

SIMCA in Crime Scene Investigations

Dr. Laura Ortiz Herrero and Dr. Luis Bartolomé Moro, University of the Basque Country (UPV/EHU)

Umetrics Webinar May 2021

Outline

Introduction

- Importance of determining the time frame at the crime scene.
- Potential of chemometrics and current status of its implementation to forensic dating.
- Objective.

Methodology

- Proposed chemometric based dating methodology.

Case study

- Questioned documents field.

Conclusions

- General conclusions drawn from the thesis work.



When did an event **occur**?

When was a trace **left** on the crime scene?



How old is an evidence?

In what order did events **occur**?



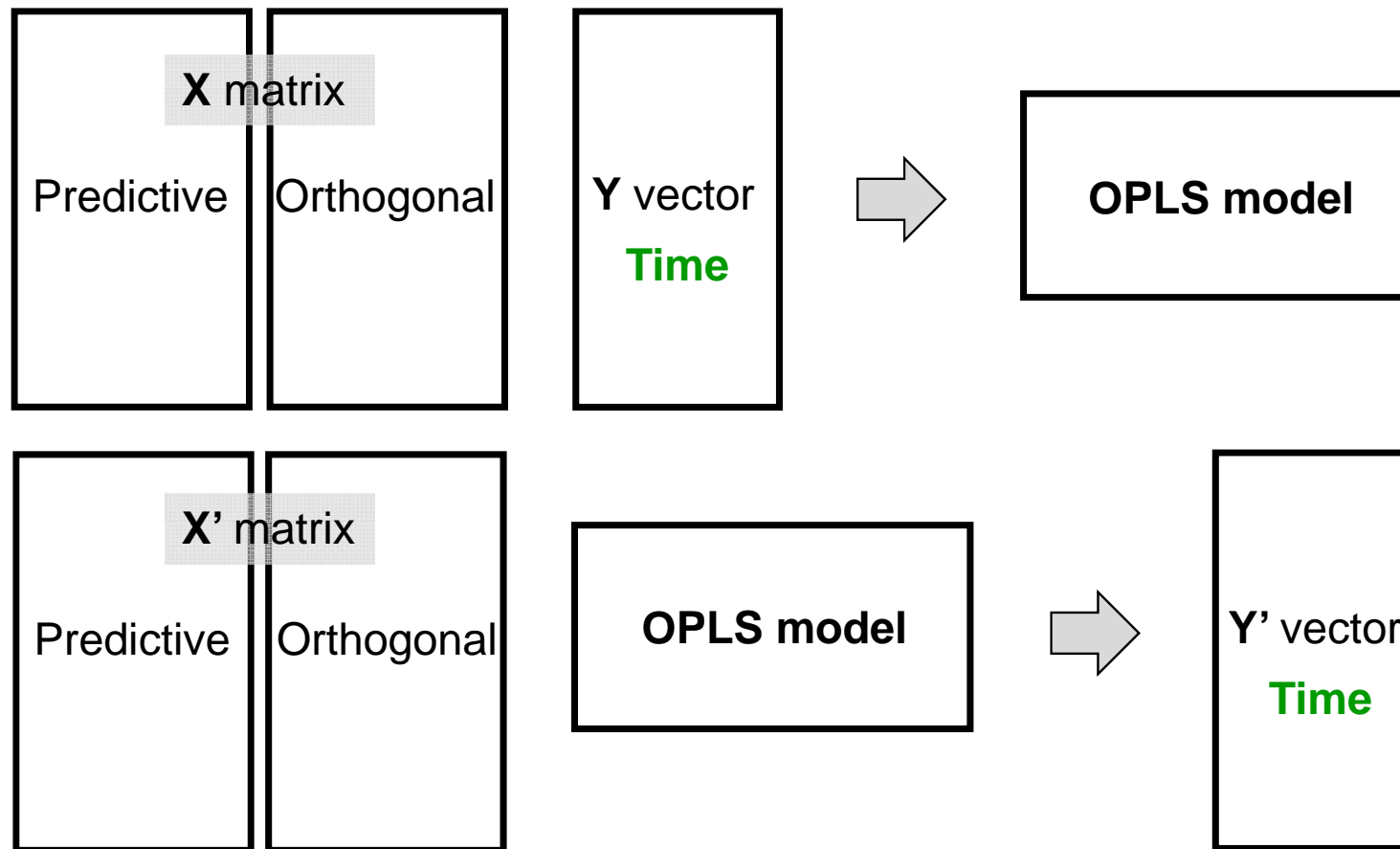
Time has been largely **unexplored** due to the complexity of the **overall challenge**.



combined with **chemometrics**

- ✓ Improves **handling** and **interpretation** of high dimensional data.
- ✓ Provides additional **information**.
- ✓ Achieves objective and meaningful **results**.
- ✓ **Fast** in the time domain.
- ✓ Reduces human **error**.

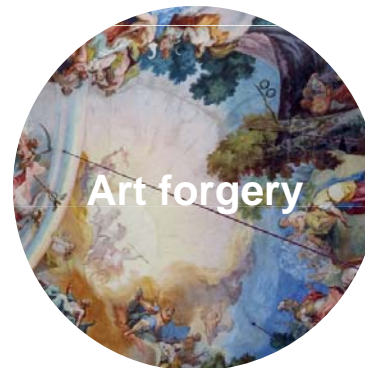
Partial least squares regression (PLSR) and its **extension OPLSR**.



Chemometrics has **recently** been **introduced** to forensic dating.

- × **Potential applicability** poorly understood and underexploited.
- × Lack of **statistical expertise** for its optimal application.
- × Limited implementation of **standardised** chemometric-based **methodologies** in forensic laboratories.

The **aim** of this thesis has been to **highlight the usefulness of multivariate regression methods** in the **forensic field** through the **development and validation of dating methodologies** in which non-destructive and micro-destructive techniques have been applied together with the (O)PLSR method.



Introduction

Methodology

Questioned documents

Conclusions

Journal of Analytical and Applied Pyrolysis 131 (2018) 9–16



Direct and indirect approaches based on paper analysis by Py-GC/MS for estimating the age of documents

L. Ortiz-Herrero^{a,*}, M.E. Blanco^a, C. García-Ruiz^{b,c}, L. Bartolomé^d



Microchemical Journal 140 (2018) 158–166



DATUVINK pilot study: A potential non-invasive methodology for dating ballpoint pen inks using multivariate chemometrics based on their UV–vis–NIR reflectance spectra

L. Ortiz-Herrero^{a,*}, L. Bartolomé^b, I. Durán^a, I. Velasco^a, M.L. Alonso^a, M.I. Maguregui^c, M. Ezcurrea^d



Chemometrics and Intelligent Laboratory Systems 207 (2020) 104187



A novel, non-invasive, multi-purpose and comprehensive method to date inks in real handwritten documents based on the monitoring of the dye ageing processes

L. Ortiz-Herrero^{a,*}, A.C. de Almeida Assis^b, L. Bartolomé^c, M.L. Alonso^a, M.I. Maguregui^d, R.M. Alonso^a, J.S. Seixas de Melo^e



Talanta 205 (2019) 120114



OPLS multivariate regression of FTIR-ATR spectra of acrylic paints for age estimation in contemporary artworks

