

Small Ambient Cabinet Model: TS032, TS035, TS037, TS050

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	44H × 24W × 24D in	111.8H × 61W × 61.6D cm
Internal Dimensions	33H × 23W × 18D in	83.8H × 58.4W × 45.7D cm
External Space	44H × 24W × 48D in Place 3" from back wall and 21" from front wall	111.8H × 61W × 12.2D cm Place 76mm from back wall and 534mm from front wall
Volume	7.9ft³	224L
Weight	150lb	69kg
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	IPXO	
Sound Level	37dB	
Communications	Ethernet-based LAN DSL Option Cellular Option	
RFID Frequency	USA 902-928 MHz EU 865-868 MHz	



