

6 8 H - 0 0 0 9 . x m l

L ▼ * o p e n

→ * s e n s o r h e a d = 6 4 0 0

▶ l i g h t

▶ c o m p s

▶ p r o m p t s

▶ c o n s t a n t s



◆ E d i t



◆ R e v e r t

◆ D o n e