

68H-0009.xml

▼ open

→ ▼ sensorhead = 6850

└─ address = "172.24.81.62"

▶ light

▶ comps

▶ prompts



◆ Edit



◆ Done