

View cultures and assess

```
graph TD; A[View cultures and assess] --> B[Centrifuge 150g for 5 minutes (pH of medium must be acidic)]; B --> C[Add 10-20% conditioned media to fresh media]; C --> D[Take sample of cells]; D --> E[Calculate cells/ml and reseed according to recommended density]; E --> F[Repeat every 2-3 days];
```

Centrifuge 150g for 5 minutes (pH of medium must be acidic)

Add 10-20% conditioned media to fresh media

Take sample of cells

Calculate cells/ml and reseed according to recommended density

Repeat every 2-3 days