

# **Spare Parts** Manual

**Cardinal 8V**

**PN: 12008809-01**

## SPARE PARTS

<b>Cabinet Assembly</b>	<b>7-2</b>
<hr/>	
<b>Main Assembly</b>	<b>7-3</b>

**Spare part numbers and spare parts are subject to change in case of technical upgrades or supplier change. New specifications and part numbers are announced through technical service bulletins.**

### Cabinet Assembly

50149856 Outer Lid Assy'

20050543 Window

50146222 Gasket

50146001 Housing

20059405 Overlay Cardinal 8V

50147717 Front Panel

50148242 Nut

50150654 Collar (1pc)

50147732 Axis

50148240 Spring

50148656 Bushing

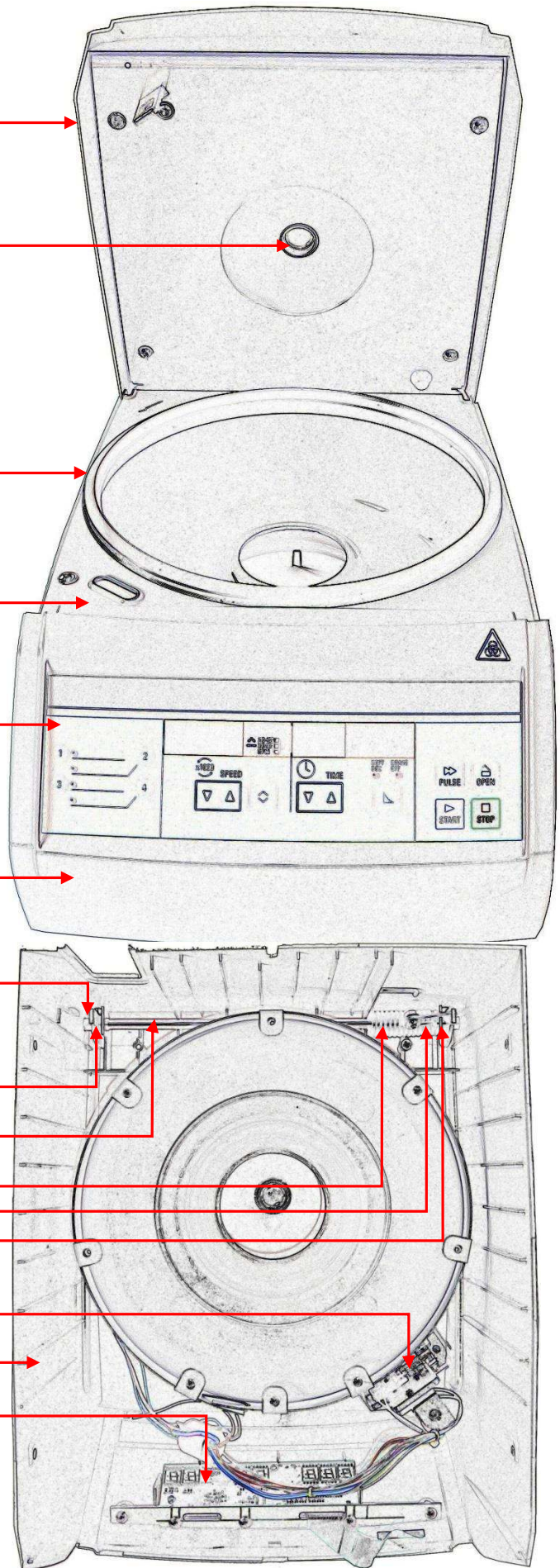
50150654 Collar (1pc)

20901162 Lid Lock Assy'

50146001 Housing

20150336 Display PCB

50104102 Display Ribbon Cable



### Main Assembly

