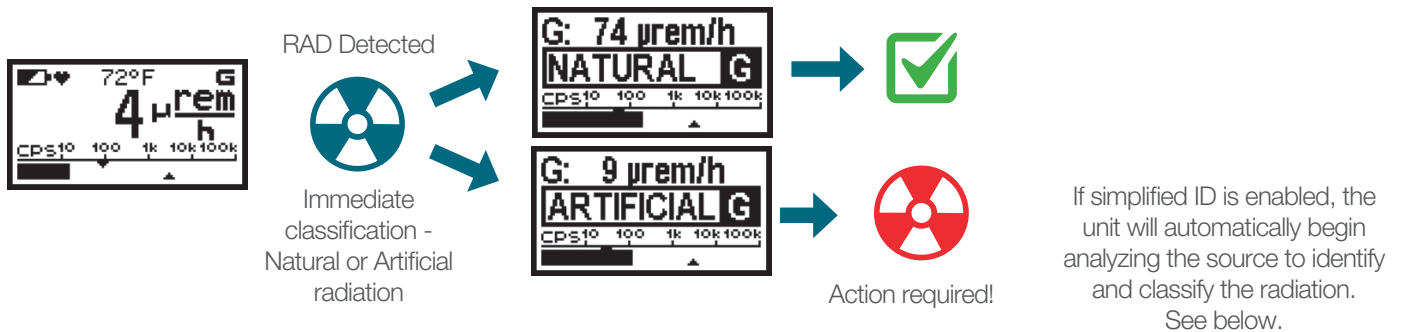


RadEye SPRD-GN - Quick Start Guide



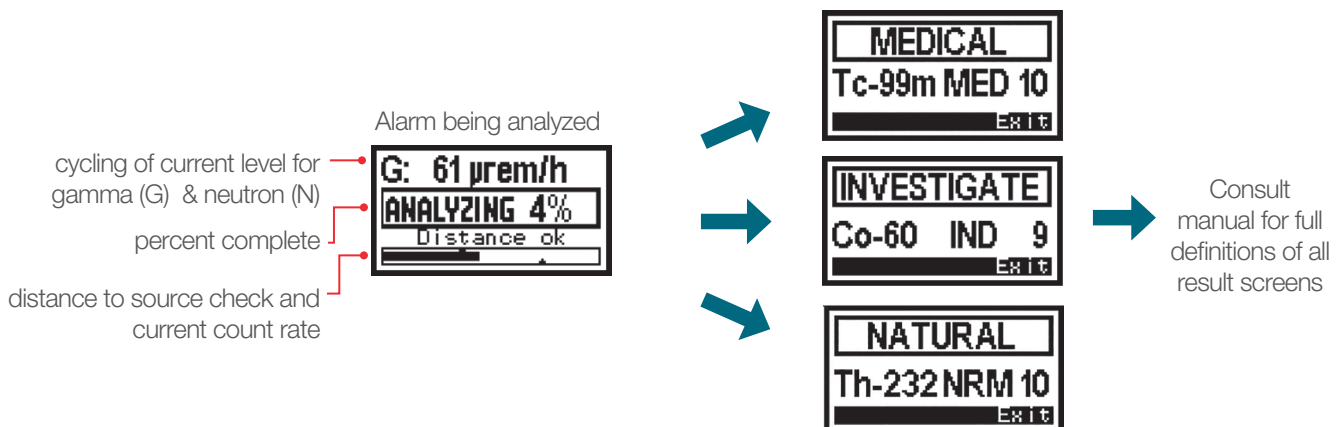
Ratemeter

Can be configured to display dose rate or count rate in the main display



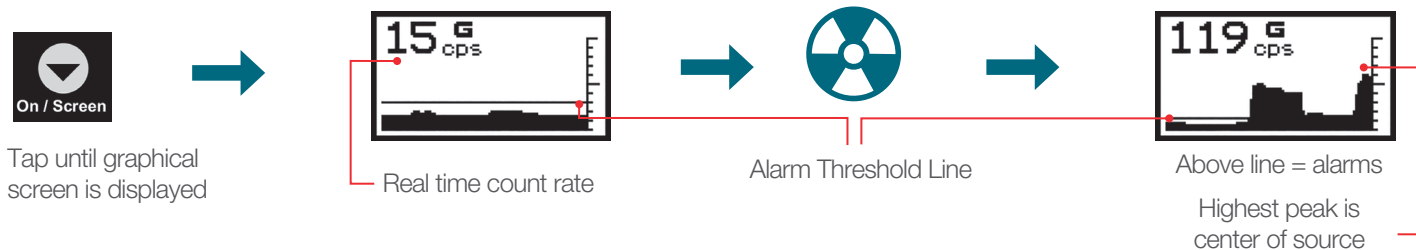
Identification and Classification

Automatically classify and ID gamma radiation after receiving an alarm



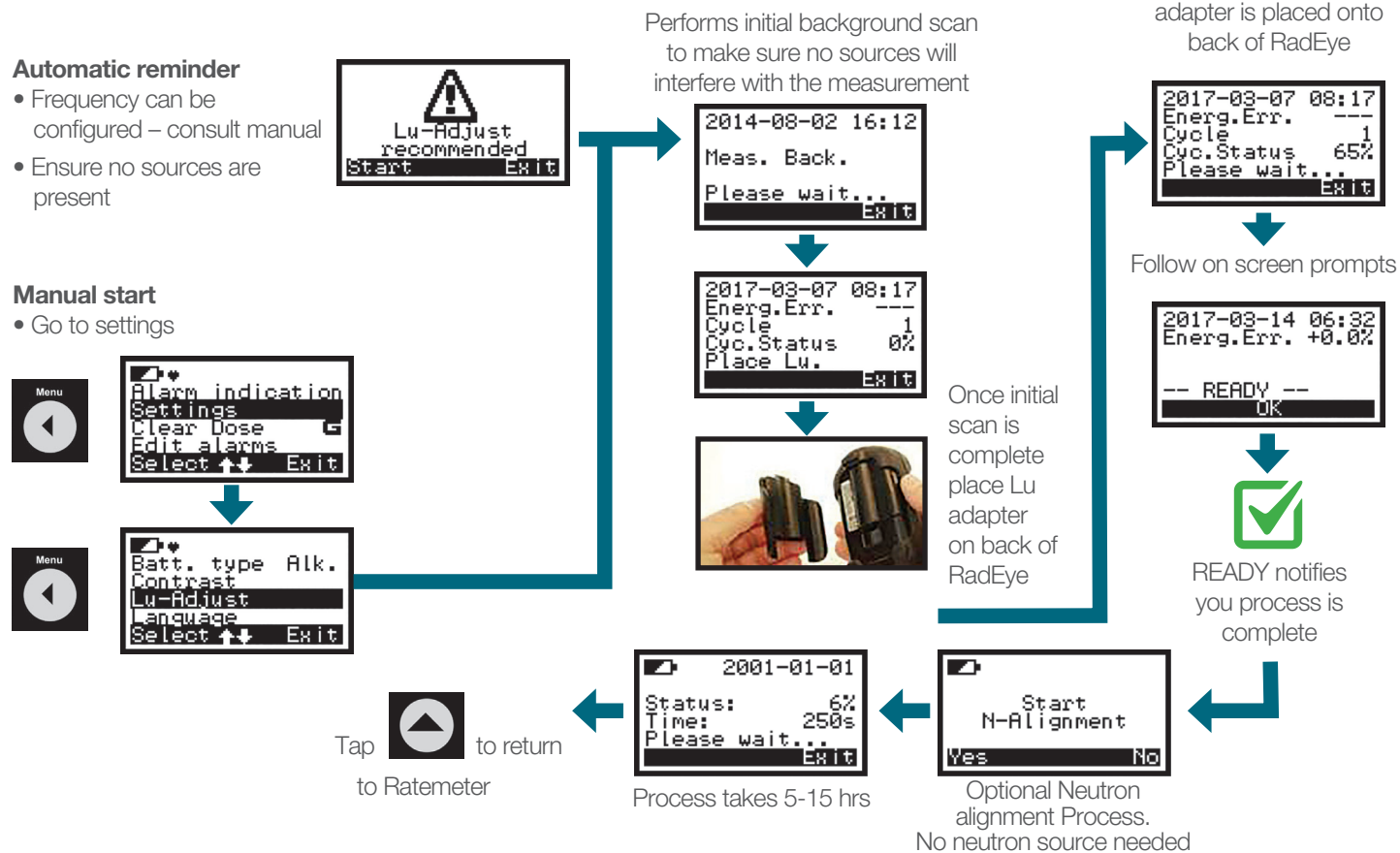
