

Stem Cells and Differentiation

Rebecca Quelch¹ and Stefan Przyborski^{1,2}

¹Department of Biosciences, Durham University, South Road, Durham, UK.

²ReproCELL Europe Ltd, NET Park Incubator, Sedgfield, UK.

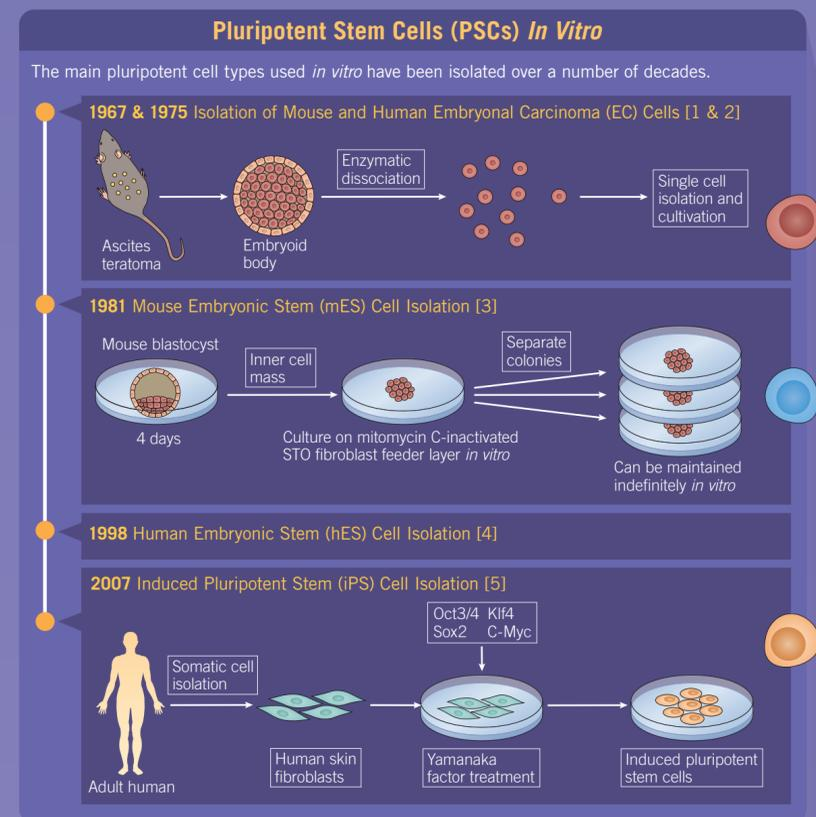
Email: s.a.przyborski@durham.ac.uk and r.h.quelch@durham.ac.uk

TOCRIS
a biotechne brand

Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals

www.tocris.com

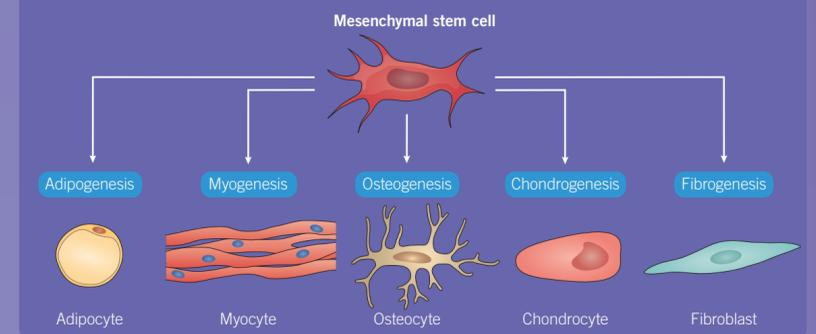
Stem cells are defined by their ability to self-renew and propensity to differentiate into functional cell types. Pluripotent Stem Cells (PSCs) differentiate into cells of all three germ layers (endoderm, ectoderm and mesoderm), whereas multipotent stem cells are more limited in their differentiation potential. The unique abilities of these cell types make them attractive tools for a wide range of applications, from regenerative medicine to drug toxicity screening. There are a large number of protocols for the maintenance of this pluripotent state, as well as for the subsequent directed differentiation of these cells *in vitro* to form specialized cell types.



