3 (forkhead box D3).⁵ Appropriate culture

conditions and qualified reagents also play an integral part in retaining the pluripotency of ES cells.¹²

MOUSE ES *IN VITRO* DIFFERENTIATION INTO EMBRYOID BODIES

Upon withdrawal of LIF and stromal contact, mouse ES cells will spontaneously differentiate into complex, three dimensional cell aggregates called embryoid bodies (EBs). Differentiation within EBs occurs in a well-defined temporal manner with the initial formation of all three embryonic germ layers (ectoderm, mesoderm, endoderm) followed by further differentiation to terminally differentiated cell types, similar to *in vivo* embryogenesis.^{12,13} Although a variety of cell types form during spontaneous differentiation, a heterogeneous mixture of cells are present and not all cell lineages (i.e. endoderm-derived) have been observed.

ES *IN VITRO* DIFFERENTIATION INTO DIFFERENT TISSUE CELL TYPES

Protocols have been developed to enrich and further expand the repertoire of tissue cell types generated following *in vitro* differentiation of mouse ES cells. A summary is presented in Table 1 (page 191). An enrichment strategy, termed genetic selection has been used to express specific fate-determining genes (e.g., *HOXB4* transcription factor implicated in definitive hematopoiesis¹⁴) or drug resistance markers under the control of lineage-specific promoters¹⁵⁻¹⁷ in ES cells to generate relatively homogenous populations of mature cell types. For example, Marchetti *et al.* used the vascular endothelium specific promoter Tie-1 linked to an antibiotic resistant gene to derive a purified population of endothelial cells from ES cells. These cells could be selectively expanded in culture and following transplantation were incorporated into the neovasculature of nude mice.¹⁵

Selection of desired cell types can also be achieved through cell sorting on the basis of one or more specific cell surface antigens. Yamashita *et al.* used this technique to isolate ES cell-derived Flk1⁺ hematopoietic progenitors by fluorescent activated cell sorting (FACS). These cells were shown to act as vascular precursor cells capable of differentiating *in vitro* and *in vivo* into both endothelial and mural cells of the vasculature.¹⁸ Therefore, cell enrichment strategies also facilitate the characterization of precursor cells, providing new insights into the cellular origins of cells in vascular development.

Selective media and culture conditions are also widely used to select a desired cell population and have been instrumental in the enrichment of neural cell types.¹⁹⁻²¹ Brustle *et al.* used this approach to generate a population of multipotent progenitor cells that express glial precursor markers. Mouse ES cells were cultured in a medium that favors survival of neural precursors followed by



StemCell Technologies
www.stemcell.com

In North America
Tel: 604.877.0713
Fax: 604.877.0704
Toll Free Tel: 1.800.667.0322
Toll Free Fax: 1.800.567.2899
e-mail: info@stemcell.com
e-mail: orders@stemcell.com

In the United Kingdom
Tel: +44.(0).20.7537.7565
Fax: +44.(0).20.7515.5408
Toll Free within United Kingdom:
Tel: 0800.731.27.14
Fax: 0800.731.27.13
e-mail: info@stemcellgb.com
www.stemcellgb.com

In Europe
Tel: +33.(0).4.76.04.75.30
Fax: +33.(0).4.76.18.99.63
e-mail: info@stemcellfrance.com

the sequential addition of various growth factor combinations [fibroblast growth factor-2 (FGF-2), FGF-2 and epidermal growth factor (EGF), FGF-2 and platelet-derived growth factor (PDGF)], resulting in cells that could be maintained for several generations in culture. Upon growth factor withdrawal, these cells differentiated into either of two specific lineages, oligodendrocytes or astrocytes.²¹ Specific combinations of cytokines have also been used to promote the growth of other mature cell types including mast cells²², hematopoietic cells²³ and dendritic cells.²⁴ For instance, using a combination of granulocyte macrophage colony-stimulating factor (GM-CSF) and interleukin-3, Fairchild and colleagues generated long-term, stable cultures of dendritic cells from differentiating ES cells.²⁴ Chemical inducers, such as the Vitamin A derivative-retinoic acid, have also been widely used to enhance neural differentiation during ES *in vitro* differentiation.^{25,26}

