

## OmicLink™ Expression Clone (CMV Promoter)

