

68H-0009.xml



*

light

source="Automatic"



*

par_in="Measured in Source"



*

sensor=""



*

cal=0.00



*

transm=0.90



/



Revert



Done