

# Small Freezer

Model: **TS036, TS038**

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	41.5H x 24W x 24D in	105.4H x 61W x 61D 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 21" from front wall	105.4H x 61W x 119.5D cm Place 76mm from back wall and 534mm from front wall
Volume	5.5ft <sup>3</sup>	156L
Weight	150lb	69kg
Electrical Input	120VAC, 60Hz, 1.00A	
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	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	37dB	
Communications	Ethernet-based LAN DSL Option Cellular Option	
RFID Frequency	USA 902–928 MHz	

