

Large Cabinet Spec Sheet

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	80H x 40W x 27D in	203.2H x 102W x 68.6D cm
Internal Dimensions	66.5H x 33W x 24D in	166.4H x 84W x 61D cm
External Space	Place 3" from back wall	Place 76mm from back wall
Volume	30 ft ³	850 L
Weight	150 lbs	69 kg
Electrical Input	120VAC, 60Hz, 0.70A 240VAC, 50Hz, 0.35A	
Electrical Output	120VAC, 60Hz, 10A (10A only if connected device has its own power protection) 240VAC, 50Hz, 5A (5A only if connected device has its own power protection)	
Power Consumption	33W max	
Battery	12VDC, 1.3A, Sealed lead acid, qty 2	
Security Device	Strike Latch Lock	
Access Control	EM4102/EM4200 Proximity Reader	
Environmental Protection	IPX0	
Sound Level	37dB	
Communications	Ethernet based LAN DSL Option Cellular Option	
RFID Frequency	USA - 902-928 MHz EU 865-868 MHz	