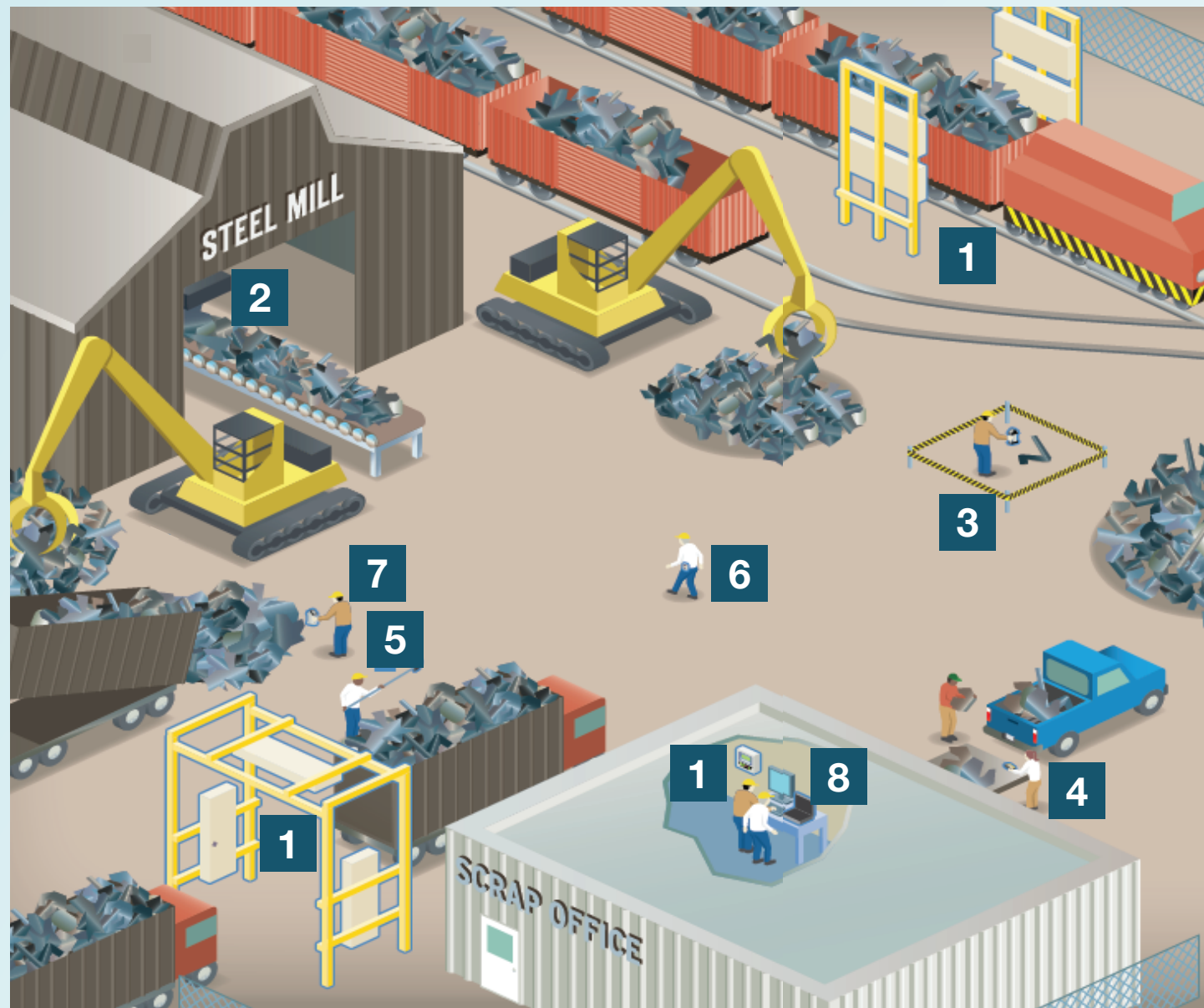


8 Reasons metal processors need radiation detection



- Radioactive sources can frequently show up at recycling facilities.
- These sources threaten works and can end up in recycled material, threatening consumers.
- Multiple points of inspection are needed – from vehicles entering to various steps in the process.

We have identified **8 key places** to consider when looking to protect your facility and the public from radiation threats.

- 1 Truck and Railcar Monitoring** Use Radiation Portal Monitoring systems for truck and railcar monitoring. **ASM IV Series monitoring system**
- 2 Charge Bucket, Conveyor, Platform Scales or Dust Collection** Ensure process monitoring systems are configurable for conveyor, platform scales or system dust collection. **ASM IV Charge Bucket monitoring system**
- 3 Secondary Inspection** Use the sensitive and fast responding **RadEye NBR High Sensitivity Gamma Radiation Monitor** to verify the radioactive find and assess if of natural or artificial (man-made) origin. Alternatively, or in addition use the ruggedized **RIIDEyeX Handheld Radioactive Isotope Identifier** to provide fast and real time identification and analysis.
- 4 Combination Detection** Belt-worn radioactive gamma isotope identification and neutron indication. **RadEye SPRD Spectroscopic Personal Radiation Detector**
- 5 Search and Find Applications** Portable Personal Radiation Detectors provide sensitive and fast detection of gamma radiation with accurate dose rate measurements. **RadEye PRD4 Personal Radiation Detector**
- 6 Gamma Neutron Paging** Use a monitor that combines gamma sensitivity and energy compensated dose rate measurement with separate, high sensitivity neutron response and alarm threshold. **RadEye GN+ Gamma Radiation monitor**
- 7 Accurate Identification of Source** Quickly locate radioactive materials with the most sensitive hand-held instrument featuring fast discrimination between man-made artificial sources and natural radiation. **RadEye NBR**
- 8 Documentation** Utilize software for documenting scans performed. **RadEye Safety Kit and Viewpoint Enterprise Remote Monitoring System**

Find out more at thermofisher.com/radiationmeasurement