

LI-6400XT Photosynthesis System

OPEN 6.8.0a

LI-6850: connected

/ User 0% full simulator

Fri Apr 27 2012 11:43:36 12.46V

Home
Menu

Config
Menu

Calib
Menu

New
Measurements

Utility
Menu