

Bio-Plex™ Protein Array System High Throughput Fluidics (HTF) System

For detailed instructions on the installation, set-up, and operation of the Bio-Plex HTF system, please refer to the Bio-Plex hardware manual.