

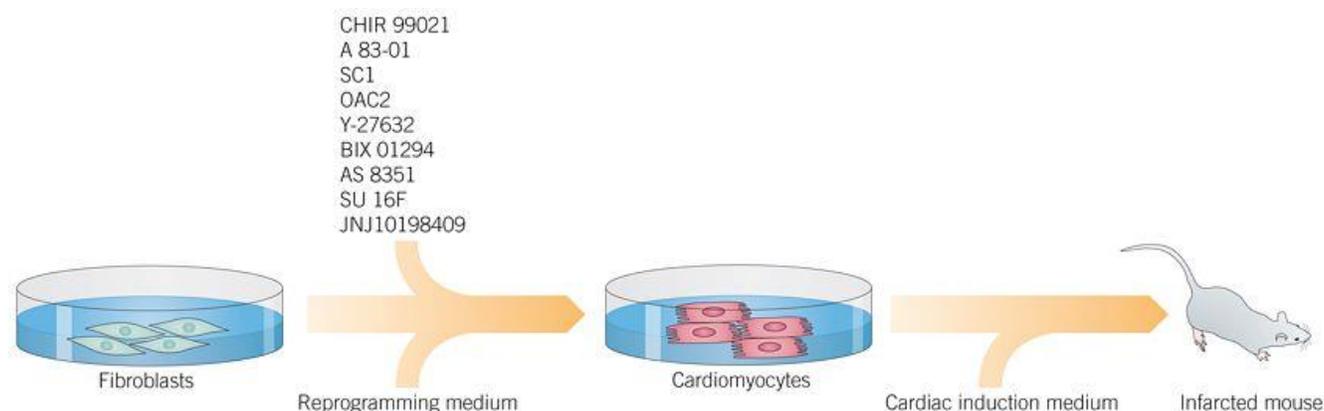
## Converting Fibroblasts into Cardiomyocytes (9C Cocktail)

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

### In Brief

Cao *et al.* converted human fibroblasts into functional cardiomyocytes using a 9C cocktail. The 9C-treated fibroblasts were transplanted into an infarcted mouse heart and converted to cardiomyocyte-like cells.

Human foreskin fibroblasts (HFF) were treated with 9C for 6 days and CIM for 5 days then transplanted into mice with infarcted hearts. When transplanted, 9C-treated fibroblasts were efficiently converted into cardiomyocyte-like cells.



### Cocktails

9C Medium		Cardiac Induction Medium	
<a href="#">CHIR 99021</a> (Cat.No. 4423)	10 µM	<a href="#">CHIR 99021</a> (Cat.No. 4423)	12 µM
<a href="#">A 83-01</a> (Cat.No. 2939)	1 µM	BMP-4	25 ng/ml
<a href="#">SC1</a> (Cat.No. 4433)	1 µM	Activin A	10 ng/ml
<a href="#">OAC-2</a> (Cat.No. 6066)	5 µM	VEGF	10 ng/ml
<a href="#">Y-27632</a> (Cat.No. 1254)	10 µM		

**9C Medium**

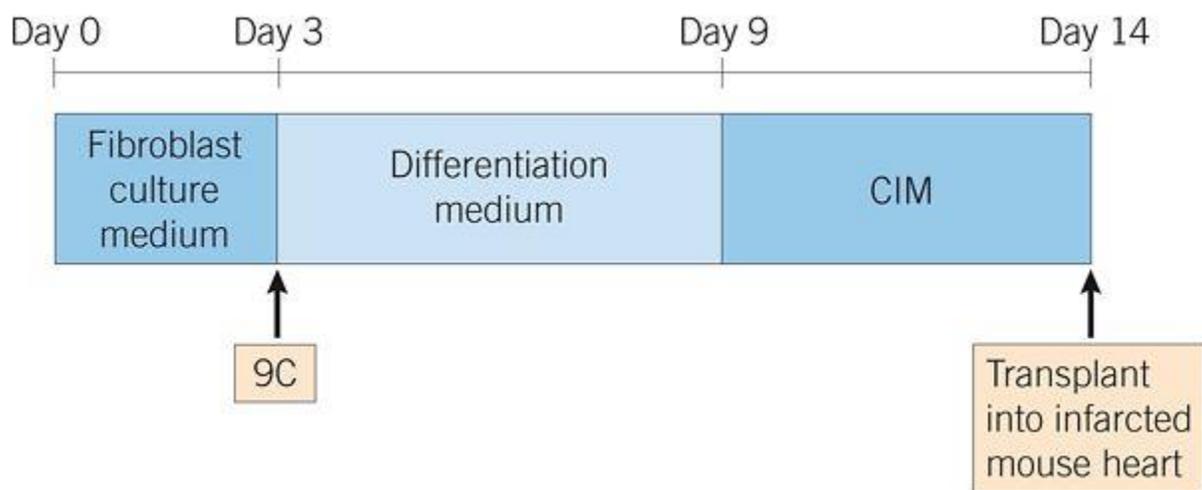
**Cardiac Induction Medium**

[BIX 01294](#) (Cat.No. 3364) 1 µM

[AS 8351](#) (Cat.No. 6044) 1 µM

[SU 16f](#) (Cat.No. 3304) 5 µM

JNJ 10198409 0.1 µM



**Reference**

**Cao et al.** (2016) Conversion of human fibroblasts into functional cardiomyocytes by small molecules. *Science* **352** 6290. PMID: [27127239](#)