

# Countertop RFID Refrigerator

Model: **TS098**

## Benefits of the TS098-Countertop RFID Refrigerator

- Maintain temperature range of +1°C to +10°C
- Prevent stock-outs; monitor expiry of products
- Customize product layout and branding
- Easily installable on a standard countertop
- Protect critical inventory through secure access
- Provide regulatory compliance
- Integrate with third-party temperature monitoring system

Attribute	Specification
External dimensions	23in H x 17.25in W x 19.25in D
Internal dimensions	16in H x 14in W x 8in D
External clearance	23in H x 25in W x 35in D
Unit volume	1.0ft <sup>3</sup>
Unit weight	64lb
Shelf weight capacity	10lb per shelf
Door	Solid, Right-side hinge
Operating temperature range	+1°C to +10°C/+5°C setpoint
Temperature uniformity	1.5°C
Temperature stability	+/- 1.0°C
Refrigerant	R600a (SNAP)
Cooling method	Forced-air-cooled
Defrost type	Auto
Security device	Electronic lock
Access control	LF/HF
<b>Communications</b>	
RFID frequency range	902MHz to 928MHz
Connection	Ethernet, optional cellular
<b>Power</b>	
Electrical power and frequency	120VAC, 60Hz, 2A
Power consumption	1.1kWhr/day
RFID battery backup	24V SLA 1.3Ah

