

## OmicsLink<sup>™</sup> Expression Clone (CMV Promoter)