Induction of endoderm-derived cells during ES *in vitro* differentiation and particularly, insulin-secreting cells have been more difficult to obtain. Soria *et al.* applied genetic selection using the human insulin promoter fused to an antibiotic resistance gene to select cells which initiate synthesis of insulin.¹⁶ Lumelsky *et al.* demonstrated the formation of insulin producing pancreatic-like islet structures from ES cells based on methods known to promote the generation of neural cell types from ES cells.²⁷ However the use of nestin as a dependable marker for evaluating pancreatic development is somewhat controversial,²⁸ and further studies are required to conclusively show that nestin-derived cells can or cannot generate cells of the pancreatic lineage. Blyszczuk and coworkers have recently described a multistep process for differentiation of insulin-secreting cells without the selection and induction of nestin-positive cells.²⁹

A recent report demonstrates the production of another cell type from the endodermal germ layer. Ali *et al.* have demonstrated that ES cells can differentiate into Type II alveolar epithelial cells, which are needed for proper gas exchange in the lungs.³⁰ These preliminary results are very promising and may provide a model system to evaluate the possibility of using ES-derived cells to regenerate damaged lung tissue.

Angiogenesis is the formation of new blood vessels by the sprouting, splitting, and remodeling of existing vessels, and is of importance in many areas of research, including the studies of solid tumor growth and metastasis in cancer biology, wound healing, inflammation-related diseases, and cardiology. An ES-based system has been established for the study of angiogenesis.³¹ Feraud *et al.* plated undifferentiated mouse ES cells in semi-solid medium containing angiogenic cytokines. Three days later, the EBs have developed branching sprouts from their centers, indicative of sprouting angiogenesis.

Promotion of tissue-specific *in vitro* differentiation has been achieved by co-culture of ES cells on different stromal cell lines. Numerous studies have used the OP9 stromal cell line for the generation of lymphoid cells (B and NK cells).³²⁻³⁵ The OP9 cell line is derived from the calvaria of op/op mice that have a spontaneous mutation in the macrophage colony-stimulating factor (M-CSF) gene. This is significant because M-CSF promotes macrophage proliferation and inhibits ES cell differentiation to lymphoid cells *in vitro*.³² Kawasaki and colleagues have also identified a stromal cell line, Pa6, which expresses neural inducing activity (termed SDIA for Stromal-cell Derived Inducing Activity). Co-culture of Pa6 and ES cells results in efficient neural differentiation and derivation of a high proportion (30%) of transplantable, dopamine-producing neurons.³⁶

FUNCTIONAL STUDIES OF TRANSPLANTED MOUSE ES AND ES-DERIVED CELLS

Many studies show that transplanted ES and ES-derived cells are able to survive, integrate and, to some measurable extent, function within the host tissue. Bjorklund *et al.* transplanted small numbers of cells from disrupted day 4 EBs into a rat model of Parkinson's Disease, and obtained dopaminergic (DA) neurons which can reverse some symptoms of this disease.³⁷ However, although 56% of the animals showed surviving grafts containing DA neurons, 20% had lethal tumor-like masses called teratomas.³⁷ Another study also demonstrated "proof of concept" that ES cells can be directed *in vivo* to produce hepatocytes in response to endogenous signaling in adult animals.³⁷

Before ES cells can be used therapeutically, methods must be developed to avoid formation of teratomas *in vivo*. The use of committed, terminally differentiated cells for transplantation, rather than ES cells themselves may reduce the incidence of tumors. For example, Yin *et al.* used a gene targeting approach to integrate the green fluorescent protein (GFP) gene into the locus of a fetal liver-specific protein, alpha-fetoprotein, in undifferentiated ES cells.³⁹ Following *in vitro* differentiation, GFP-positive cells were isolated by FACS and injected into haptoglobin-deficient mice. The authors detected low levels of haptoglobin suggesting that ES-derived cells are able to mediate liver specific functions *in vivo*. More importantly, in over 100 mice transplanted with the ES-derived differentiated hepatocytes, teratomas were not observed.³⁹ Other investigators using ES-derived oligodendrocytes^{21,26,40} and DA neurons^{36,41} report no evidence of teratomas, although long-term studies are needed to confirm these results. McKay and co-workers enhanced the generation of ES-derived DA neurons using growth factor manipulation and expression of the transcription factor nuclear receptor related-1 (Nurr1), which has a role in differentiation of midbrain precursors to DA neurons. Transplanted neural cells were shown to integrate and remain functional for up to 8 weeks, with no evidence of teratoma formation.⁴¹

