

Large Biomedical Freezer Spec Sheet

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	Place 3" from back wall	Place 76mm from back wall
Volume	8.3ft ³	91 L
Weight	200 lbs	91 kg
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	180W max	
Battery	12VDC, 1.3A, Sealed lead acid, qty 2	
Security Device	Electromagnetic Lock	
Access Control	EM4102/EM4200 Proximity Reader	
Environmental Protection	IPXO	
Sound Level	37dB	
Communications	Ethernet based LAN DSL Option Cellular Option	
RFID Frequency	USA - 902-928 MHz	

