

## Hybridoma antibody production

