

Large Cabinet

Model: TS030, TS030-P

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	81H x 44W x 30D in Place 3" from back wall and 21" from front wall	206H x 90W x 103D cm Place 76mm from back wall and 534mm from front wall
Volume	30ft ³	850L
Weight	200lb	91kg
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	Electromagnetic Lock x2	
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	

