

## Differentiation of Hematopoietic Progenitors from hPSCs

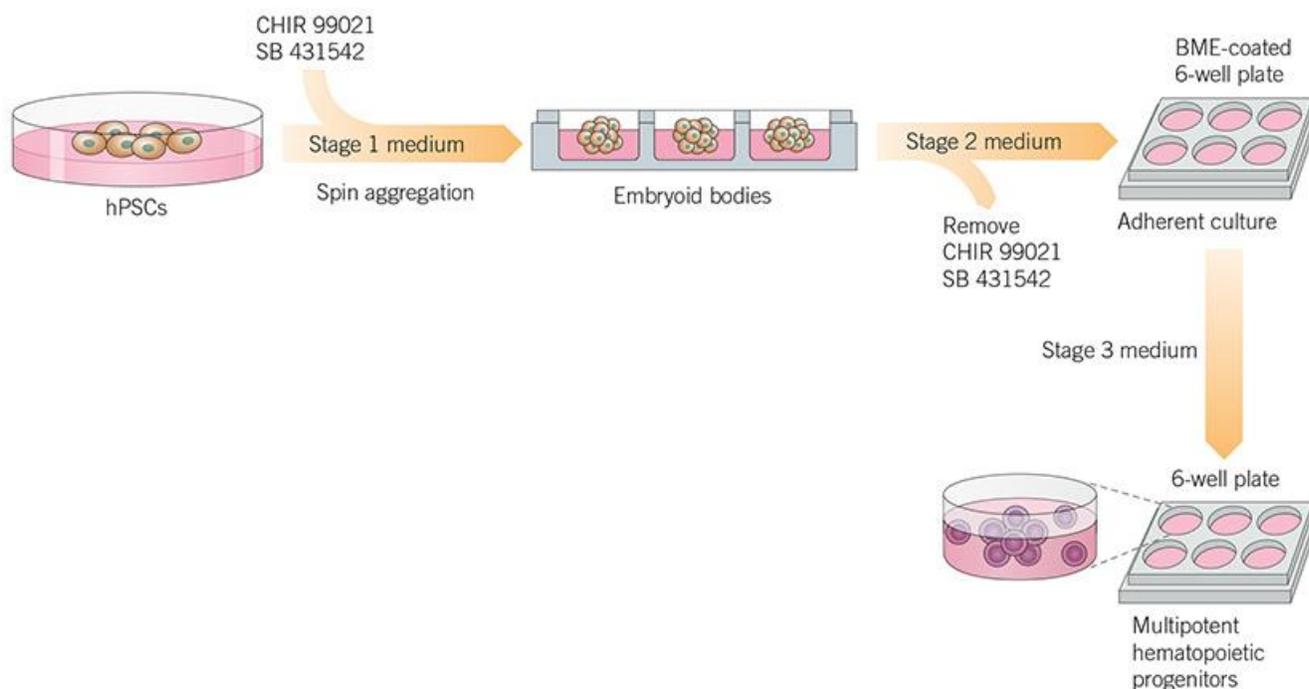
*This is intended as a guide only; for full experimental details please read the reference provided.*

### In Brief

Nafria *et al.* describe a 3-stage procedure for the [differentiation](#) of human pluripotent stem cells (hPSCs) to multipotent hematopoietic progenitors in 14+ days using a cocktail of small molecules and growth factors.

