

Spare Parts Manual

Cardinal 8HC

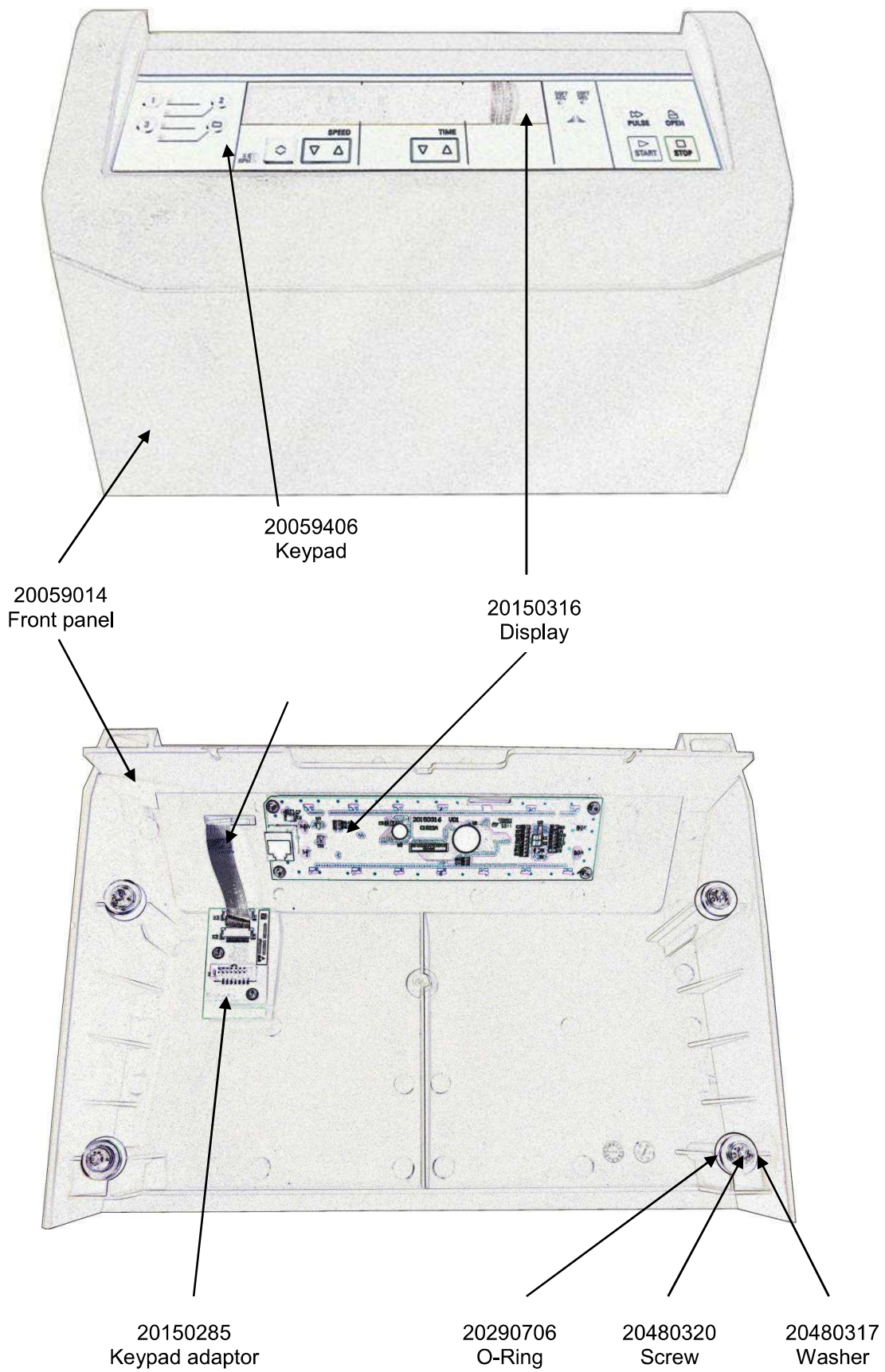
PN: 12007215-01

SPARE PARTS

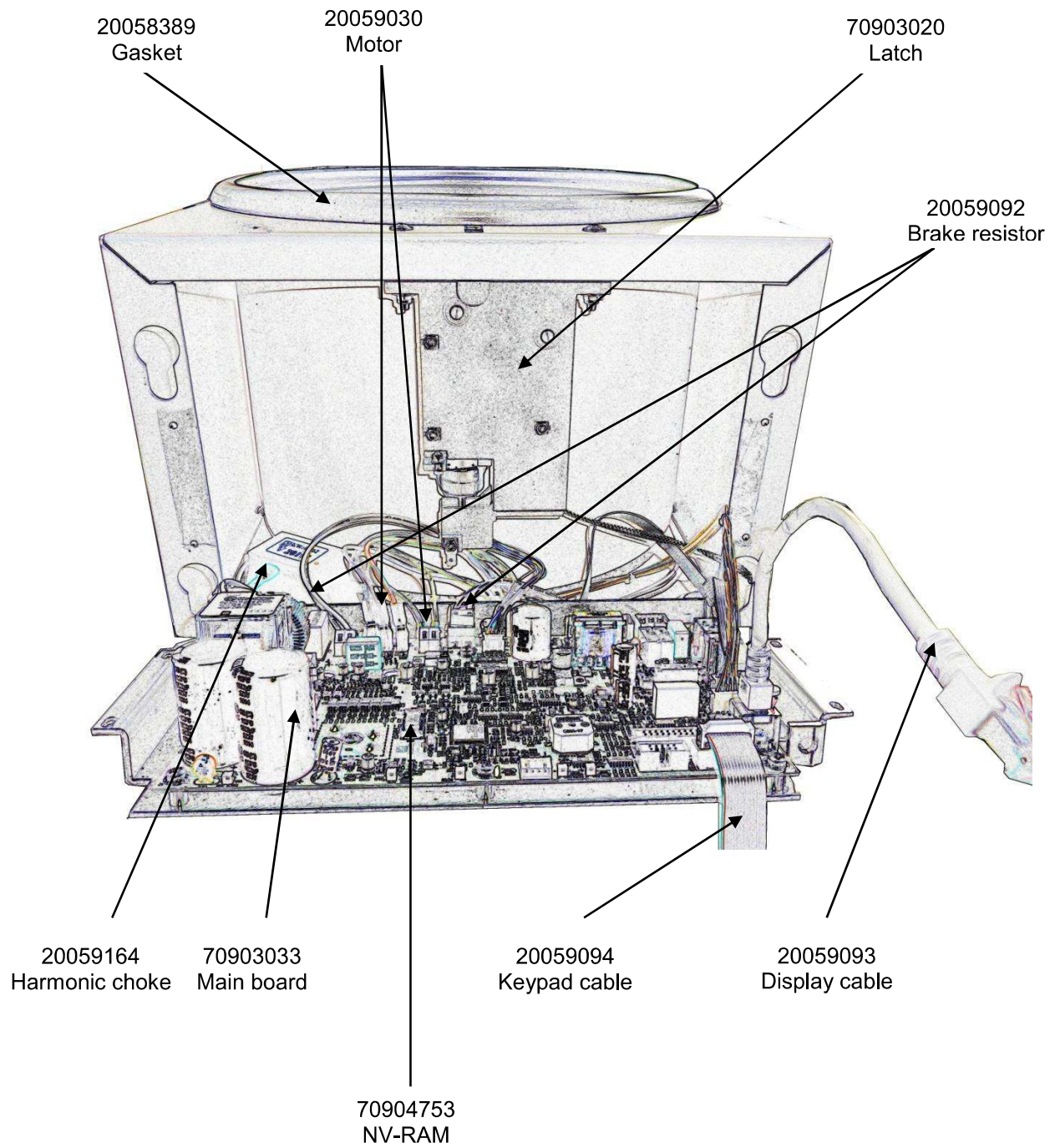
Front Panel	7-2
Front View	7-3
Back View	7-4
Top View	7-5
Drive	7-6

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.

