

applied biosystems



## Quick reference guide: RapidHIT ID System

DNA results in 90 minutes

**ThermoFisher**  
SCIENTIFIC

# Applied Biosystems™ RapidHIT™ ID System investigative lead workflow

1

Obtain a sample



Buccal/cheek swab



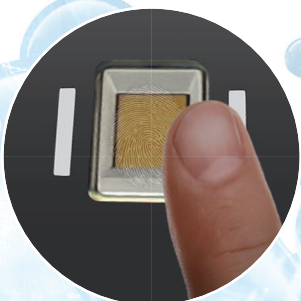
Crime scene  
samples

2

Sign in (3 options)



Facial  
recognition



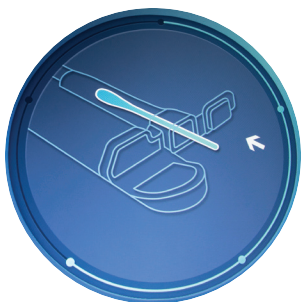
Fingerprint



Six-digit pin

3

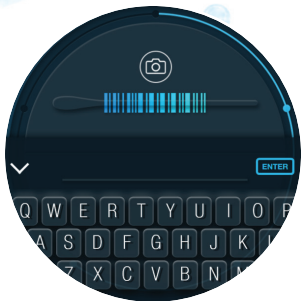
Insert cartridge and enter  
sample identification



Insert the sample into  
a sample cartridge



Instrument will read the RFID from the RFID chip



Scan barcode or manually  
enter sample name

4

Process sample on the instrument



For Applied Biosystems™ RapidINTEL™ Plus Cartridge, select instrument protocol



Confirm sample and run information and start run



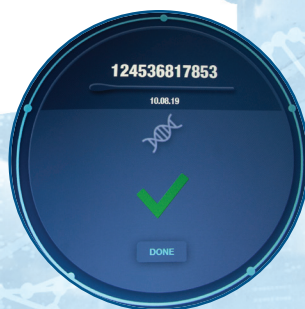
Results will be generated in ~90 minutes

5

Remove the sample cartridge



Discard or store the sample cartridge based upon standard operating procedures



A green checkmark indicates that a profile was generated, yellow indicates that the profile needs review by the lab, and red indicates there was no result



Review results in Applied Biosystems™ RapidLINK™ Software



## Two sample cartridges for various sample types

Self-contained cartridge technology uses standard laboratory steps to process samples.

### Applied Biosystems™ RapidINTEL™ Plus Sample Cartridge



### Applied Biosystems™ RapidHIT™ ID ACE GlobalFiler™ Express Sample Cartridge



### Workflow process at a glance

Collect sample

Process sample on location  
in 90 minutes or less

Get hit/no-hit  
confirmation report



Find out more at [thermofisher.com/rapidDNA](https://thermofisher.com/rapidDNA)

**ThermoFisher**  
SCIENTIFIC