

Troubleshooting Checklist



Steps:

1. State the problem

The Problem is _____

2. List All Possible Explanations

1. _____
2. _____
3. _____
4. _____
5. _____
6. _____
7. _____
8. _____
9. _____
10. _____

3. Collect the Data

Controls: _____
Storage: _____
Equipment: _____
Procedure: _____
Other: _____

4. Eliminate Explanations (Cross out in section 2)

5. Check with Experimentation (List your tests)

Exp 1: _____
Exp 2: _____
Exp 3: _____
Exp 4: _____
Exp 5: _____

6. Identify the Cause