Another important consideration for evaluating the efficacy of ES-derived cells is determination of the functional durability of the transplanted cells. One group found that a large proportion of mice receiving ES-derived insulin-producing cells developed hyperglycemia 12 weeks later, implying that the implanted cells had a limited *in vivo* life span.¹⁶ Interestingly, Kyba *et al.* described an overexpression strategy with the transcription factor HOXB4, that has been implicated in self-renewal of definitive hematopoietic stem cells, combined with co-culture on the OP9 stromal layer to generate ES-derived hematopoietic stem cells. These cells were able to engraft lethally irradiated adults and contribute to long-term, multi-lineage hematopoiesis in primary and secondary recipients demonstrating long-term functionality of these cells.¹⁴ This study is intriguing as it illustrates that ES-derived somatic stem cells, rather than terminally differentiated cells, may provide a suitable cell source for transplantation.

In summary, significant progress has been made in controlling the *in vitro* differentiation of ES cells into a wide variety of cell types. This provides a powerful model system to investigate molecular and cellular processes involved in lineage-specification and embryogenesis. Protocols based on the identification of genes and cell surface markers for somatic stem cells or lineage precursors may further improve the efficacy and safety of these strategies to provide functional cells for regenerative medicine.

TABLE 1

Differentiated cell types derived from mouse embryonic stem cells *in vitro*

Embryonic Germ Layer	Cell Type	Reference(s)
Ectoderm (external layer)	neurons	19, 25
	dopaminergic neurons	20, 37, 41
	motor neurons	42
	oligodendrocytes	21, 26, 40, 43
	astrocytes	44
	epithelial cells	45
Mesoderm (middle layer)	adipocytes	46
	cardiomyocytes	47-50
	chondrocytes	51
	hematopoietic cells	23, 52 - 54
	hematopoietic stem cells	14
	dendritic cells	24
	mast cells	22
	B cells	32, 33, 35
	natural killer cells	33, 34
	lymphoid precursors	55
	endothelial cells	15, 18, 31
Endoderm (internal layer)	osteoblasts	56
	striated muscle cells	57
	smooth muscle cells	18
	macrophages	58
	pancreatic-like islets	27
	insulin-producing cells	16, 29
	alveolar epithelium	30
	hepatocytes	59

Embryonic stem cells differentiate *in vitro* into each of the three embryonic germ layers (ectoderm, mesoderm, endoderm) in a manner that recapitulates *in vivo* embryogenesis. These cells can subsequently be induced to differentiate into the majority of cell types of the adult mouse.

REFERENCES:

- Evans MJ, Kaufman MH: Establishment in culture of pluripotential cells from mouse embryos. *Nature* 292:154-156, 1981
- Martin GR: Isolation of a pluripotent cell line from early mouse embryos cultured in medium conditioned by teratocarcinoma stem cells. *Proc Natl Acad Sci USA* 78: 7634-7638, 1981
- Bradley A, Evans M, Kaufman MH, Robertson E: Formation of germ-line chimaeras from embryo-derived teratocarcinoma cell lines. *Nature* 309: 255-256, 1984
- Williams RL, Hilton DJ, Pease S, Willson TA, Stewart CL, Gearing DP, Wagner EF, Metcalf D, Nicola NA, Gough NM: Myeloid leukaemia inhibitory factor maintains the developmental potential of embryonic stem cells. *Nature* 336: 684-687, 1988
- Chambers I, Smith A: Self-renewal of teratocarcinoma and embryonic stem cells. *Oncogene* 23: 7150-60, 2004
- Ying QL, Nichols J, Chambers I, Smith A: BMP induction of Id proteins suppresses differentiation and sustains embryonic stem cell self-renewal in collaboration with STAT3. *Cell* 115:281-92, 2003
- Feng Z, Srivastava AS, Mishra R, Carrier E: A regulatory role of Wnt signaling pathway in the hematopoietic differentiation of murine embryonic stem cells. *Biochem Biophys Res Commun* 324:1333-9, 2004
- Sato N, Meijer L, Skaltsounis L, Greengard P, Brivanlou AH: Maintenance of pluripotency in human and mouse embryonic stem cells through activation of Wnt signaling by a pharmacological GSK-3-specific inhibitor. *Nat Med* 10: 55-63, 2004
- Nichols J, Zevnik B, Anastassiadis K, Niwa H, Klewe-Nebenius D, Chambers I, Scholer H, Smith A: Formation of pluripotent stem cells in the mammalian embryo depends on the POU transcription factor Oct4. *Cell* 95:379-91, 1998

