

# Large Freezer

Model: **TS078, TS031, TS042, TS043**

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	71H x 24W x 27D in	180.4H x 61W x 66D cm
Internal Dimensions	49H x 19.5W x 15D in	124.5H x 49.5W x 38D cm
External Space	71H x 24W x 50D in Place 3" from back wall and 21" from front wall	180.4H x 61W x 127D cm Place 76mm from back wall and 534mm from front wall
Volume	8.3ft <sup>3</sup>	235L
Weight	200lb	91kg
Electrical Input	120VAC, 60Hz, 1.00A 240VAC, 50Hz, 0.50A	
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	180W 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 EU 865–868 MHz	

