

ULT Freezer

Model: **TS057**

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	78H x 38W x 37D in	198H x 96.5W x 94D cm
Internal Dimensions	52H x 23W x 19D in	132H x 58.4W x 48.3D cm
External Space	83H x 52W x 81D in Place 6" from back wall Allow min. 34" for door clearance	210.81H x 132W x 205.7D cm Place 15cm from back wall Allow min. 86cm for door clearance
Volume	13ft ³	368L
Weight	725lb	329kg
Temperature Range	-85°C to -50°C	
Refrigerant	CFC/HCFC-free environmentally safe refrigerant mixtures/non-flammable in both stages	
Cooling Method	2-stage cascade system, CO ₂ and LN ₂ options available	
Electrical Output	120VAC, 60Hz, 16FLA 230VAC, 60Hz, 12FLA	
Power Consumption	1.9KW max	
Battery	12VDC, 1.3A, Sealed Lead Acid, qty 2	
Security Device	Electromagnetic Lock	
Access Control	LF Proximity Reader	
Environmental Protection	IP41, RoHS-compliant	
Sound Level	37dB	
Communications	Ethernet-based LAN WiFi Option Cellular Option	
RFID Frequency	USA 902-928 MHz	