- Mitsui K, Tokuzawa Y, Itoh H, Segawa K, Murakami M, Takahashi K, Maruyama M, Maeda M, Yamanaka S: The homeoprotein Nanog is required for maintenance of pluripotency in mouse epiblast and ES cells. *Cell* 113:631-42, 2003
- Chambers I, Colby D, Robertson M, Nichols J, Lee S, Tweedie S, Smith A: Functional expression cloning of Nanog, a pluripotency sustaining factor in embryonic stem cells. *Cell* 113:643-55, 2003
- Rathjen J, Rathjen PD: Mouse ES cells: experimental exploitation of pluripotent differentiation potential. *Curr Opin Genet Dev* 11: 587-594, 2000
- Weiss MJ, Orkin SH: *In vitro* differentiation of mouse embryonic stem cells. New approaches to old problems. *J Clin Invest* 97: 591-595, 1996
- Kyba M, Perlingeiro RC, Daley GQ: HoxB4 confers definitive lymphoid-myeloid engraftment potential on embryonic stem cell and yolk sac hematopoietic progenitors. *Cell* 109: 29-37, 2002
- Marchetti S, Gimond C, Iljin K, Bourcier C, Alitalo K, Pouyssegur J, Pages G: Endothelial cells genetically selected from differentiating mouse embryonic stem cells incorporate at sites of neovascularization *in vivo*. *J Cell Sci* 115: 2075-2085, 2002
- Soria B, Roche E, Berna G, Leon-Quinto T, Reig JA, Martin F: Insulin-secreting cells derived from embryonic stem cells normalize glycemia in streptozotocin-induced diabetic mice. *Diabetes* 49: 157-162, 2000
- Klug MG, Soonpaa MH, Koh GY, Field LJ: Genetically selected cardiomyocytes from differentiating embryonic stem cells form stable intracardiac grafts. *J Clin Invest* 98: 216-224, 1996
- Yamashita J, Itoh H, Hirashima M, Ogawa M, Nishikawa S, Yurugi T, Naito M, Nakao K: Flk1-positive cells derived from embryonic stem cells serve as vascular progenitors. *Nature* 408: 92-96, 2000
- Okabe S, Forsberg-Nilsson K, Spiro AC, Segal M, McKay RD: Development of neuronal precursor cells and functional postmitotic neurons from embryonic stem cells *in vitro*. *Mech Dev* 59: 89-102, 1996
- Lee SH, Lumelsky N, Studer L, Auerbach JM, McKay RD: Efficient generation of midbrain and hindbrain neurons from mouse embryonic stem cells. *Nat Biotechnol* 18: 675-679, 2000
- Brustle O, Jones KN, Learish RD, Karram K, Choudhary K, Wiestler OD, Duncan ID, McKay RD: Embryonic stem cell-derived glial precursors: a source of myelinating transplants. *Science* 285: 754-756, 1999
- Tsai M, Wedemeyer J, Ganiatsas S, Tam SY, Zon LI, Galli SJ: *In vivo* immunological function of mast cells derived from embryonic stem cells: an approach for the rapid analysis of even embryonic lethal mutations in adult mice *in vivo*. *Proc Natl Acad Sci USA* 97: 9186-9190, 2000
- Wiles MV, Keller G: Multiple hematopoietic lineages develop from embryonic stem (ES) cells in culture. *Development* 111: 259-267, 1991
- Fairchild PJ, Brook FA, Gardner RL, Graca L, Strong V, Tone Y, Tone M, Nolan KF, Waldmann H: Directed differentiation of dendritic cells from mouse embryonic stem cells. *Curr Biol* 10: 1515-1518, 2000
- Bain G, Kitchens D, Yao M, Huettner JE, Gottlieb DI: Embryonic stem cells express neuronal properties *in vitro*. *Dev Biol* 168: 342-357, 1995
- McDonald JW, Liu XZ, Qu Y, Liu S, Mickey SK, Turetsky D, Gottlieb DI, Choi DW: Transplanted embryonic stem cells survive, differentiate and promote recovery in injured rat spinal cord. *Nat Med* 5: 1410-1412, 1999
- Lumelsky N, Blondel O, Laeng P, Velasco I, Ravin R, McKay R: Differentiation of embryonic stem cells to insulin-secreting structures similar to pancreatic islets. *Science* 292: 1389-1394, 2001
- Selander L, Edlund H: Nestin is expressed in mesenchymal and not epithelial cells of the developing mouse pancreas. *Mech Dev* 113: 189-192, 2002
- Blyszczuk P, Asbrand C, Rozzo A, Kania G, St-Onge L, Rupnik M, Wobus AM: Embryonic stem cells differentiate into insulin-producing cells without selection of nestin-expressing cells. *Int J Dev Biol* 48:1095-104, 2004
- Ali NN, Edgar AJ, Samadikuchaksaraei A, Timson CM, Romanska HM, Polak JM, Bishop AE: Derivation of Type II alveolar epithelial cells from mouse embryonic stem cells. *Tissue Eng* 8: 541-550, 2002
- Feraud O, Cao Y, Vittet D: Embryonic stem cell-derived embryoid bodies development in collagen gels recapitulates sprouting angiogenesis. *Lab Invest* 81:1669-81, 2001