L. Ortiz-Herrero^{a,*}, I. Cardaba^b, S. Setien^a, L. Bartolomé^c, M.L. Alonso^a, M.I. Maguregui^b



Polymer Degradation and Stability 179 (2020) 109263



Extension study of a statistical age prediction model for acrylic paints

L. Ortiz-Herrero^{a,*}, I. Cardaba^b, L. Bartolomé^c, M.L. Alonso^a, M.I. Maguregui^b



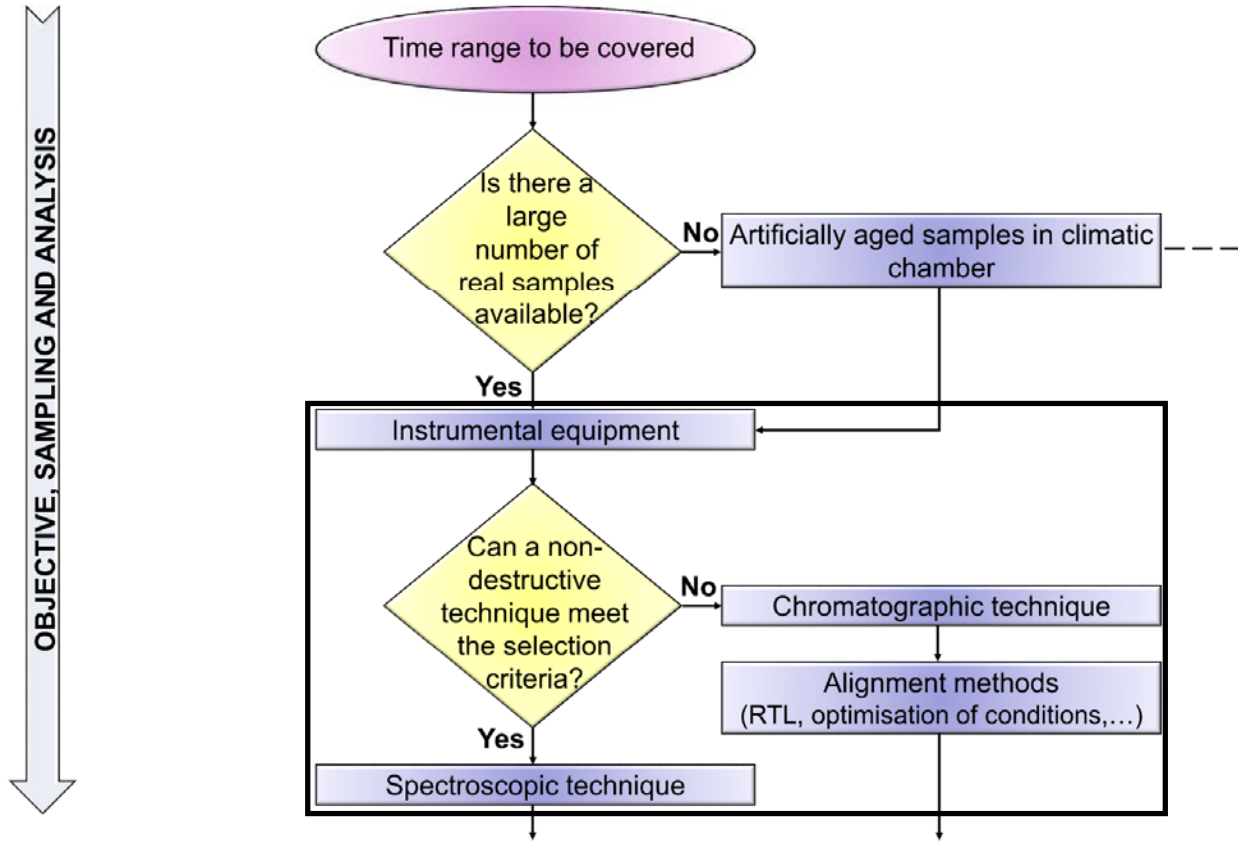
Trends in Analytical Chemistry 141 (2021) 116278



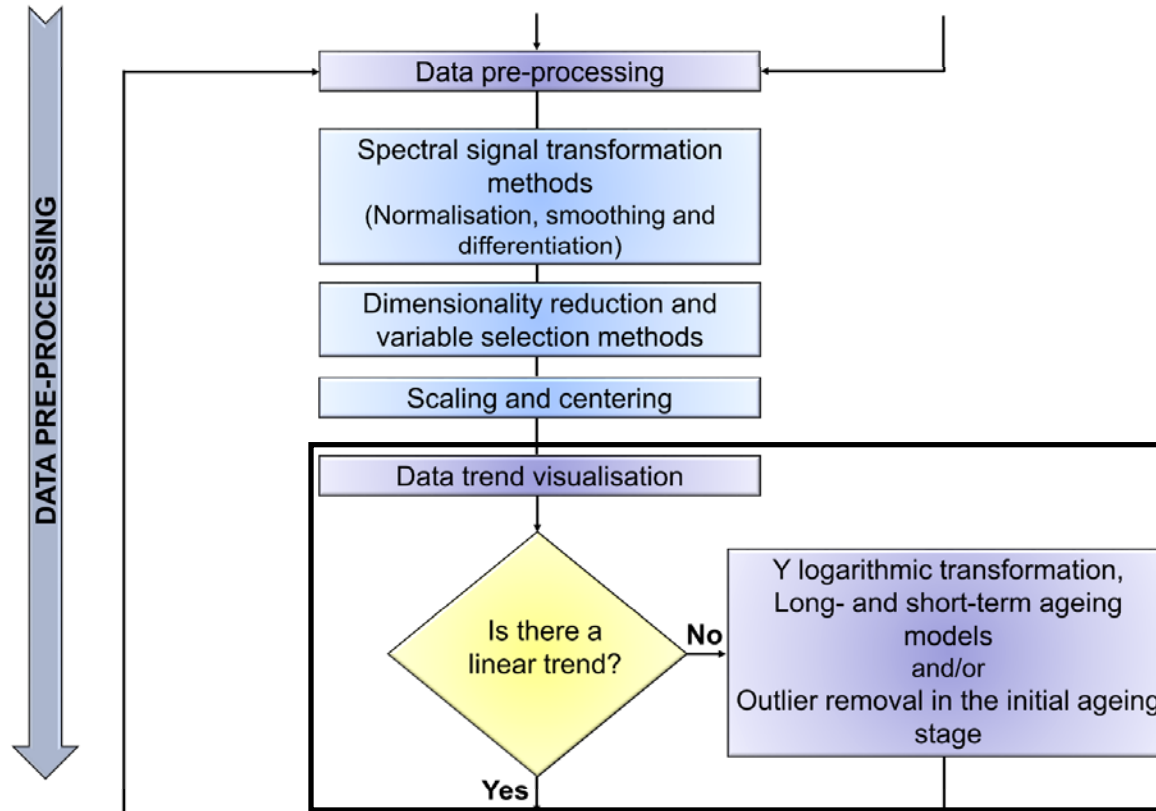
Multivariate (O)PLS regression methods in forensic dating

