

# ULT Under Counter Freezer

Model: **TS079**

Attribute	Specification (Imperial)	Specification (Metric)
External Dimensions	32.7H x 35.4W x 22.3D in	83.1H x 89.9W x 56.6D cm
Internal Dimensions	25.2H x 18.9W x 13D in	64H x 48W x 33D cm
External Space	32.7H x 40W x 50D in	83.1H x 101.6W x 127D cm
Volume	3.6ft <sup>3</sup>	102L
Weight	265lb	120kg
Maximum Shelf Weight	44lb	20kg
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 <sub>2</sub> and LN <sub>2</sub> options available	
Electrical Output	120VAC, 60Hz, 5A	
Power Consumption	367W 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	33.4dB	
Communications	Ethernet-based LAN WiFi Option Cellular Option	
RFID Frequency	USA 902-928 MHz	

