

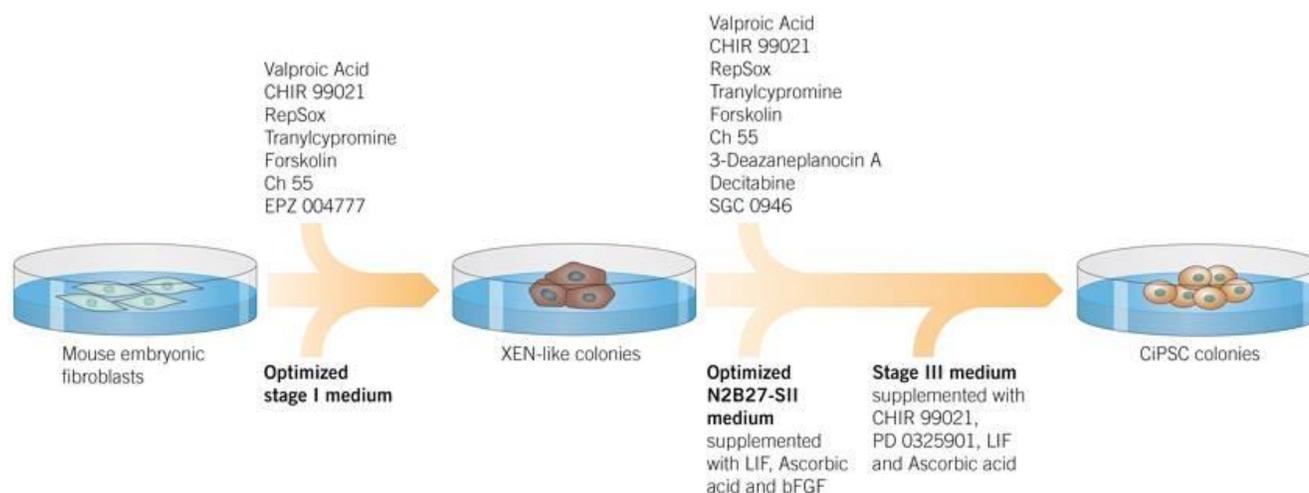
Highly Efficient Generation of CiPSCs from MEFs

This is intended as a guide only; for full experimental details please read the reference provided.

In Brief

Zhao et al. describe a protocol for the highly efficient reprogramming of mouse embryonic fibroblasts (MEFs) to chemically-induced pluripotent stem cells (CiPSCs) in 16-20 days using a small molecule cocktail.

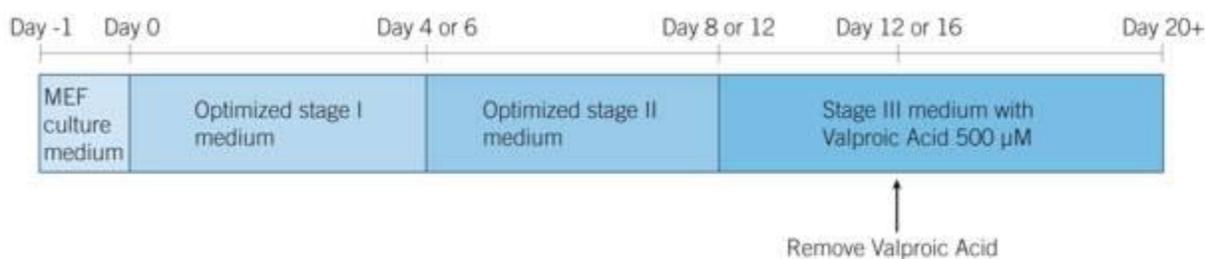
Briefly, MEFs were grown in MEF culture medium. On day 0 the medium was changed to optimized stage I medium. On day 4-6 small extraembryonic endoderm (XEN)-like colonies appeared and stage I medium was exchanged for optimized N2B27-SII medium. Cells were cultured for a further 4-6 days, then on day 8 or 12 N2B27-SII medium was exchanged for stage III medium containing 500 µM Valproic Acid (VPA). VPA was removed on day 12 or 16, and cells were cultured for a further 4-10 days in stage III medium, when CiPSC colonies emerged.



Cocktails

Optimized Stage I Medium (FBS/KSR-based medium)		Optimized N2B27-SII Medium (N2B27-based medium)		Stage III Medium (N2B27-based medium)	
Valproic acid (Cat.No. 2815)	100 µM	Valproic acid (Cat.No. 2815)	1 mM	CHIR 99021 (Cat.No. 4423)	3 µM
CHIR 99021 (Cat.No. 4423)	40 µM	CHIR 99021 (Cat.No. 4423)	10 µM	PD 0325901 (Cat.No. 4192)	1 µM

Optimized Stage I Medium (FBS/KSR-based medium)		Optimized N2B27-SII Medium (N2B27-based medium)		Stage III Medium (N2B27-based medium)	
RepSox (Cat.No. 3742)	10 µM	RepSox (Cat.No. 3742)	10 µM	LIF	10 ng/ml
Tranylcypromine (Cat.No. 3852)	5 µM	Tranylcypromine (Cat.No. 3852)	5 µM	L-Ascorbic Acid (Cat.No. 4055)	50 µg/ml
Forskolin (Cat.No. 1099)	10 µM	Forskolin (Cat.No. 1099)	10 µM		
Ch 55 (Cat.No. 2020)	1 µM	Ch 55 (Cat.No. 2020)	1 µM		
EPZ 004777 (Cat.No. 5567)	5 µM	3-Deazaneplanocin A (Cat.No. 4703)	0.05 µM		
		Decitabine (Cat.No. 2624)	0.5 µM		
		SGC 0946 (Cat.No. 4541)	5 µM		
		L-Ascorbic Acid (Cat.No. 4055)	50 µg/ml		
		LIF	10 ng/ml		
		bFGF	25 ng/ml		



Reference

Zhao *et al.* (2018) Single-cell RNA-seq reveals dynamic early embryonic-like programs during chemical reprogramming. *Cell Stem Cell*. **23** 31 PMID: [29937202](#)