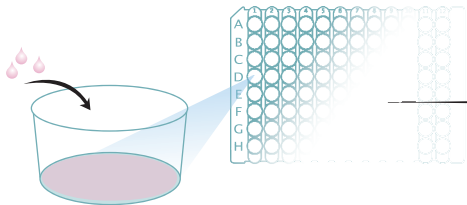


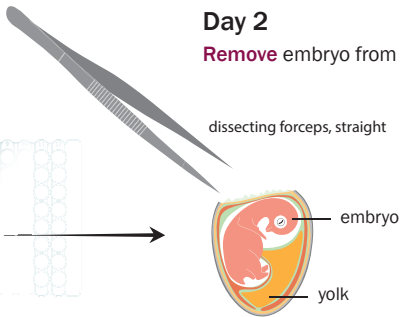
## Day 1

**Prepare** cell culture plates by coating with Mouse Laminin I or Bovine Fibronectin Protein.

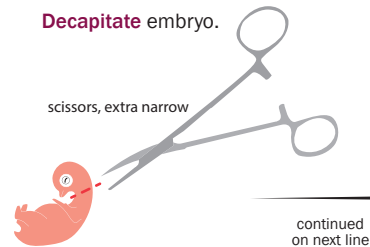


## Day 2

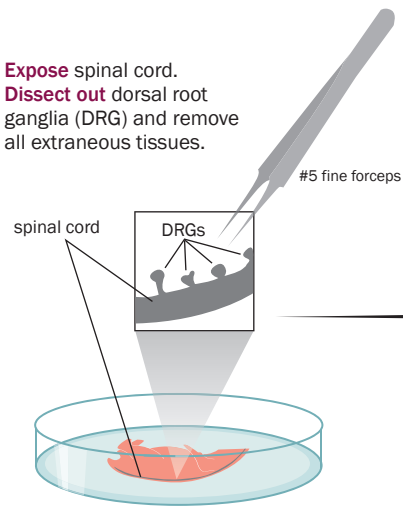
**Remove** embryo from egg .



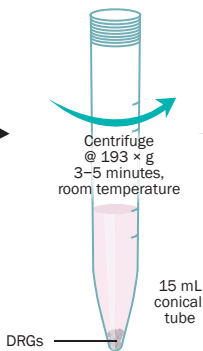
**Decapitate** embryo.



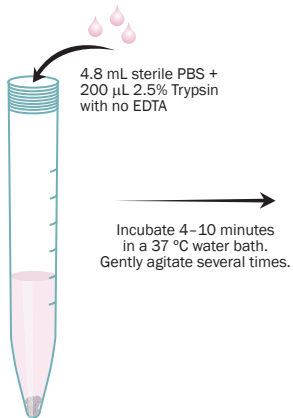
**Expose** spinal cord.  
**Dissect out** dorsal root ganglia (DRG) and remove all extraneous tissues.



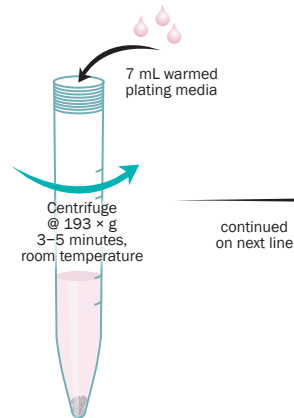
**Collect** cleaned DRGs in cold dissection media.  
**Centrifuge. Decant** supernatant.



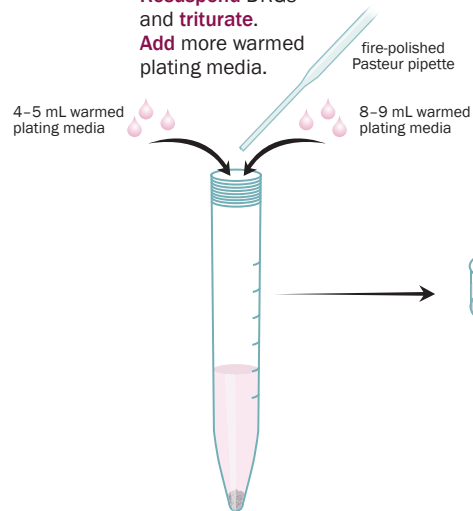
**Resuspend** DRGs.  
**Mix** by gentle agitation.



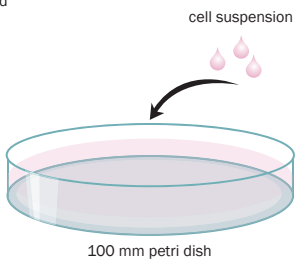
**Add** warmed plating media.  
**Centrifuge. Decant** supernatant.



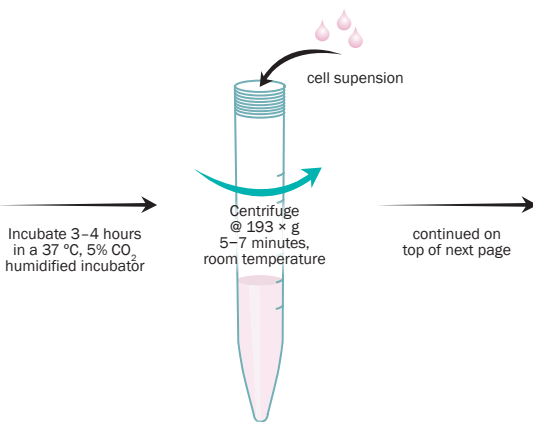
**Resuspend** DRGs and triturate.  
**Add** more warmed plating media.



**Transfer** cell suspension.



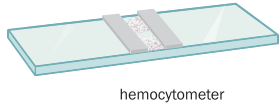
**Transfer** DRG neurons.  
**Centrifuge. Decant** supernatant.



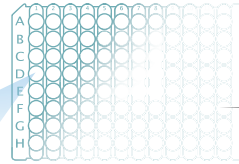
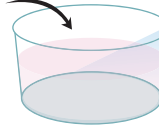
**Resuspend** DRGs neurons  
in 2–5 mL warmed culture media.  
**Count** cells.

**Reconstitute** DRG neurons with warmed  
culture media. **Seed** neurons onto  
Laminin I/Fibronectin-coated cell culture plates.

continued from  
previous page



100  $\mu$ L cell suspension



continued  
on next line

**Culture** DRG neurons for desired  
amount of time. **Exchange** media  
every 3–4 days.

culture media