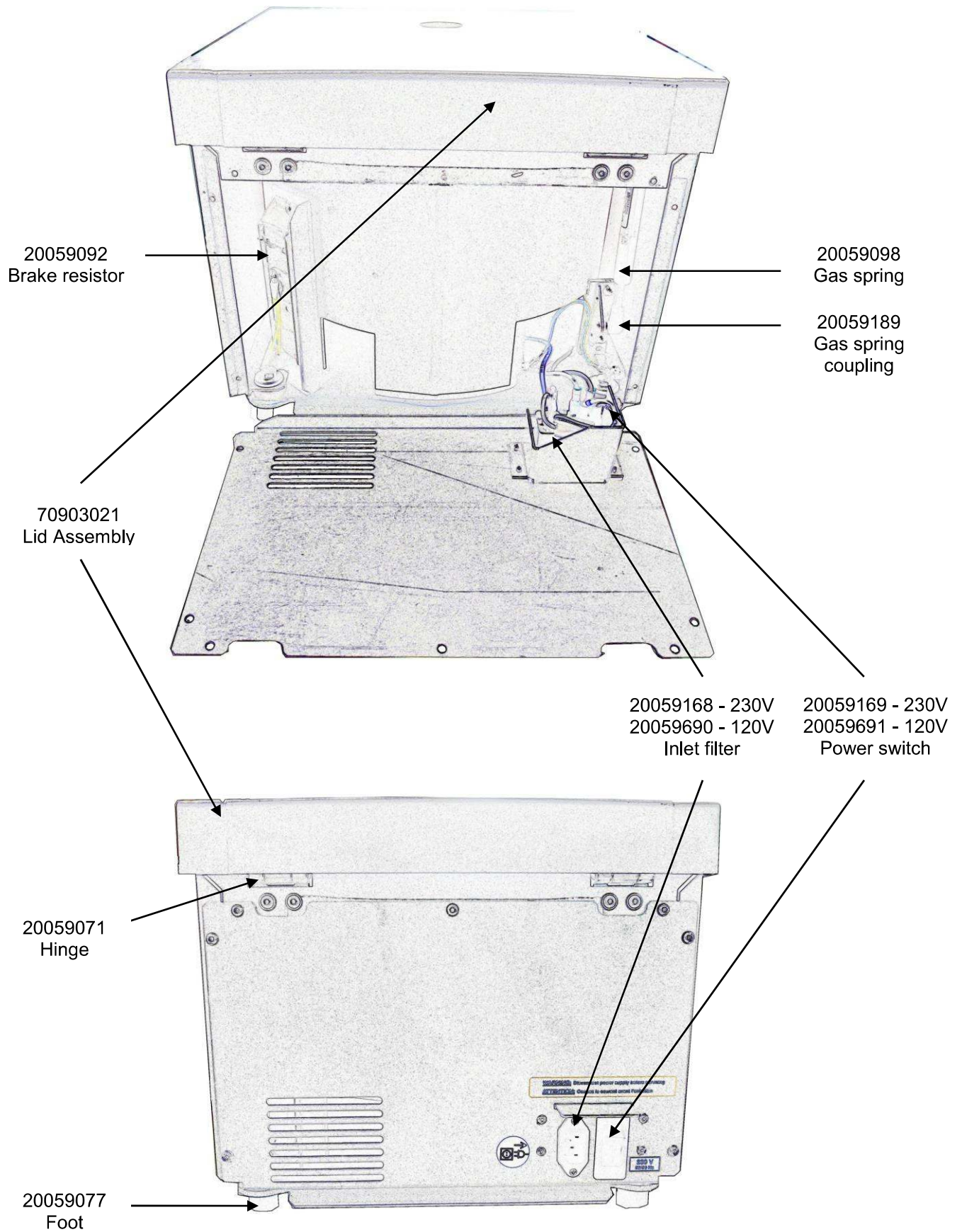
Front panel