L. Ortiz-Herrero^{a,*}, M.I. Maguregui^b, L. Bartolomé^c

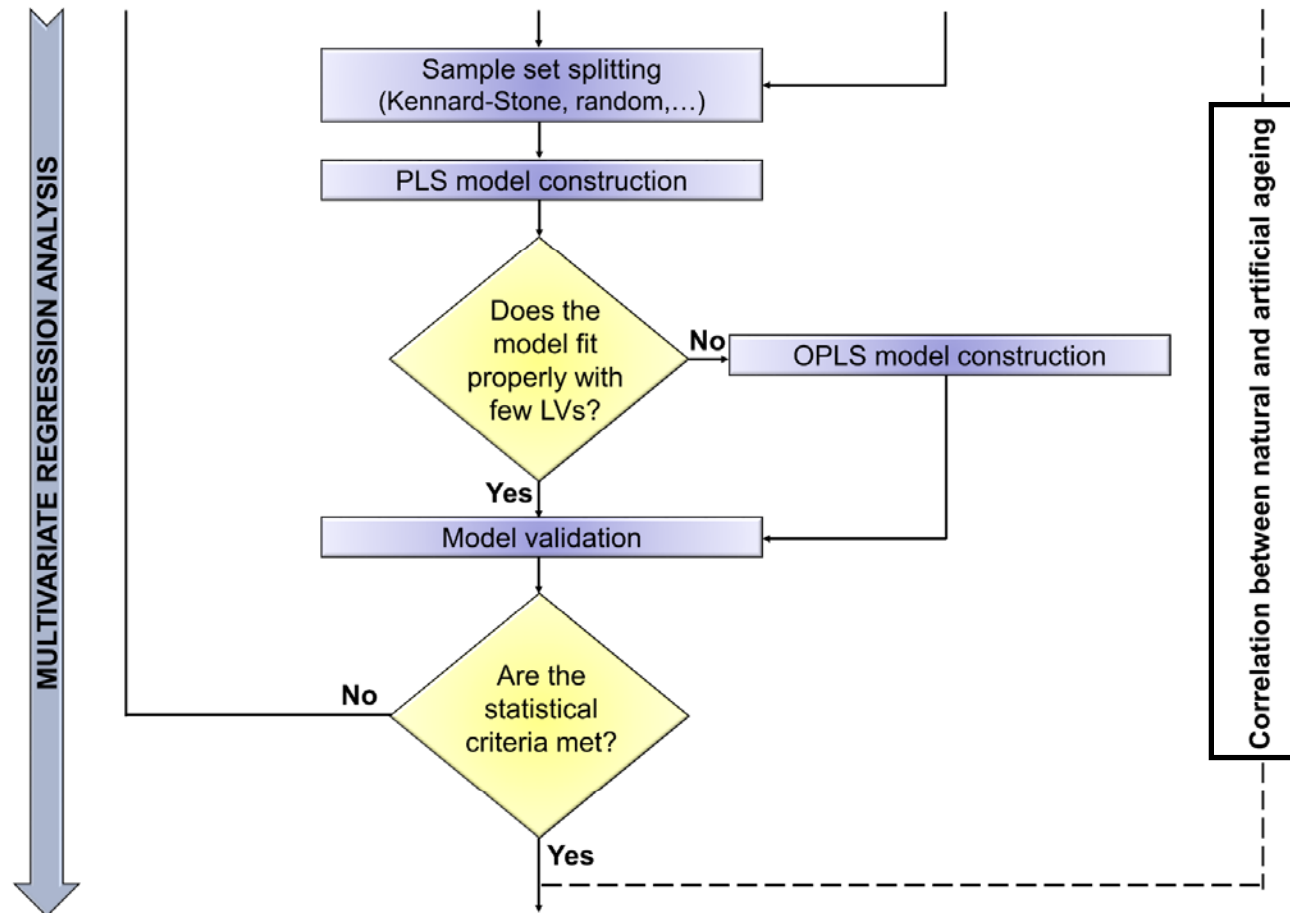


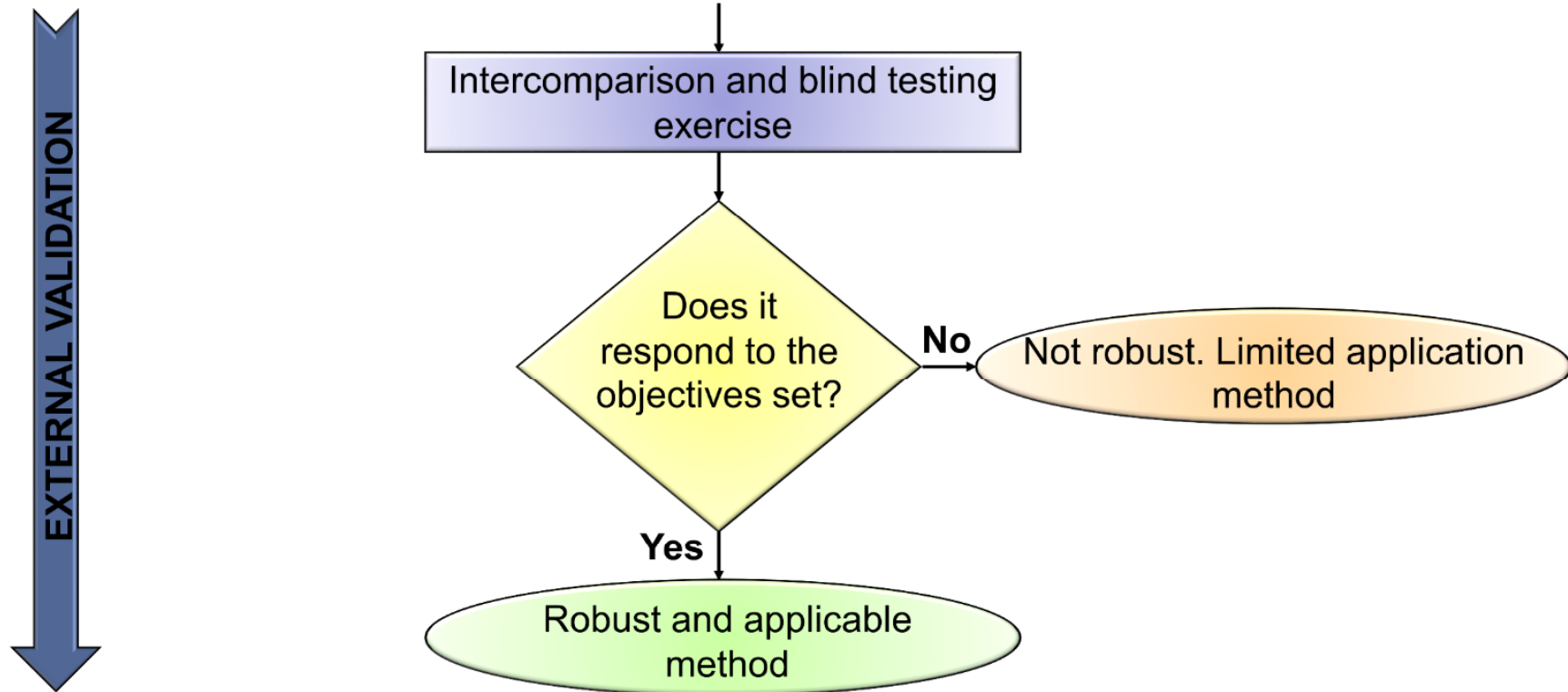


SIMCA 15.0.2 Umetrics® software



SIMCA 15.0.2 Umetrics® software





Contents lists available at [ScienceDirect](https://www.sciencedirect.com)

Chemometrics and Intelligent Laboratory Systems

journal homepage: www.elsevier.com/locate/chemometrics

A novel, non-invasive, multi-purpose and comprehensive method to date inks in real handwritten documents based on the monitoring of the dye ageing processes



L. Ortiz-Herrero^{a,*}, A.C. de Almeida Assis^b, L. Bartolomé^c, M.L. Alonso^a, M.I. Maguregui^d, R.M. Alonso^a, J.S. Seixas de Melo^e

^a Analytical Chemistry Department, Faculty of Science and Technology, University of the Basque Country (UPV/EHU), Barrio Sarriena S/N, 48940, Leioa, Bizkaia, Spain

^b Scientific Police Laboratory, Judiciary Police, Rua Gomes Freire, 174, 1169-007, Lisbon, Portugal

^c Advances Research Facilities (SGIker), Martina Casiano Technology Platform, University of the Basque Country (UPV/EHU), Barrio Sarriena S/N, 48940, Leioa, Bizkaia, Spain

^d Painting Department, Faculty of Fine Arts, University of the Basque Country (UPV/EHU), Barrio Sarriena S/N, 48940, Leioa, Bizkaia, Spain

^e Coimbra Chemistry Centre, Department of Chemistry, University of Coimbra, Rua Larga, 3004-535, Coimbra, Portugal



Requirements to be met by the **methodologies** to be developed:



Drawbacks encountered in the development of these methodologies:



Objective

- ✓ Method for **dating handwritten documents**.
- ✓ **Accurate** prediction of the ink stroke.
- ✓ Applied to the **largest number** of pen **inks** and to a **wide time range**.
- ✓ **Non-invasive** measurement by Vis-MSP.
- ✓ **Reliable** results with 95% confidence.
- ✓ **Minimal influence** of paper **support** and preservation **conditions**.

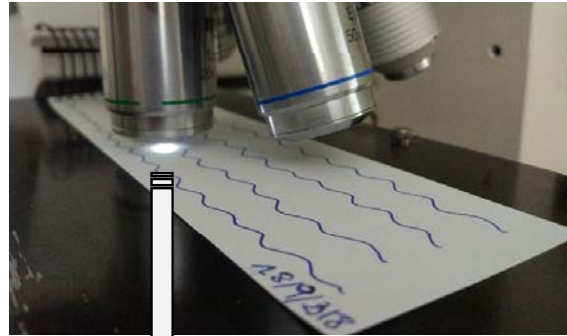
Ink sampling

Pen brand	Pen type	No. of samples	Temporary range of application (mo.)
Bic® Cristal Medium France	Ballpoint	25	25
Bic® Cristal Medium USA	Ballpoint	30	27
Staedtler® Stick 430 M	Ballpoint	27	23
Paper Mate® Flex Grip Elite 1.4 M	Ballpoint	20	25
Paper Mate® 0.7mm	Liquid (gel ink)	20	25
Pilot® Super Grip M	Ballpoint	25	25
Pilot® G-1	Liquid (gel ink)	20	25
Faber-Castell® Medium	Ballpoint	28	23
Uni-Ball® Signo Broad	Liquid (gel ink)	23	25
Uni-Ball® Jetstream Sport	Ballpoint	21	25
Inoxcrom® TC Ball Stainless Steel Tip M	Liquid (gel ink)	22	25

