# **Gold Purity Analysis: Portable XRF vs. Fire Assay**

Fire assay is a method widely used to determine the purity of gold in jewelry. We are often asked the question: "how does portable XRF compare to fire assay?" The answer is that the accuracy of Thermo Scientific™ Niton™ DXL and XL2 Precious Metal Analyzers is comparable to the accuracy of fire assay and is typically within 1/8 of a karat of the fire assay results. In addition to being a non-destructive method, portable XRF is a much more comprehensive analysis, yet simple, fast, and does not require any hazardous chemicals.



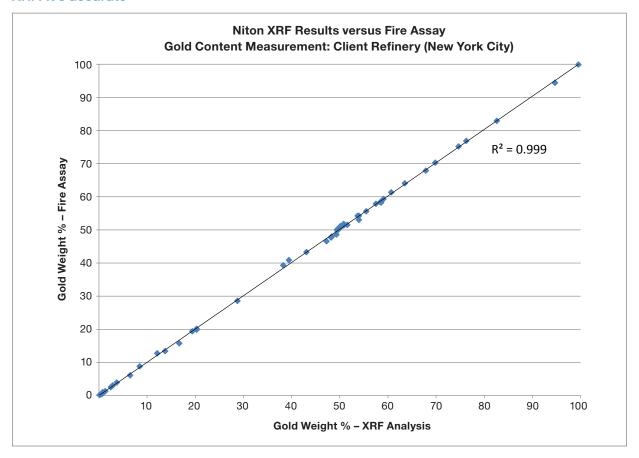
Niton DXL countertop analyzer

# XRF vs. Fire Assay

Feature	XRF	Fire Assay	XRF Advantages
Level of complexity	Simple	Complex	<ul><li>Easy to use</li><li>Requires less lab space</li></ul>
Analysis duration	3-10 seconds	2-3 hours	<ul><li>High throughput</li><li>Customer satisfaction</li></ul>
Eliminate the need for a skilled operator	✓	Х	Training costs
Eliminate the need for harsh chemicals	✓	Х	<ul><li>Employee safety</li><li>Environmental compliance</li><li>Cost of disposal</li></ul>
Portable	✓	Х	Transfer between locations
Non-destructive	✓	Х	<ul><li>Customer satisfaction</li><li>Value retention of jewelry</li></ul>
Number of elements analyzed	22 (incl. Au, Ag, Ni, W)	1 (Au)	Comprehensive information Indication about fraudulent merchandise



## XRF: It's accurate



# **Fire Assay**



# XRF: It's simple



## thermoscientific.com/niton

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

### Americas Boston, MA USA

+1 978 670 7460 niton@thermofisher.com

#### Europe, Middle East, Africa & South Asia Munich, Germany +49 89 3681 380 niton.eur@thermofisher.com

Asia Pacific New Territories, Hong Kong +852 2885 4613 niton.asia@thermofisher.com

In addition to these offices, Thermo Fisher Scientific maintains a network of representative organizations throughout the world.

