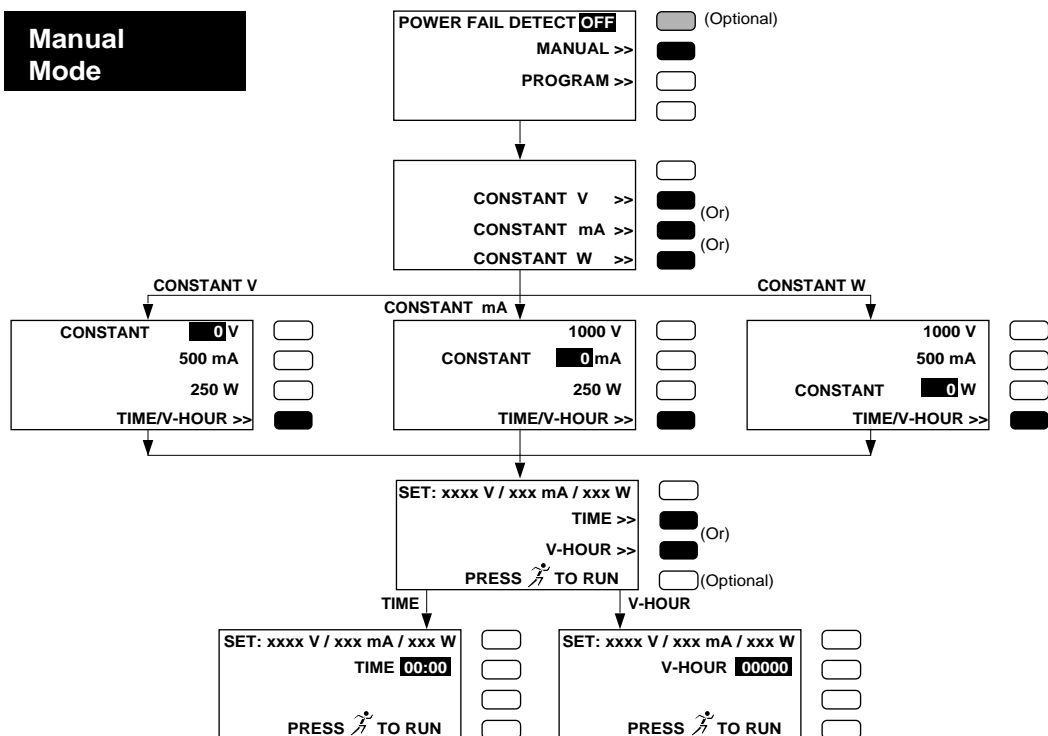


POWER/PAC 1000

QUICK REFERENCE GUIDE

Manual Mode



**Program Mode:
Creating a Method**

