

Refrigerator Spec Sheet

| 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 54.3W x 40.6D cm |
| External Space | Place 3" from back wall | Place 76mm from back wall |
| Volume | 5.5ft ³ | 156 L |
| Weight | 150 lbs | 69 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 | 150W 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 | |