Differentiation

Adult Stem Cells (ASCs)

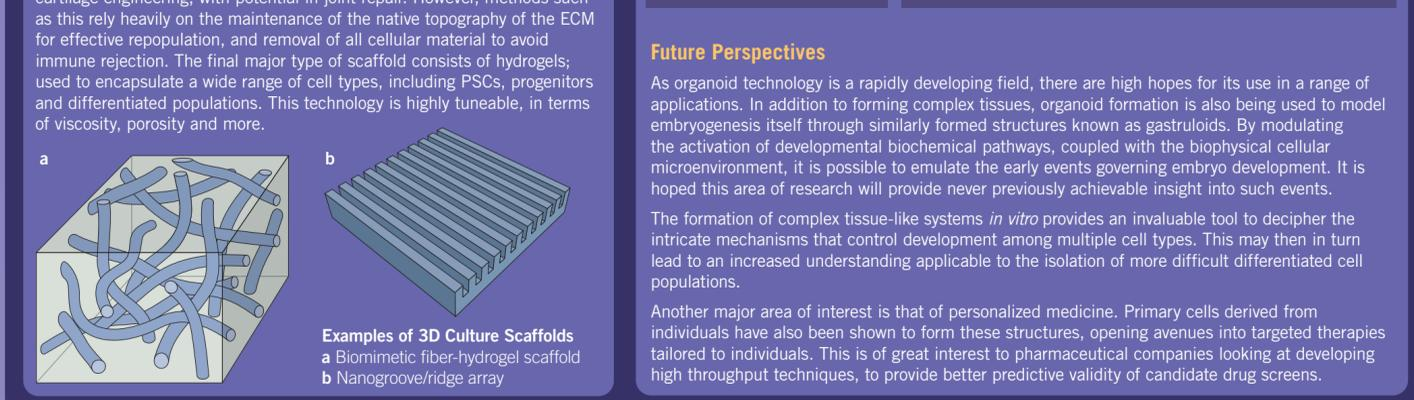
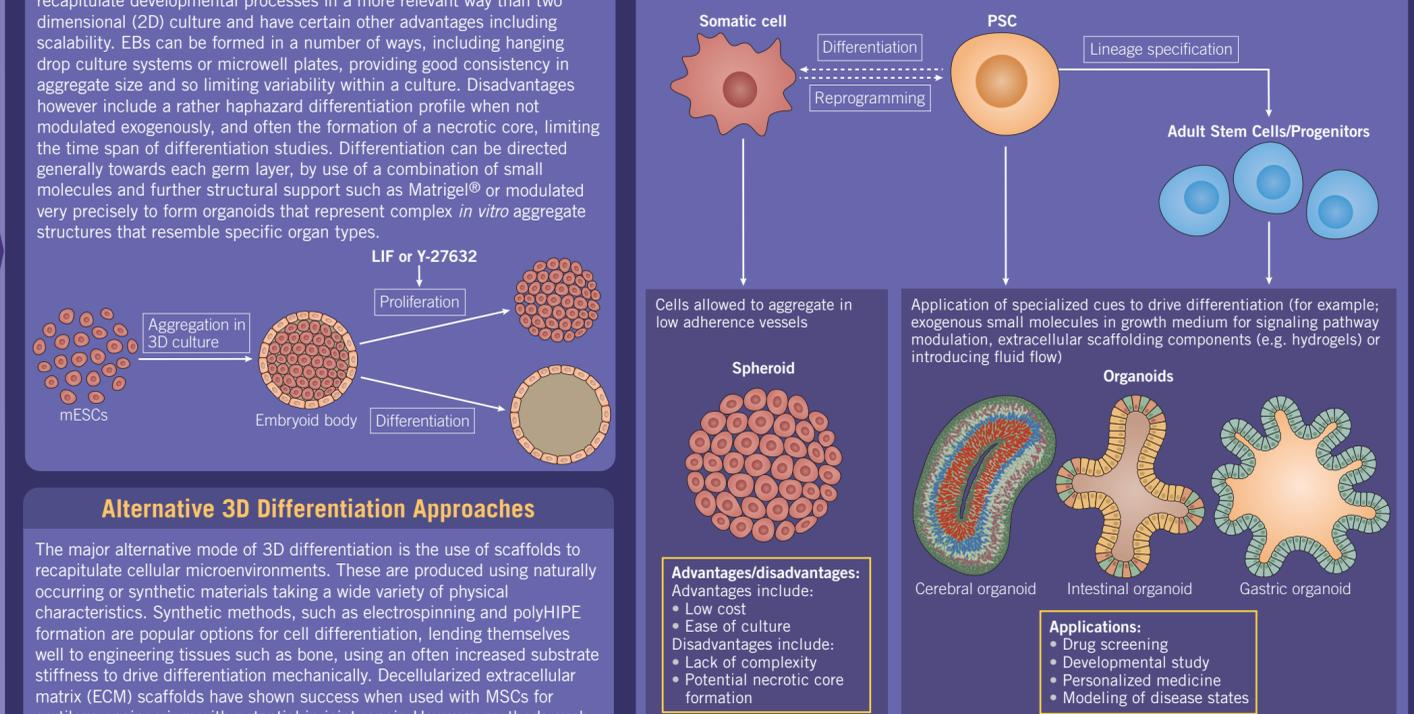
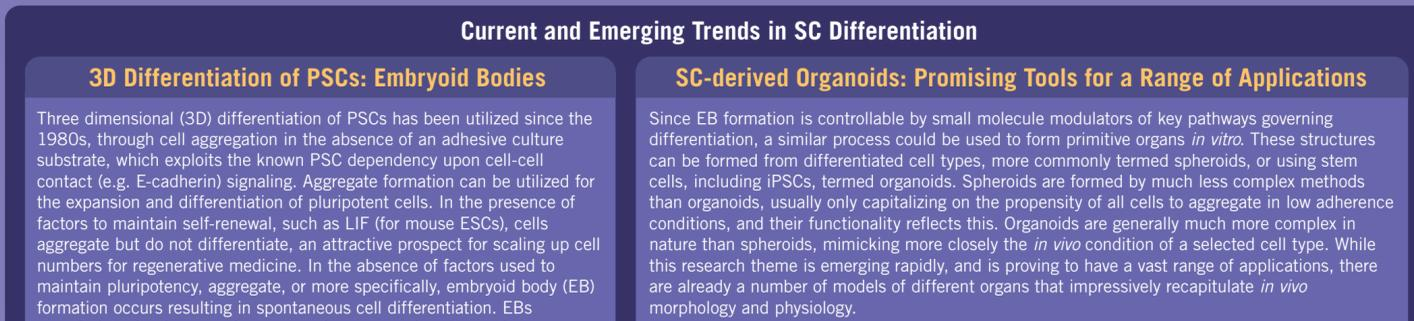
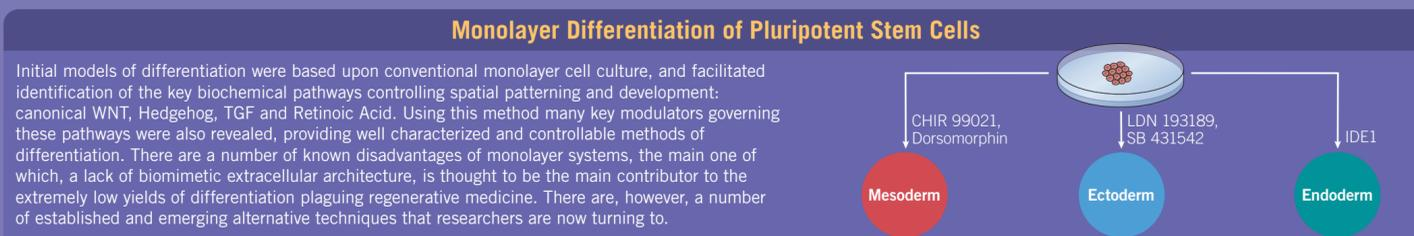
It is thought that most organs contain resident adult (or somatic) stem cell populations, i.e. groups of cells that are able to self-renew and differentiate into a limited number of cell types to repopulate an organ for maintenance of functionality throughout normal cell turnover or upon challenge/injury. Mesenchymal stem cells (MSCs) are a very well characterized somatic stem cell type, due to their ease of isolation from connective tissue and relatively large capacity for differentiation. Recently, the receptor LGR5 has gained provenance as a marker for a range of more specific organ adult cell types, such as in the intestine and liver. Such markers are important in the development of organoid technology.