11 writing tools from **7 brands**

Max. **2.5 natural years**

Vis-MSP analysis



Spectrophotometer J&M TIDAS® MSP-800

Interpolation yes

1 nm resolution

Single scan, 3 accumulations

1 pixel bunching

400-800 nm range, reflectance acquisition mode

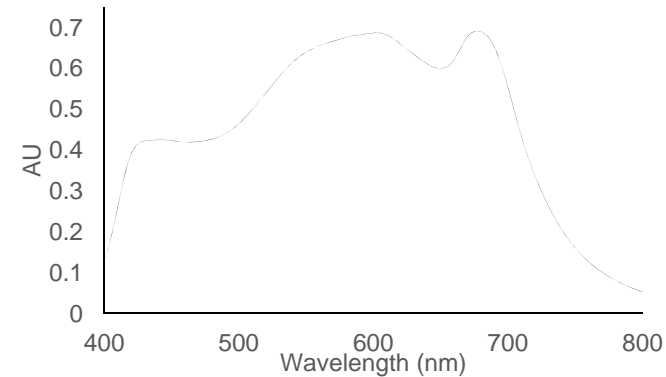
Zeiss® Axiotech 100 Microscope

220.0 x 127.0 μm diaphragm

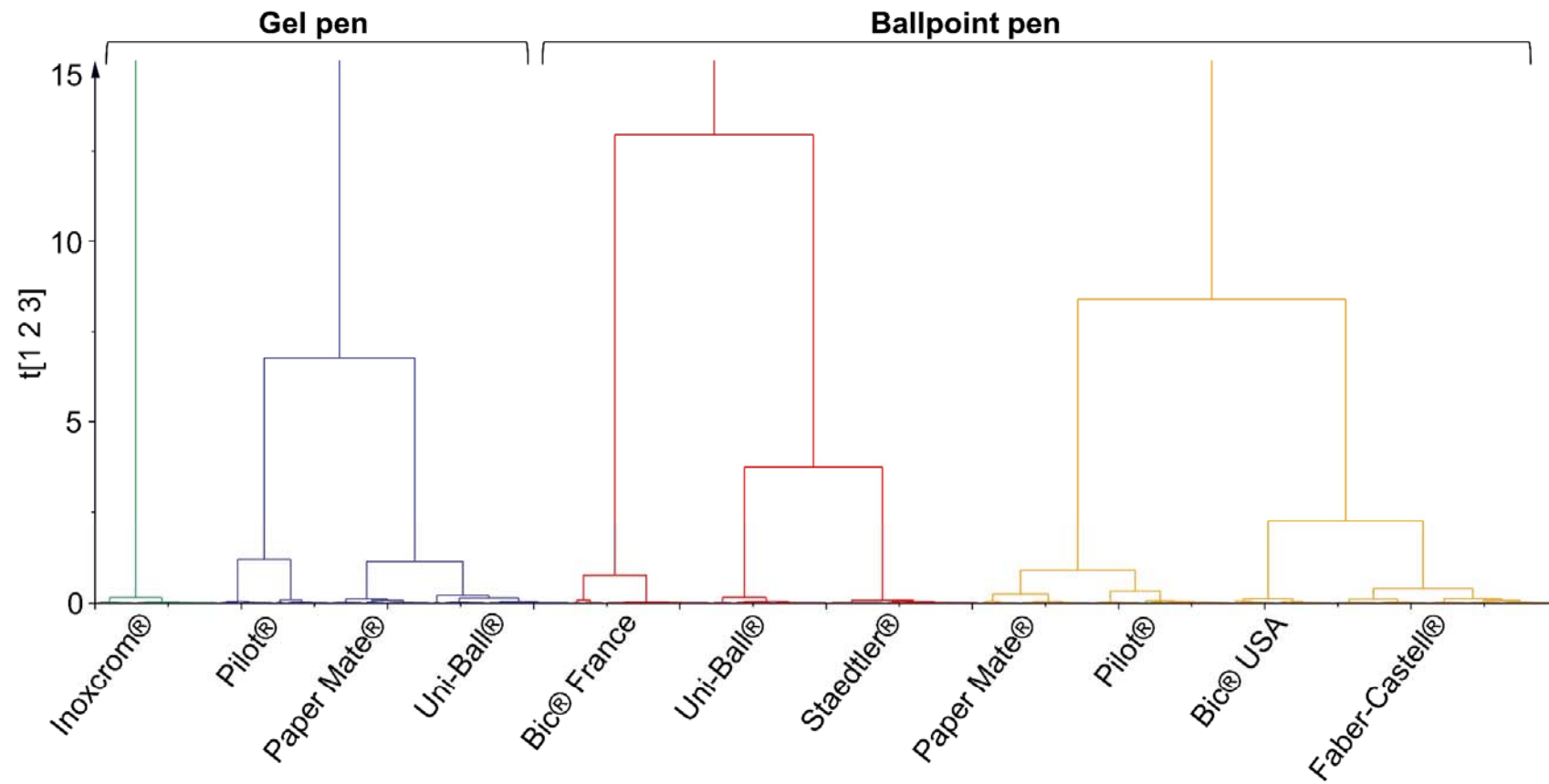
640 x 80 image resolution

20x objective magnification

10x light intensity



Pre-classification by HCA

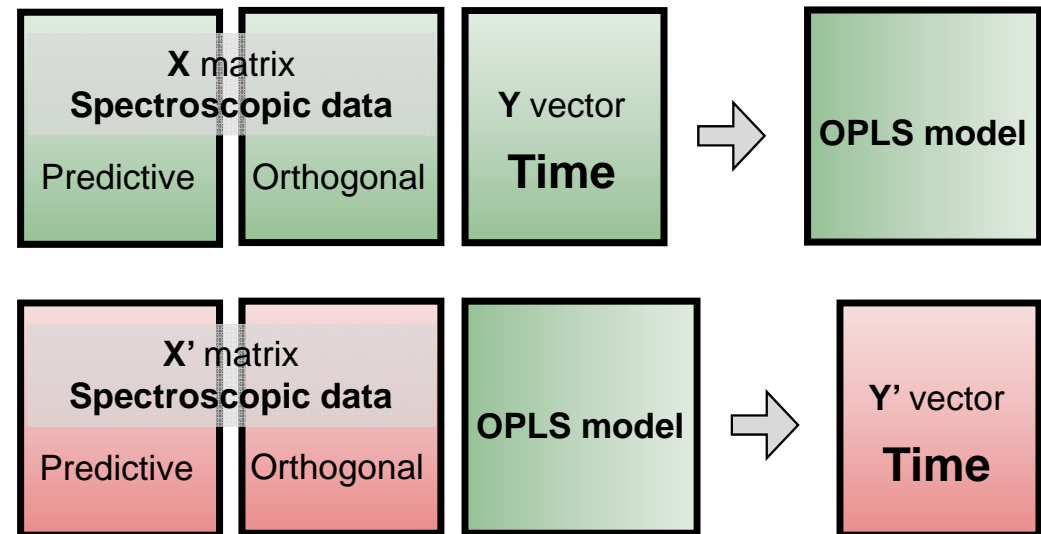


OPLS model building and validation

Ink type	Pen brand
Ballpoint	Bic® France
	Bic® USA
	Faber-Castell®
	Staedtler®
	Pilot®
	Paper Mate®
	Uni-Ball®
	Paper Mate®
	Uni-Ball®
	Pilot®
Gel	Inoxcrom®

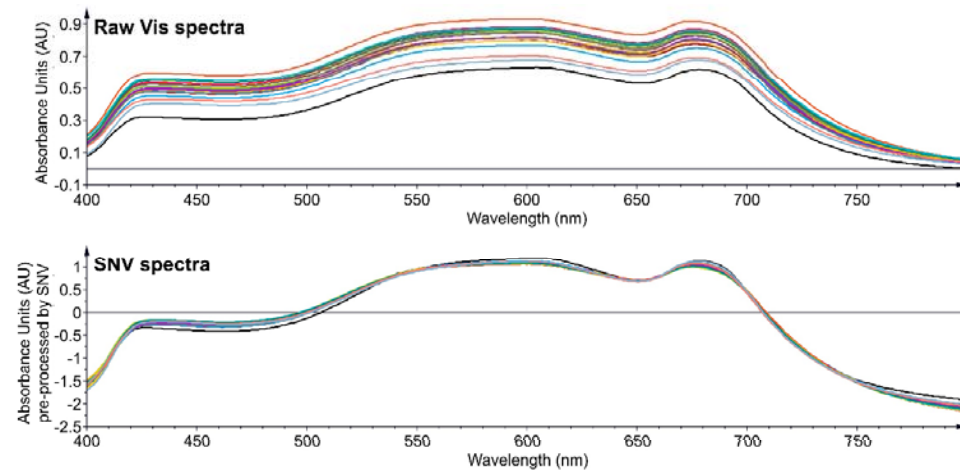
SIMCA V. 15.0.2 Umetrics® software Stone algorithm, 80% training and 20% test

One OPLS model per writer (n=31) Rapidminer® Studio V. 9.8 software



OPLS model building and validation

Ink type	Pen brand	No. calibration/validation samples	Pre-treatment (scale/filter)
Ballpoint	Bic® France	20/5	UV and SNV
	Bic® USA	23/7	Centring and SNV
	Faber-Castell®	21/7	UV and MSC
	Staedtler®	20/7	UV and 1st derivative
	Pilot®	20/5	UV and SNV
	Paper Mate®	15/5	UV and 1st derivative
	Uni-Ball®	16/5	UV and MSC
Gel	Paper Mate®	15/5	UV and SNV
	Uni-Ball®	18/5	UV and 1st derivative
	Pilot®	15/5	UV and MSC
	Inoxcrom®	17/5	UV and MSC



Remove systematic **non-time-related** variation in X

OPLS model building and validation

Model fitting Internal predictive testing
by cross-validation

Ink type	Pen brand	No. calibration/validation samples	Pre-treatment (scale/filter)	LV	R ² Y	Q ²	RMSEE (days)	RMSECV (days)
Ballpoint	Bic® France	20/5	UV and SNV	4	0.87	0.74	1.2	1.3
	Bic® USA	23/7	Centring and SNV	4	0.86	0.74	1.2	1.3
	Faber-Castell®	21/7	UV and MSC	2	0.71	0.52	1.4	1.4
	Staedtler®	20/7	UV and 1st derivative	6	0.95	0.83	1.2	1.3
	Pilot®	20/5	UV and SNV	1	0.65	0.55	1.4	1.4
	Paper Mate®	15/5	UV and 1st derivative	3	0.78	0.53	1.4	1.5
	Uni-Ball®	16/5	UV and MSC	3	0.82	0.63	1.3	1.3
Gel	Paper Mate®	15/5	UV and SNV	3	0.70	0.41	1.2	1.3
	Uni-Ball®	18/5	UV and 1st derivative	2	0.58	0.41	1.5	1.5
	Pilot®	15/5	UV and MSC	1	0.60	0.45	1.2	1.3
	Inoxcrom®	17/5	UV and MSC	3	0.50	0.31	1.4	1.4

