

# Pathatrix® system—generic bead beating protocol