Finder Mode

Quickly pinpoint radiation sources using simple graphical display of source strength



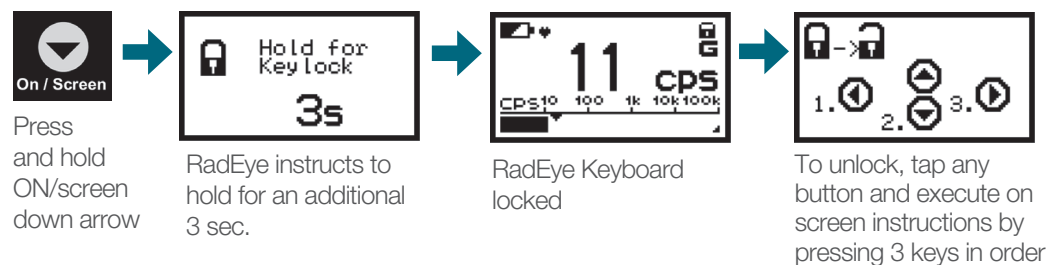
Optimization with Lu Test Adapter

Fast, field optimization to keep RadEye operating at its best

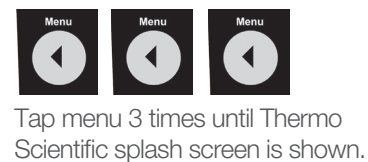


Locking Keyboard

Lock keyboard to prevent accidental operation while in a pocket



Power Off



ThermoFisher
SCIENTIFIC