OPLS model building and validation

$$E\% = \left(\frac{y_{real} - y_{predicted}}{y_{real}} \right) \times 100$$

External predictive testing



Ink type	Pen brand	No. calibration/validation samples	Pre-treatment (scale/filter)	LV	R ² Y	Q ²	RMSEE (days)	RMSECV (days)	RMSEP (days)	Accuracy error (%)
Ballpoint	Bic® France	20/5	UV and SNV	4	0.87	0.74	1.2	1.3	1.1	10
	Bic® USA	23/7	Centring and SNV	4	0.86	0.74	1.2	1.3	1.1	10
	Faber-Castell®	21/7	UV and MSC	2	0.71	0.52	1.4	1.4	1.2	15
	Staedtler®	20/7	UV and 1st derivative	6	0.95	0.83	1.2	1.3	1.2	15
	Pilot®	20/5	UV and SNV	1	0.65	0.55	1.4	1.4	1.2	12
	Paper Mate®	15/5	UV and 1st derivative	3	0.78	0.53	1.4	1.5	1.3	18
	Uni-Ball®	16/5	UV and MSC	3	0.82	0.63	1.3	1.3	1.3	22
Gel	Paper Mate®	15/5	UV and SNV	3	0.70	0.41	1.2	1.3	1.3	13
	Uni-Ball®	18/5	UV and 1st derivative	2	0.58	0.41	1.5	1.5	1.4	25
	Pilot®	15/5	UV and MSC	1	0.60	0.45	1.2	1.3	1.3	16
	Inoxcrom®	17/5	UV and MSC	3	0.50	0.31	1.4	1.4	1.2	15