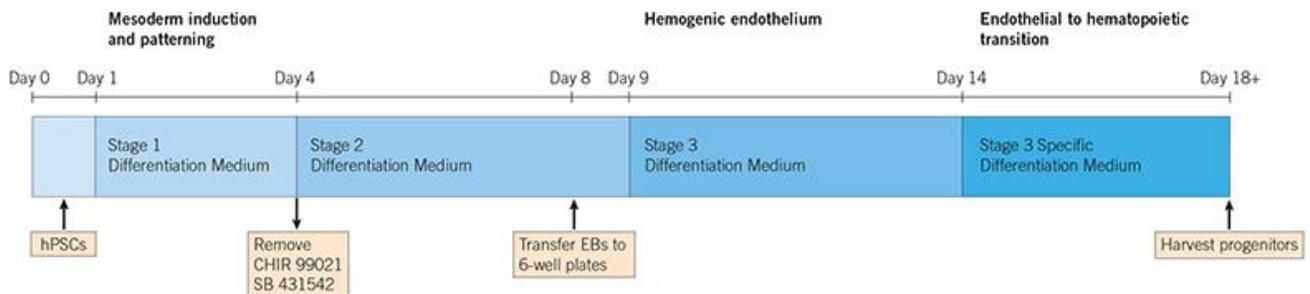
Briefly, hPSCs are co-cultured with feeder cells for 1 week, after which the hPSCs are harvested, dissociated and seeded in non-adherent round-bottom 96 well plates in stage 1 differentiation medium. On day 0 plates are centrifuged and cells aggregate forming embryoid bodies (EB; 1 per well). On day 1, [SB 431542](#) and [CHIR 99021](#) are added to the Stage 1 differentiation medium to direct the mesoderm toward the definitive hematopoietic program. On day 4 the medium is exchanged for Stage 2 medium without small molecule inhibitors. At day 8 of differentiation EBs are transferred to Basement Membrane Extract (BME, e.g. [Cultrex™](#)) coated 6-well plates containing stage 2 medium. The next day (Day 9) the medium is removed and replaced with stage 3 medium. Then around day 13-15 Stage 3 medium is replaced with Stage 3 Specific medium to promote proliferation and differentiation of hematopoietic lineage cells.

Vascular structures and hemogenic endothelium expressing the SOX17 marker are generated from day 8 to 14 and from day 14 onward, SOX17- CD34+ RUNX1C-/+ hematopoietic progenitors are generated within the SOX17+ endothelium displaying similarities to human aorta-gonad mesonephros (AGM) and adult hematopoietic stem cells.



## Cocktails

Stage 1 Differentiation Medium		Stage 2 Differentiation Medium		Stage 3 Differentiation Medium		Stage 3 Specific Medium	
STAPEL medium		STAPEL medium		STAPEL medium		STAPEL medium	
<a href="#">CHIR 99021</a>	0.5 µM	<a href="#">BMP-4</a>	20 ng/mL	<a href="#">BMP-4</a>	10 ng/mL	<a href="#">SCF</a>	100 ng/mL
Cat.No 4423		314-BP/CF		314-BP/CF		7466-SC/CF	
<a href="#">BMP-4</a>	20 ng/mL	<a href="#">VEGF</a>	50 ng/mL	<a href="#">SCF</a>	100 ng/mL	<a href="#">Flt3</a>	100 ng/mL
314-BP/CF		293-VE/CF		7466-SC/CF		308-FK/CF	
<a href="#">VEGF</a>	25 ng/mL	<a href="#">SCF</a>	50 ng/mL	<a href="#">Flt3</a>	100 ng/mL	<a href="#">TPO</a>	50 ng/mL
293-VE/CF		7466-SC/CF		308-FK/CF		288-TP/CF	
<a href="#">SCF</a>	25 ng/mL	<a href="#">IGF-2</a>	10 ng/mL	<a href="#">TPO</a>	50 ng/mL	<a href="#">IL-6</a>	25 ng/mL
7466-SC/CF		292-G2		288-TP/CF		7270-IL/CF	
<a href="#">Activin A</a>	7.5 ng/mL	<a href="#">FGF-2</a>	10 ng/mL	<a href="#">VEGF</a>	50 ng/mL	<a href="#">IL-3</a>	25 ng/mL
388-AC/CF		3718-FB		293-VE/CF		203-IL/CF	
<a href="#">FGF-2</a>	10 ng/mL			<a href="#">IL-6</a>	25 ng/mL	<a href="#">Penicillin/Streptomycin</a>	Optional
3718-FB				7270-IL/CF		B21210	
<b>From Day 1</b>				<a href="#">IL-3</a>	25 ng/mL		
				203-IL/CF			
<a href="#">CHIR 99021</a>	3 µM			<a href="#">IGF-2</a>	10 ng/mL		
Cat.No 4423				292-G2			
<a href="#">SB 431542</a>	3.8 µM			<a href="#">FGF-2</a>	10 ng/mL		
Cat.No 1614				3718-FB			
				<a href="#">Penicillin/Streptomycin</a>	Optional		
				B21210			



## Reference

Nafria *et al.* (2020) Protocol for the generation of definitive hematopoietic progenitors from human pluripotent stem cells. STAR Protoc 1 100130 PMID: [33377024](https://pubmed.ncbi.nlm.nih.gov/33377024/)