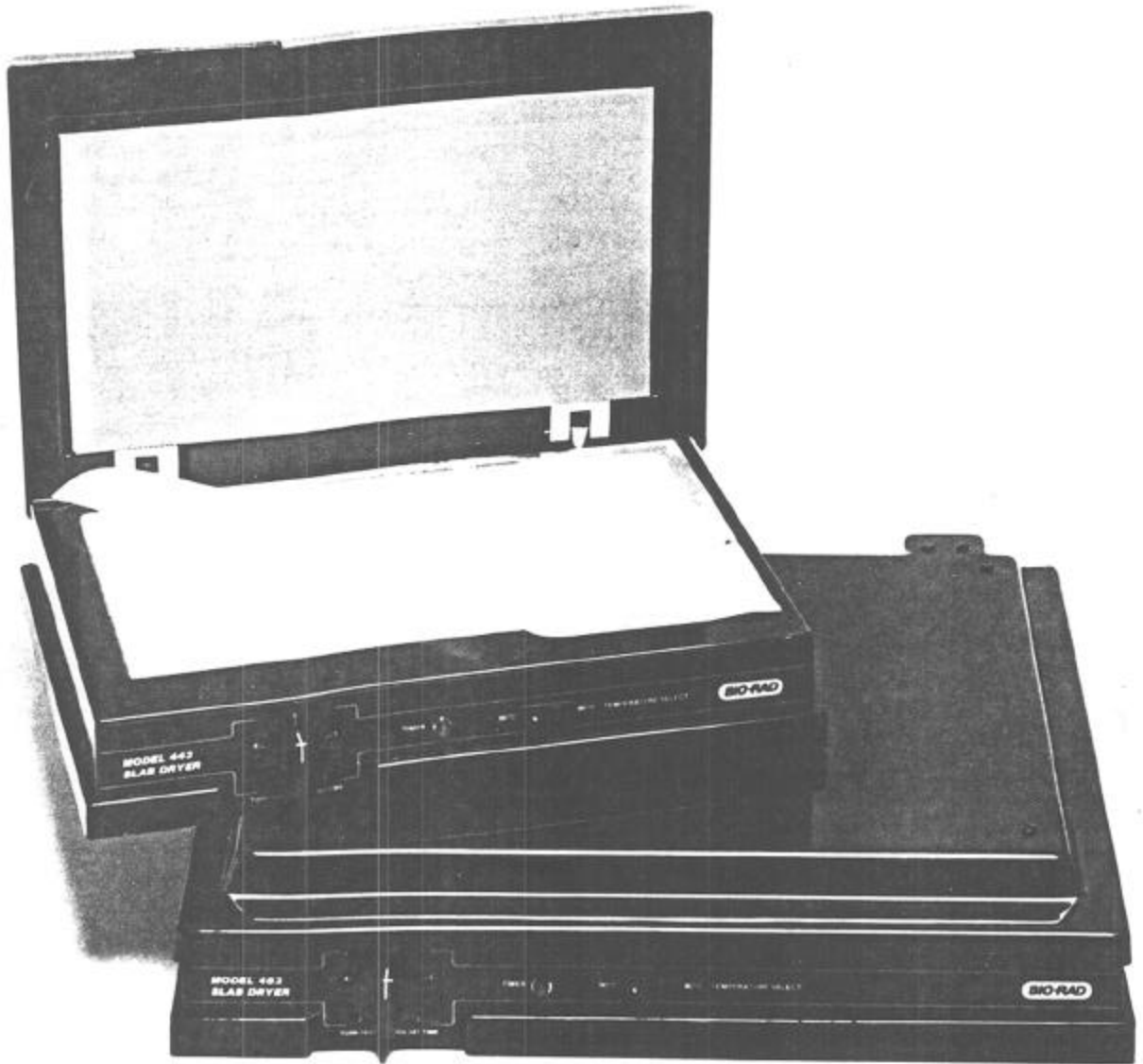


Model 443 and Model 483 Slab Dryer

Instruction Manual



BIO-RAD