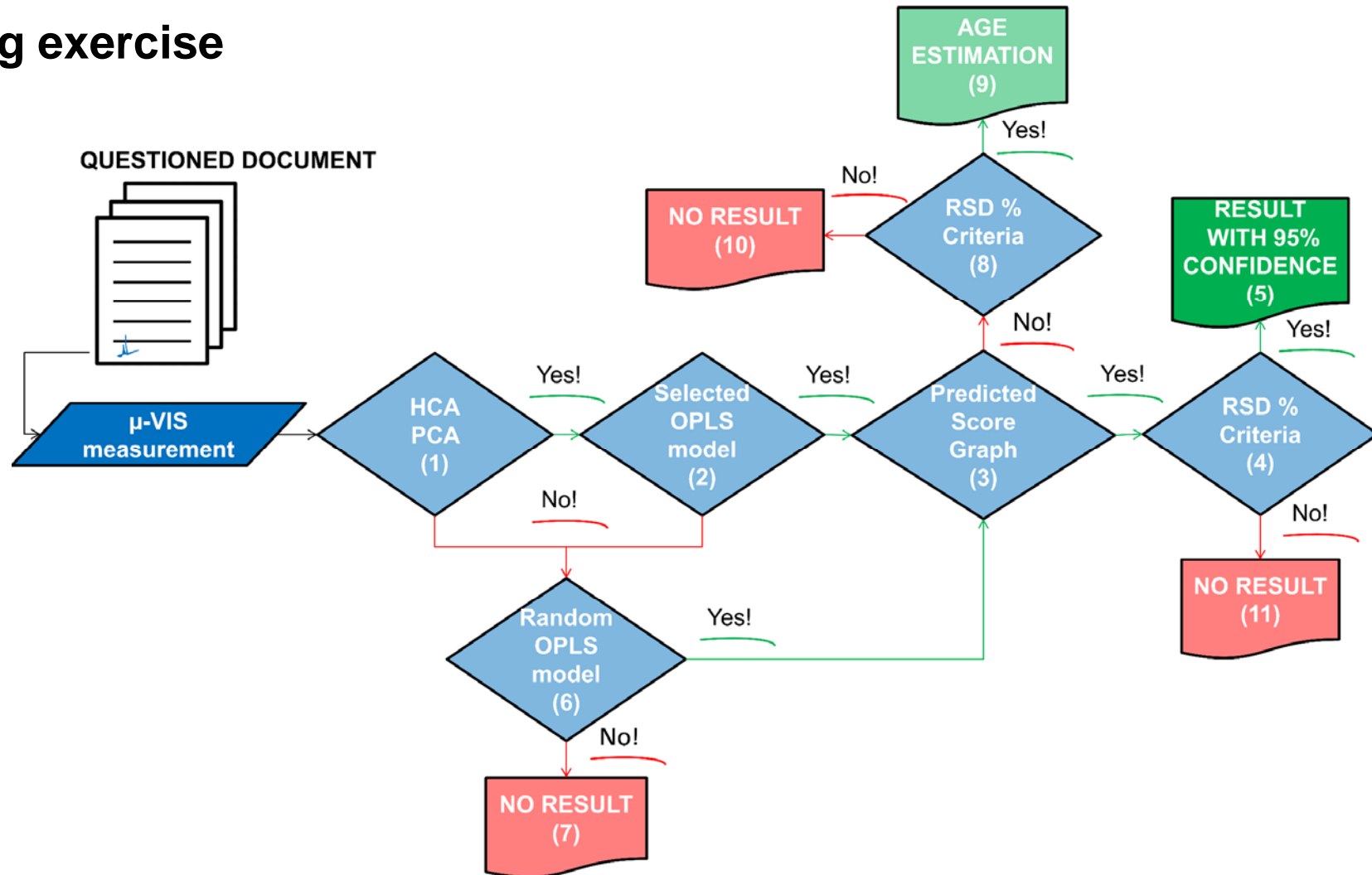
Blind testing exercise

SDG-Q7
F12018

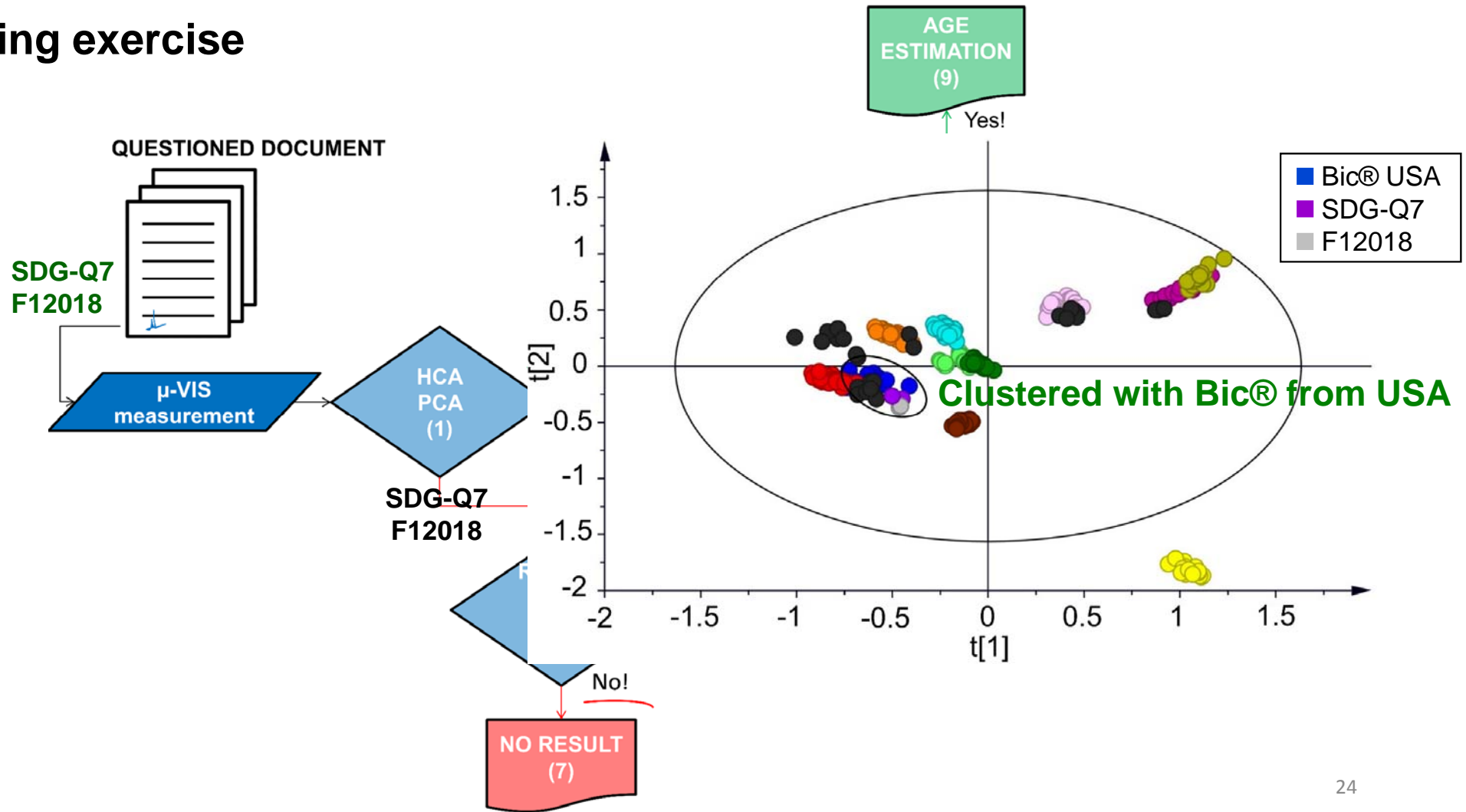
F22018
F32018
F42018

SDG-Q6
SDG-Q8

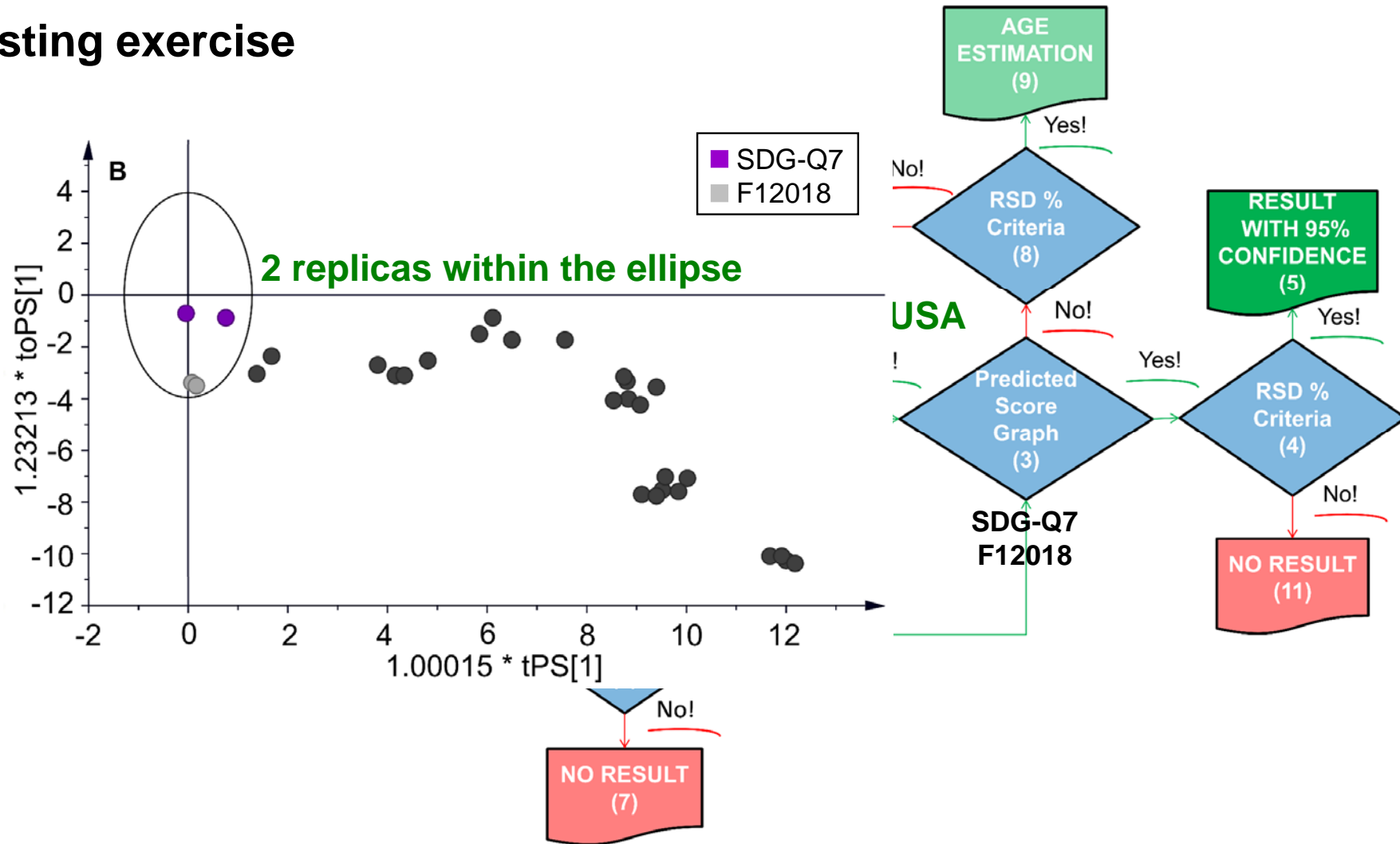
F12019
F22019
F12017
F32017
F42017
F52017