32. Nakano T, Kodama H, Honjo T: Generation of lymphohematopoietic cells from embryonic stem cells in culture. *Science* 265: 1098-1101, 1994
33. Nakayama N, Fang I, Elliott G: Natural killer and B-lymphoid potential in CD34+ cells derived from embryonic stem cells differentiated in the presence of vascular endothelial growth factor. *Blood* 91: 2283-2295, 1998
34. Lian RH, Maeda M, Lohwasser S, Delcommenne M, Nakano T, Vance RE, Raulat DH, Takei F: Orderly and nonstochastic acquisition of CD94/NKG2 receptors by developing NK cells derived from embryonic stem cells *in vitro*. *J Immunol* 168: 4980-4987, 2002
35. Cho SK, Webber TD, Carlyle JR, Nakano T, Lewis SM, Zuniga-Pflucker JC: Functional characterization of B lymphocytes generated *in vitro* from embryonic stem cells. *Proc Natl Acad Sci USA*, 96: 9797-9802, 1999
36. Kawasaki H, Mizuseki K, Nishikawa S, Kaneko S, Kuwana Y, Nakanishi S, Nishikawa SI, Sasai Y: Induction of midbrain dopaminergic neurons from ES cells by stromal cell-derived inducing activity. *Neuron* 28:31-40, 2000
37. Bjorklund LM, Sanchez-Pernaute R, Chung S, Andersson T, Chen IY, McNaught KS, Brownell AL, Jenkins BG, Wahlestedt C, Kim KS, Isacson O: Embryonic stem cells develop into functional dopaminergic neurons after transplantation in a Parkinson rat model. *Proc Natl Acad Sci USA* 99: 2344-2349, 2002
38. Choi D, Oh HJ, Chang UJ, Koo SK, Jiang JX, Hwang SY, Lee JD, Yeoh GC, Shin HS, Lee JS, Oh B: *In vivo* differentiation of mouse embryonic stem cells into hepatocytes. *Cell Transplant* 11: 359-368, 2002
39. Yin Y, Lim YK, Salto-Tellez M, Ng SC, Lin CS, Lim SK: AFP(+), ESC-derived cells engraft and differentiate into hepatocytes *in vivo*. *Stem Cells* 20: 338-346, 2002
40. Liu S, Qu Y, Stewart TJ, Howard MJ, Chakraborty S, Holekamp TF, McDonald JW: Embryonic stem cells differentiate into oligodendrocytes and myelinate in culture and after spinal cord transplantation. *Proc Natl Acad Sci USA* 97: 6126-6131, 2000
41. Kim JH, Auerbach JM, Rodriguez-Gomez JA, Velasco I, Gavin D, Lumelsky N, Lee SH, Nguyen J, Sanchez-Pernaute R, Bankiewicz K, McKay R: Dopamine neurons derived from embryonic stem cells function in an animal model of Parkinson's disease. *Nature* 418: 50-56, 2002
42. Wichterle H, Lieberam I, Porter J, Jessell T: Directed differentiation of embryonic stem cells into motor neurons. *Cell* 110: 385, 2002
43. Billon N, Jolicoeur C, Ying QL, Smith A, Raff M: Normal timing of oligodendrocyte development from genetically engineered, lineage-selectable mouse ES cells. *J Cell Sci* 115: 3657-3665, 2002
44. Fraichard A, Chassande O, Bilbaut G, Dehay C, Savatier P, Samarut J: *In vitro* differentiation of embryonic stem cells into glial cells and functional neurons. *J Cell Sci* 108: 3181-3188, 1995
45. Bagutti C, Wobus AM, Fassler R, Watt FM: Differentiation of embryonal stem cells into keratinocytes: comparison of wild-type and beta 1 integrin-deficient cells. *Dev Biol*: 179: 184-196, 1996