Maintenance of Stem Cell Phenotype *In Vitro*

The maintenance of stem cell self-renewal and an undifferentiated status is often overlooked. However stem cells occupy a continuum of developmental states (e.g. "naïve" vs. "primed" pluripotency) that can, in some cases, be defined by the culture environment. Methods of maintaining pluripotency were initially focused around growth on "feeder" layers of fibroblasts. The spotlight subsequently shifted to specialized growth medium formulations and the use of qualified serum and small molecules. LIF for mESCs and bFGF for hESCs, are most commonly used to maintain pluripotency, but there are a range of other compounds, including Y-27632, A83-01 or SB 431542, which can be used to maintain cell phenotype.

More recently focus has shifted further towards the composition of the growth substrate itself; research concerning substrate coatings (including tailored integrins) and substrate topography (hydrogels and other scaffolds) is ongoing for both the maintenance and differentiation of PSCs.



Products available from Tocris

Reprogramming
(±)-Bay K 8644, BIX 01294, CHIR 99021, DBZ, 3-Deazaneplanocin A, Kenpaullone, L-Ascorbic acid, PD 0325901, RepSox, (S)-(-)-Dimethindene, SMER 28, Thiazovivin, Trichostatin A, Valproic acid, sodium salt
Differentiation
Neural Stem Cells
DAPT, Dibutyl-γ-cAMP, Forskolin, Fluoxetine, IBMX, LDN 193189, Metformin, SAG dihydrochloride, SU 5402
Cardiomyocyte
Dorsomorphin, 1-EBIO, ISX 9, IWP 4, Wnt-C59
Mesenchymal Stem Cells
AICAR, Dexamethasone, Purmorphamine, SP 600125, Zebularine
Other
IDE1
Proliferation and Cell Viability
A 769662, A 83-01, BIO, CH 223191, Go 6983, IWP 2, LY 294002, Pluripotin, Prostaglandin E2, PD 173074, PD 98059, SB 202190, SB 203580, SB 216763, SB 431542, Troglitazone, U0126, Y-27632
Stem Cell Signaling
AMD 3100, Cyclopamine, IWP 2, Mitomycin C, XAV 939

Stem cell products available from other Bio-Techne brands

Cell Culture and Differentiation
Cultrex® Stem Cell-Qualified BME – ECM scaffolding for stem cells Cultrex® BME, Type 2 – ECM scaffolding for organoid growth N21-MAX – Media supplement to support stem cell and organoid health N-2 MAX – Media supplement to support stem cell and organoid health
Verifying Pluripotency and Stemness
Human Pluripotent Stem Cell Functional Identification Kit GloLIVE™ Human Pluripotent Stem Cell Live Cell Imaging Kit
Growth Factors and Proteins
Recombinant Human/Mouse/Rat/Canine/Equine BDNF Protein rhFGF basic (146 aa) Protein rhGDNF Protein rhFGF basic (146 aa) Protein rHLIF Protein

Abbreviations

EC, Embryonal Carcinoma; ECM, Extracellular Matrix; ES, Embryonic Stem; iPS, Induced Pluripotent Stem; AS, Adult Stem; EB, Embryoid Body; MSC, Mesenchymal Stem Cell; 2D, two dimensional; 3D, three dimensional

References

- Finch, BW and Ephrussi, B (1967) *Proc. Natl. Acad. Sci. U.S.A.* **57** 615
- Fogh, J and Tepper, G (1975) New human tumor cell lines. In *Human Tumor Cells In Vitro*. J. Fogh, (ed.) 115–159. Plenum, New York.
- Evans, HJ and Kaufman, MH (1981) *Nature* **292** 154
- Thomson, J et al (1998) *Science* **282** 1145
- Takahashi, K et al (2007) *Cell* **131** 861

Recommended Reviews

Dutta, D et al (2017) *Trends Mol. Med.* **23** 393
Clarke, K et al (2015) *Tocris scientific review series: Using small molecules to control stem cell growth and differentiation.*

For copies of this poster, please visit tocris.com
© 2018 Tocris Cookson, Ltd.
Tocris is a Bio-Techne brand