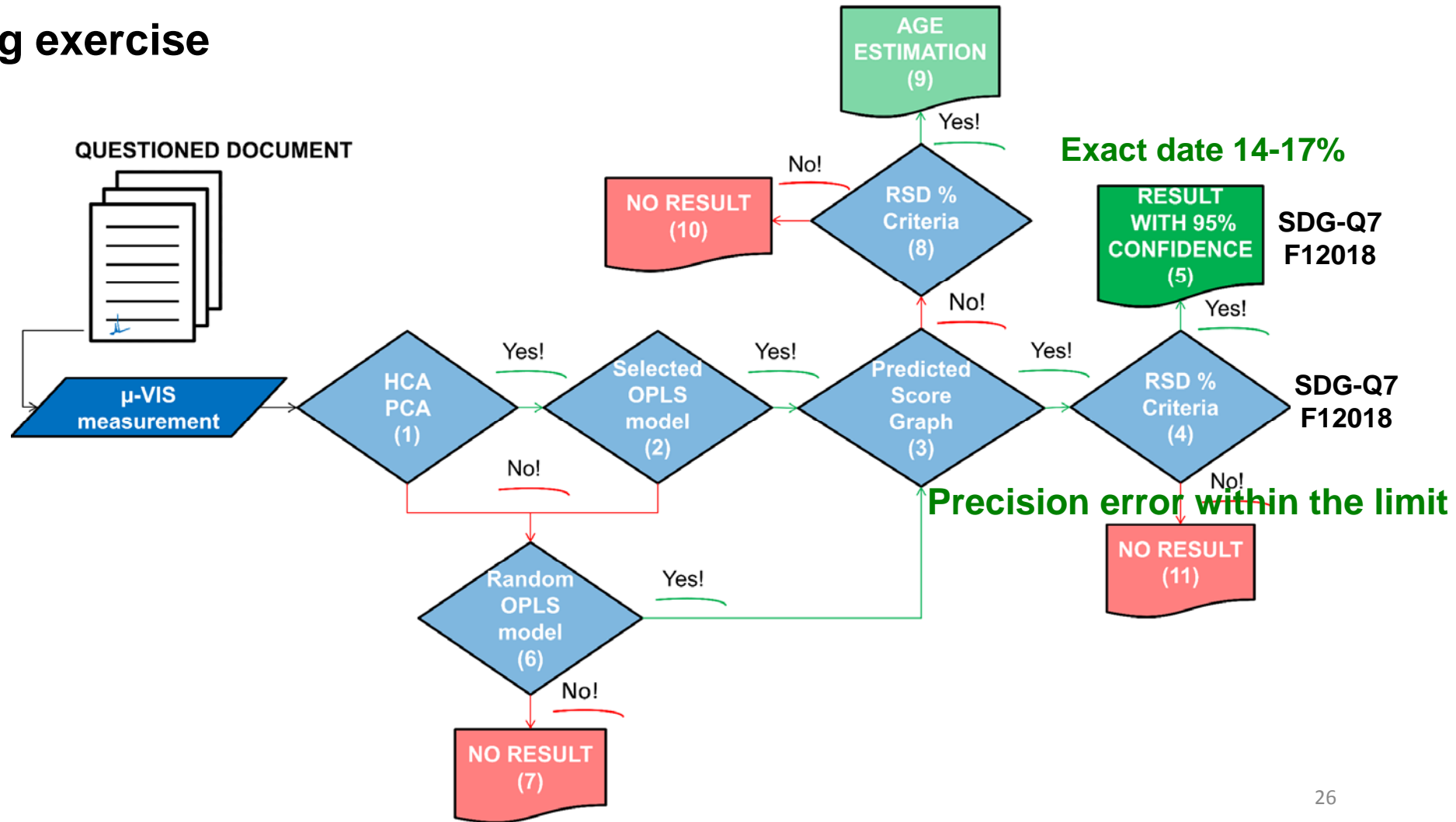
Blind testing exercise



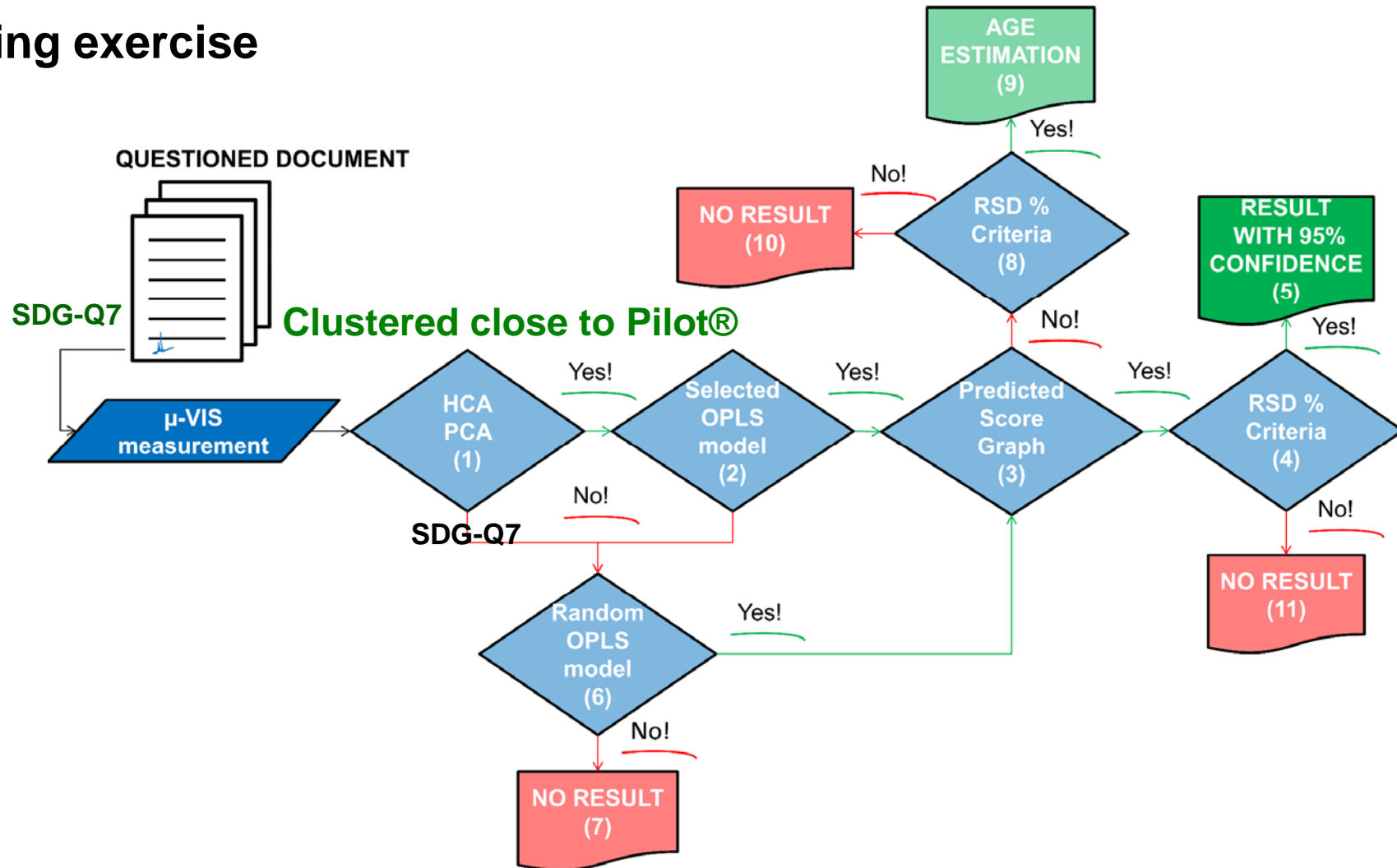
Blind testing exercise



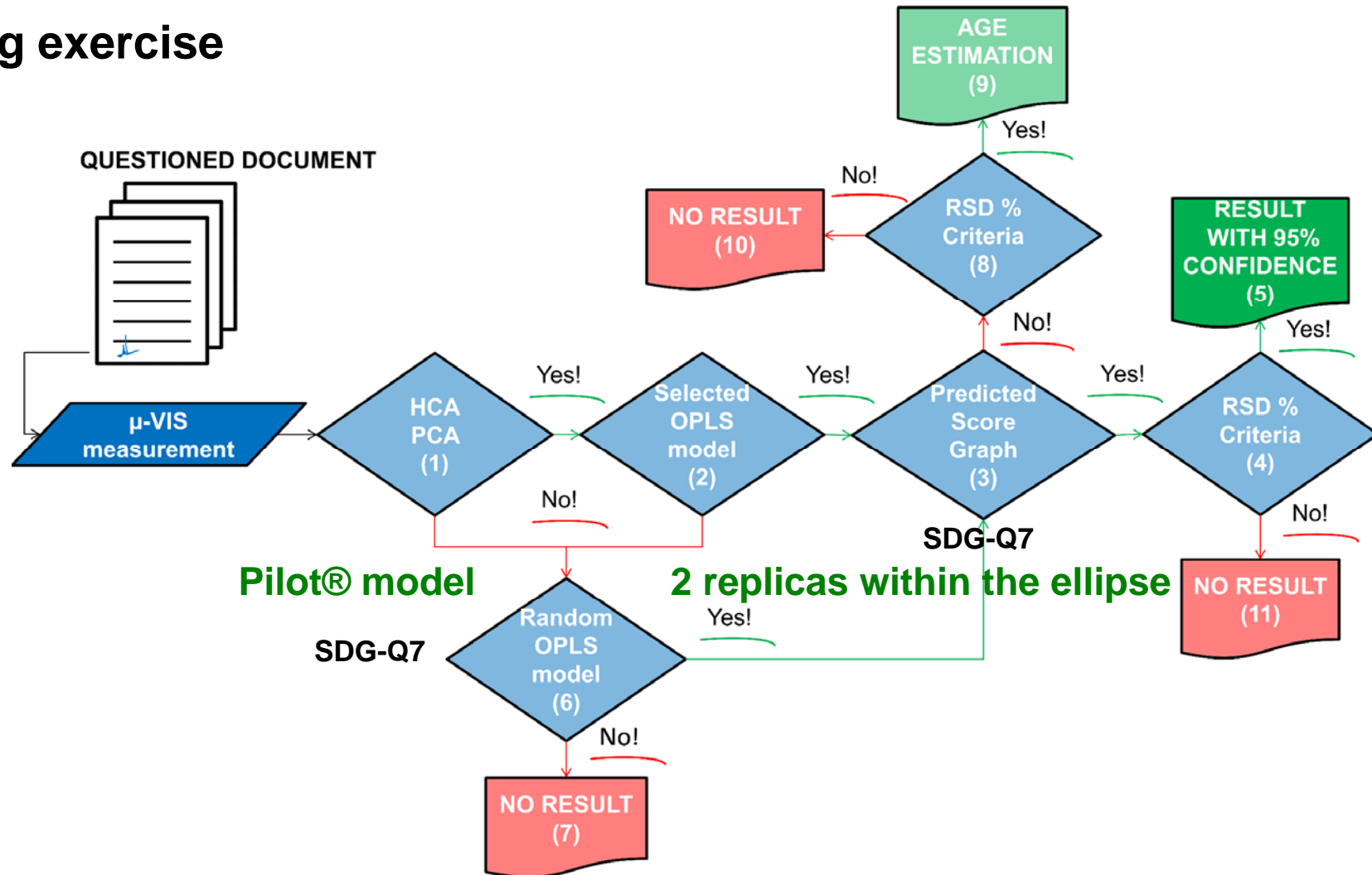
Blind testing exercise