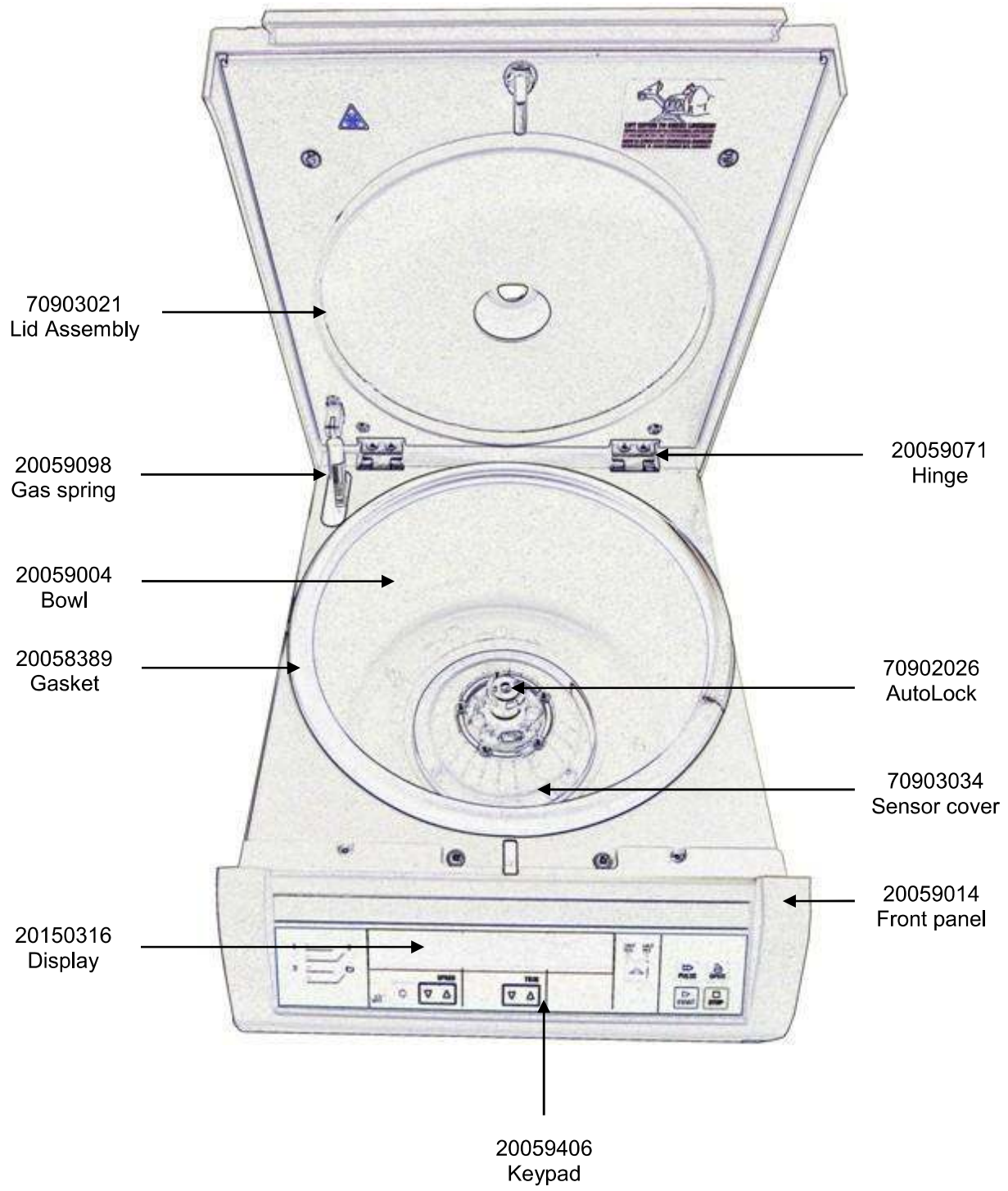
Front view



Back view



Top view



Drive

