

# Hybridoma antibody production

