

Small Auto Defrost Freezer (EU)

Model: **TS047e**

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	39.5H x 23.5W x 26D in	100.3H x 59.6W x 66D cm
Internal Dimensions	27H x 21W x 16D in	68.6H x 53.4W x 40.6D cm
External Space	41.5H x 24W x 48D in Place 3" from back wall and 23" from front wall	105.4H x 61W x 122D cm Place 76mm from back wall and 534mm from front wall
Volume	5.5ft ³	156L
Weight	130lb	59kg
Electrical Input	230VAC, 50Hz, 1.0A	
Electrical Output	230VAC, 50Hz, 5A	
Power Consumption	150W max	
Battery	12VDC, 1.3A, Sealed Lead Acid, qty 2	
Security Device	Electromagnetic Lock	
Access Control	EM4102/EM4200 Proximity Reader	
Environmental Protection	IPX0	
Sound Level	33.4dB	
Communications	Ethernet-based LAN DSL Option Cellular Option	
RFID Frequency	EU 865–868 MHz	

