

# Pathatrix® system—BAX® lysis linking step

