



## **Troubleshooting guide for Eppendorf Minispin - Z605220 and Z606235**

ERROR	CAUSE	SOLUTION
NO DISPLAY	NO MAIN POWER CONNECTION  POWER FAILURE	PLUG IN MAINS CABLE ON BOTH SIDES  CHECK MAINS FUSE OF THE PLUG/LAB
LID CANNOT BE OPENED	POWER FAILURE  ROTOR IS STILL SPINNING	USE EMERGENCY LID RELEASE  WAIT FOR ROTOR TO COME TO A STANDSTILL
CENTRIFUGE SHAKES ON ACCELERATION	ROTOR NOT LOADED SYMETRICALLY	STOP CENTRIFUGE RUN AND LOAD CENTRIFUGE SYMETRICALLY
“LID”	LID NOT CLOSED CORRECTLY	PRESS LID CLOSED. ERROR WITH LID CLOSING MECHANISM.
“INT”	POWER FAILURE DURING RUN	CHECK THE MAINS PLUG AND RESTART THE CENTRIFUGE USING START/STOP
“Err 5”	ROTOR LOADED INCORRECTLY  DEFECTIVE LID LATCH	CHECK LOADING OF ROTOR
“Err 9” TO “Err 14”	ELECTRONICS ERROR	SWITCH OFF CENTRIFUGE AND THEN SWITCH ON AGAIN
“Err 60” TO “Err 68”	DRIVE ERROR	ALLOW CENTRIFUGE TO COOL DOWN THEN RE-START