46. Dani C, Smith AG, Dessolin S, Leroy P, Staccini L, Villageois P, Darimont C, Ailhaud G: Differentiation of embryonic stem cells into adipocytes *in vitro*. *J Cell Sci* 110: 1279-1285, 1997
47. Boheler KR, Czyz J, Tweedie D, Yang HT, Anisimov SV, Wobus AM: Differentiation of pluripotent embryonic stem cells into cardiomyocytes. *Circ Res* 91:189-201, 2002
48. Wobus AM, Wallukat G, Hescheler J: Pluripotent mouse embryonic stem cells are able to differentiate into cardiomyocytes expressing chronotropic responses to adrenergic and cholinergic agents and Ca²⁺ channel blockers. *Differentiation* 48: 173-182, 1991
49. Wobus AM, Kaomei G, Shan J, Wellner MC, Rohwedel J, Ji G, Fleischmann B, Katus HA, Hescheler J, Franz WM: Retinoic acid accelerates embryonic stem cell-derived cardiac differentiation and enhances development of ventricular cardiomyocytes. *J Mol Cell Cardiol* 29: 1525-1539, 1997
50. Maltsev VA, Rohwedel J, Hescheler J, Wobus AM: Embryonic stem cells differentiate *in vitro* into cardiomyocytes representing sinusnodal, atrial and ventricular cell types. *Mech Dev* 44: 41-50, 1993
51. Kramer J, Hegert C, Guan K, Wobus AM, Muller PK, Rohwedel J: Embryonic stem cell-derived chondrogenic differentiation *in vitro*: activation by BMP-2 and BMP-4. *Mech Dev* 92: 193-205, 2000
52. Carotta S, Pilat S, Mairhofer A, Schmidt U, Dolznig H, Steinlein P, Beug H: Directed differentiation and mass cultivation of pure erythroid progenitors from mouse embryonic stem cells. *Blood* 104:1873-80, 2004
53. Nakano T, Kodama H, Honjo T: *In vitro* development of primitive and definitive erythrocytes from different precursors. *Science* 272: 722-724, 1996
54. Keller G, Kennedy M, Papayannopoulou T, Wiles MV: Hematopoietic commitment during embryonic stem cell differentiation in culture. *Mol Cell Biol* 13: 473-486, 1993
55. Gutierrez-Ramos JC, Palacios R: *In vitro* differentiation of embryonic stem cells into lymphocyte precursors able to generate T and B lymphocytes *in vivo*. *Proc Natl Acad Sci USA* 89: 9171-9175, 1992
56. Buttery LD, Bourne S, Xynos JD, Wood H, Hughes FJ, Hughes SP, Episkopou V, Polak JM: Differentiation of osteoblasts and *in vitro* bone formation from mouse embryonic stem cells. *Tissue Eng* 7: 89-99, 2001
57. Rohwedel J, Maltsev V, Bober E, Arnold HH, Hescheler J, Wobus AM: Muscle cell differentiation of embryonic stem cells reflects myogenesis *in vivo*: developmentally regulated expression of myogenic determination genes and functional expression of ionic currents. *Dev Biol* 164: 87-101, 1994
58. Lindmark H, Rosengren B, Hurt-Camejo E, Bruder CE: Gene expression profiling shows that macrophages derived from mouse embryonic stem cells is an improved *in vitro* model for studies of vascular disease. *Exp Cell Res* 300:335-44, 2004
59. Jones EA, Tosh D, Wilson DI, Lindsay S, Forrester LM: Hepatic differentiation of mouse embryonic stem cells. *Exp Cell Res* 272: 15-22, 2002