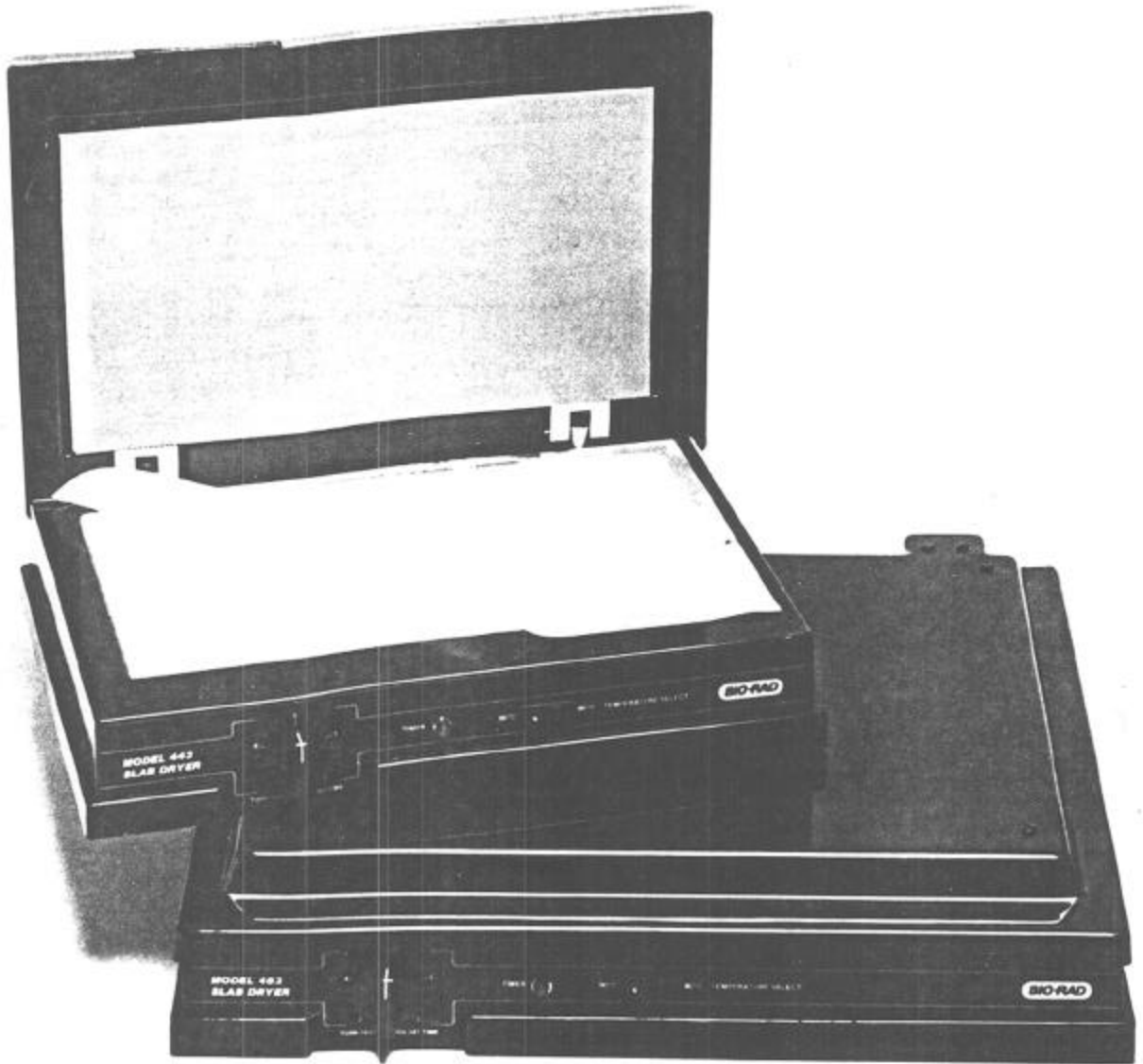


# **Model 443 and Model 483 Slab Dryer**

## **Instruction Manual**



**BIO-RAD**

