

## Small Cabinet Spec Sheet

| Attribute                | Specification (Imperial)   | Specification (Metric)    |
|--------------------------|--|---------------------------|
| External Dimensions      | 44H x 24W x 24D in   | 11.8H x 61W x 61.6D cm    |
| Internal Dimensions      | 33H x 23W x 18D in   | 83.8H x 58.4W x 45.7D cm  |
| External Space           | Place 3" from back wall  | Place 76mm from back wall |
| Volume                   | 7.9ft <sup>3</sup>   | 224 L                     |
| Weight                   | 150 lbs  | 69 kg                     |
| Electrical Input         | 120VAC, 60Hz, 0.70A<br>240VAC, 50Hz, 0.35A   |                           |
| Electrical Output        | 120VAC, 60Hz, 10A (10A only if connected device has its own power protection)<br>240VAC, 50Hz, 5A (5A only if connected device has its own power protection) |                           |
| Power Consumption        | 33W max  |                           |
| Battery                  | 12VDC, 1.3A, Sealed lead acid, qty 2   |                           |
| Security Device          | Strike Latch Lock  |                           |
| Access Control           | EM4102/EM4200 Proximity Reader   |                           |
| Environmental Protection | IPX0   |                           |
| Sound Level              | 37dB   |                           |
| Communications           | Ethernet based LAN<br>DSL Option<br>Cellular Option  |                           |
| RFID Frequency           | USA - 902-928 MHz<br>EU 865-868 MHz  |                           |

