

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

