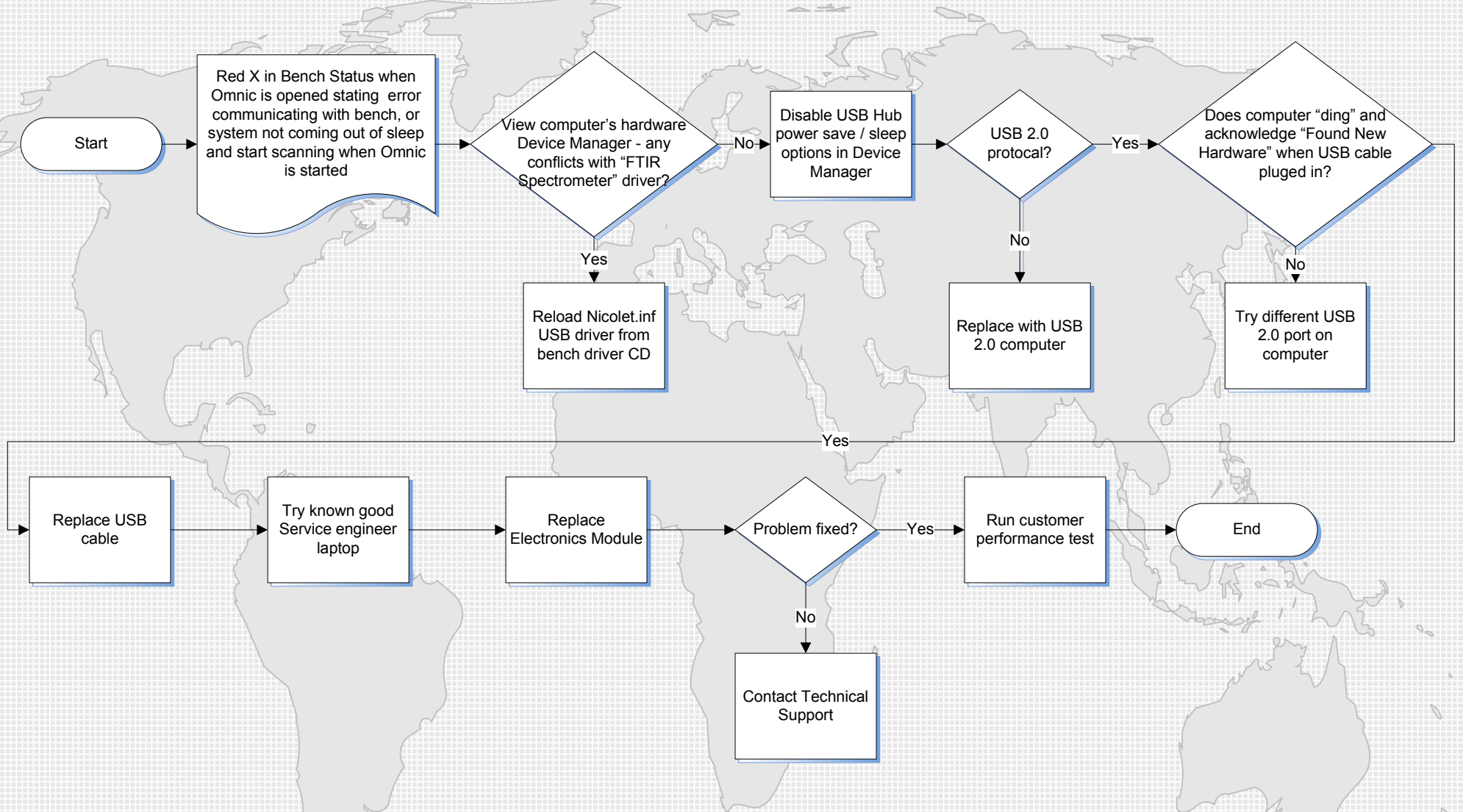


Basic troubleshooting FTIR/computer USB communication



Note

Any one of the fixes in the flow chart can resolve the USB communications problem. Once the communications problem is resolved, the interferometer should be auto aligned. Collect a customer performance test to verify instrument performance.