

Safety and security radiation monitoring

Scenario:

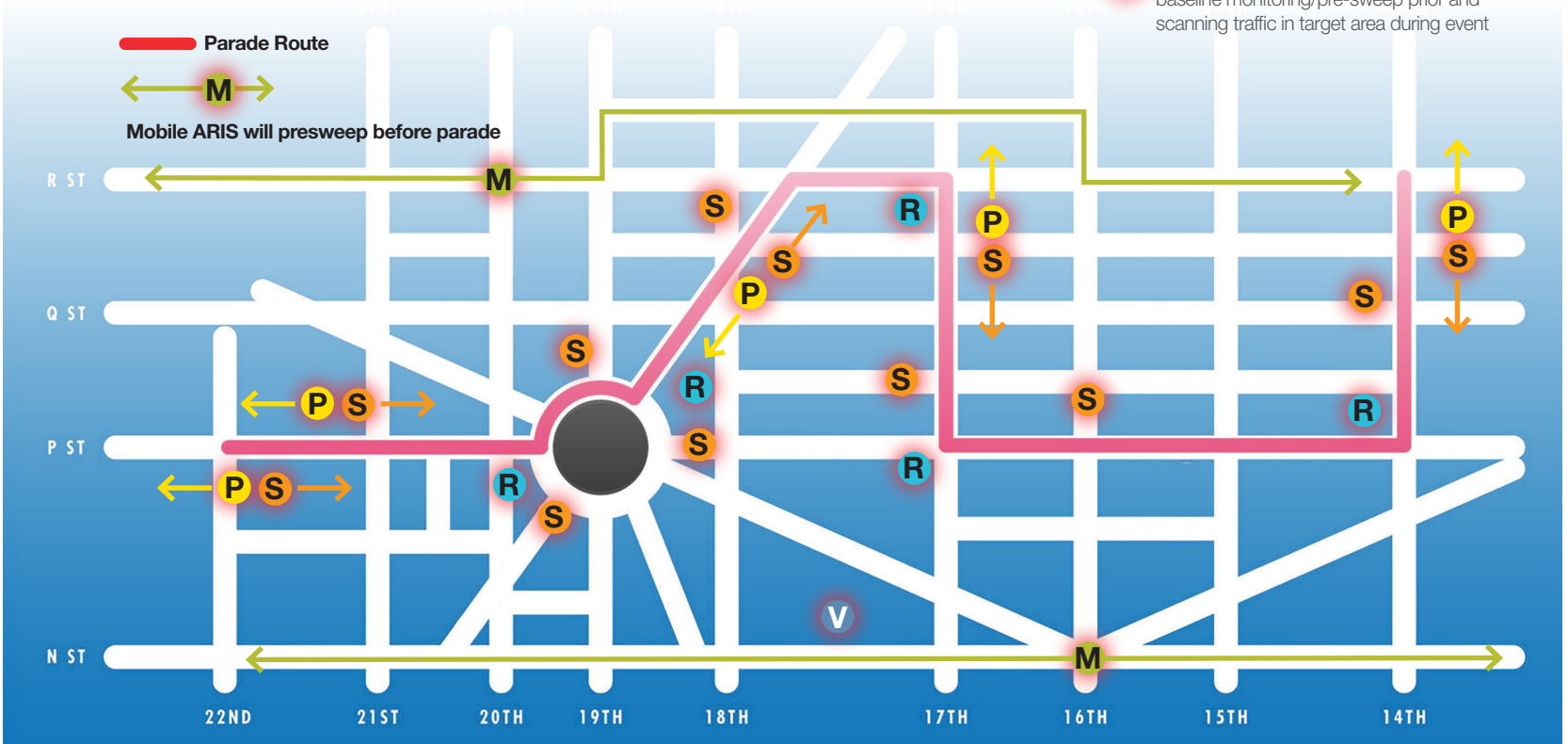
Monitor all traffic and fans along the parade route. Baseline scans prior to the event and monitor for changes during event.

Areas of concern:

- Size of area to monitor with limited checkpoints for screening and monitoring
- Fans carrying sources along route
- Sources placed in high traffic areas
- Vehicles in motion armed with devices

Key products and set up:

- P** Thermo Scientific™ PackEye™ Radiation Detection Backpack - baseline scanning prior and event time comparison
 - Start and finish lines
 - Walking the route
- R** Thermo Scientific™ RadHalo™ RDP Area Monitor - screening spectators at large funnel points
- V** Network and Remote Monitoring
 - Thermo Scientific™ ViewPoint™ Enterprise Remote Monitoring - command and control software to monitor all devices
 - SEE Multimeter Viewer - cloud based solution setup in minutes with built in plume tracker
- S** Thermo Scientific™ RadEye SPRD™ Spectroscopic Personal Radiation Detector - police officers wearing as primary means of locating source as they walk the event
- M** Thermo Scientific™ Mobile ARIS™ - baseline monitoring/pre-sweep prior and scanning traffic in target area during event



To register, please visit www.thermofisher.com/radiationsecurity