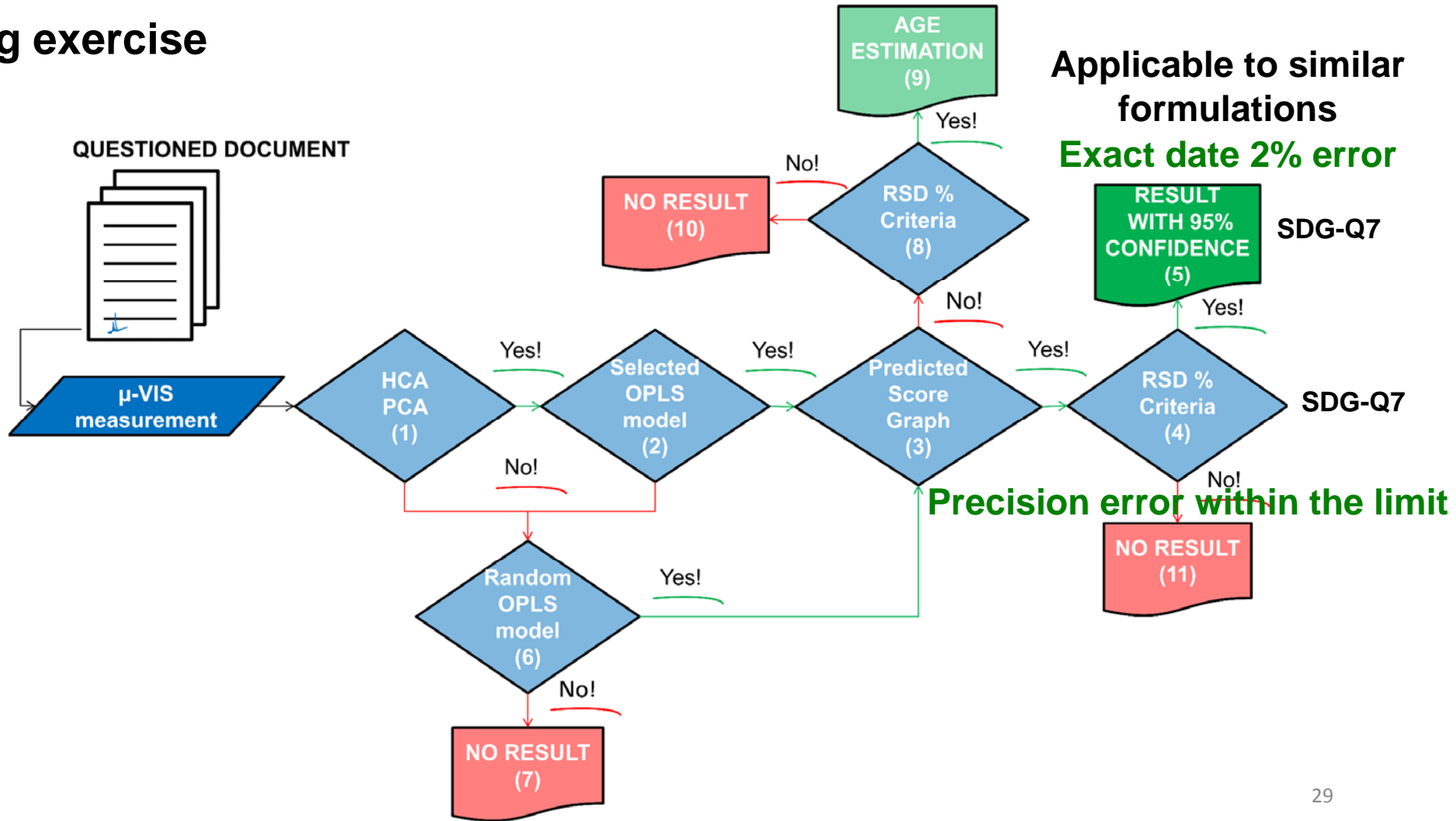
Blind testing exercise



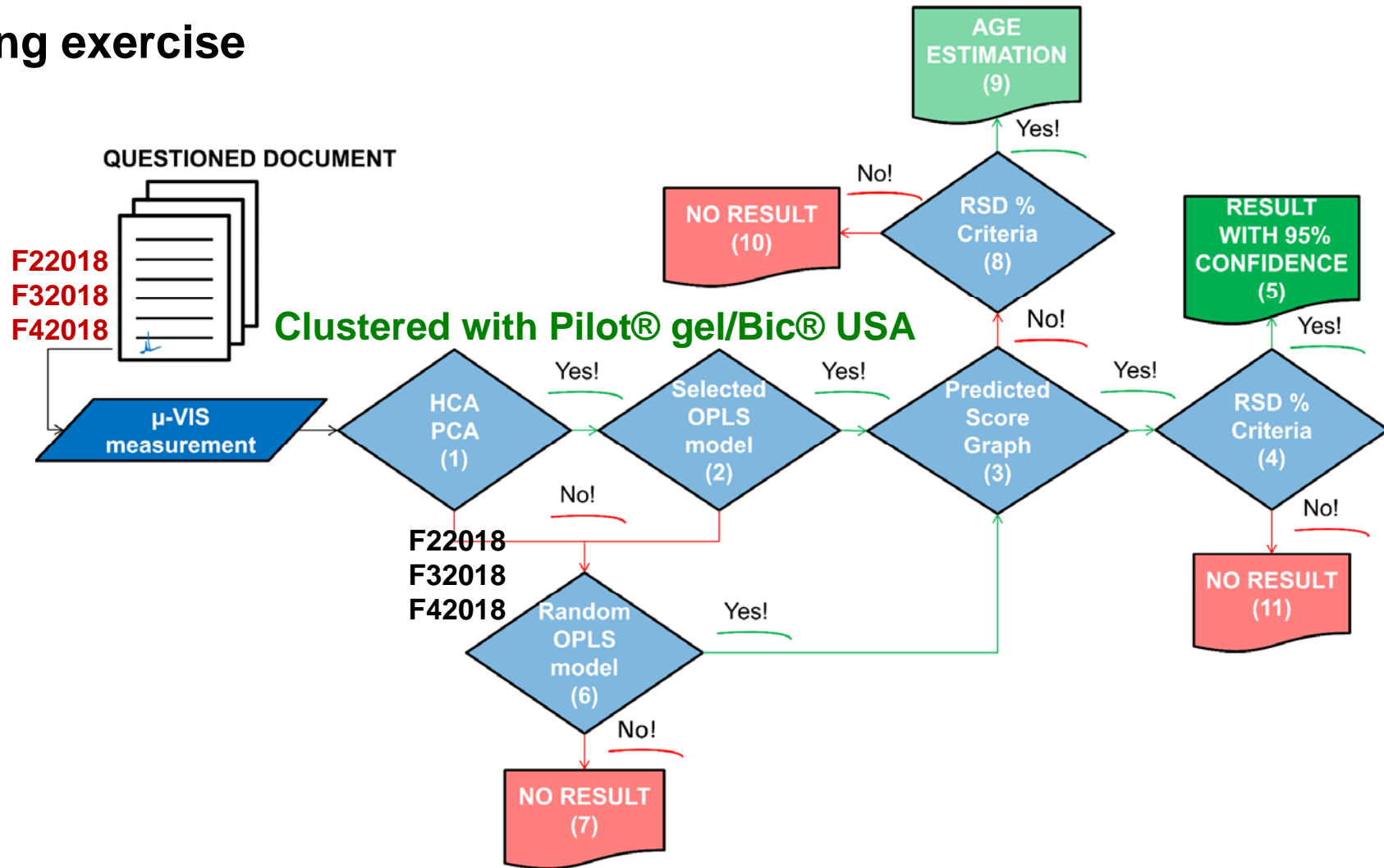
Blind testing exercise



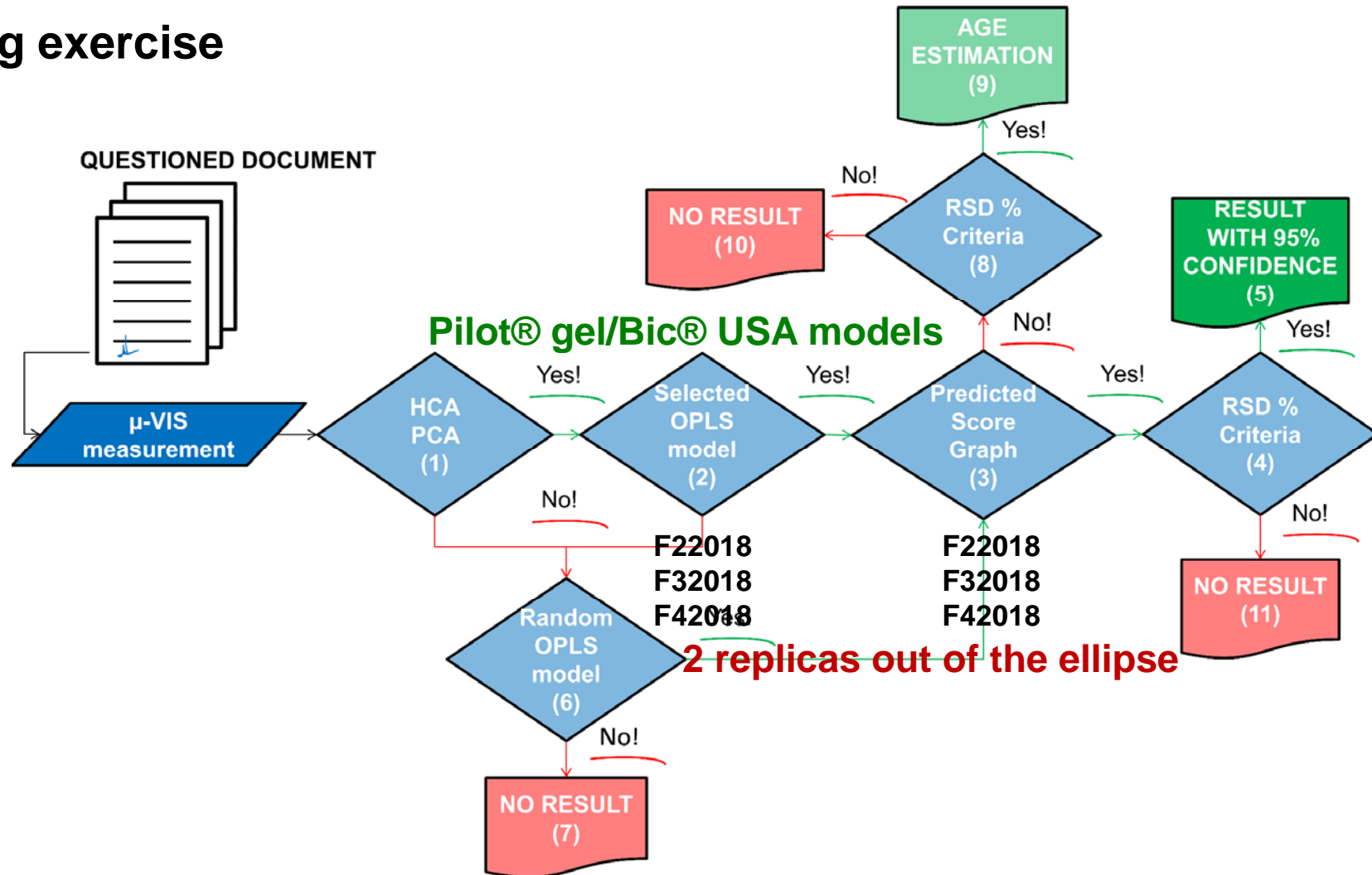
Blind testing exercise



Blind testing exercise

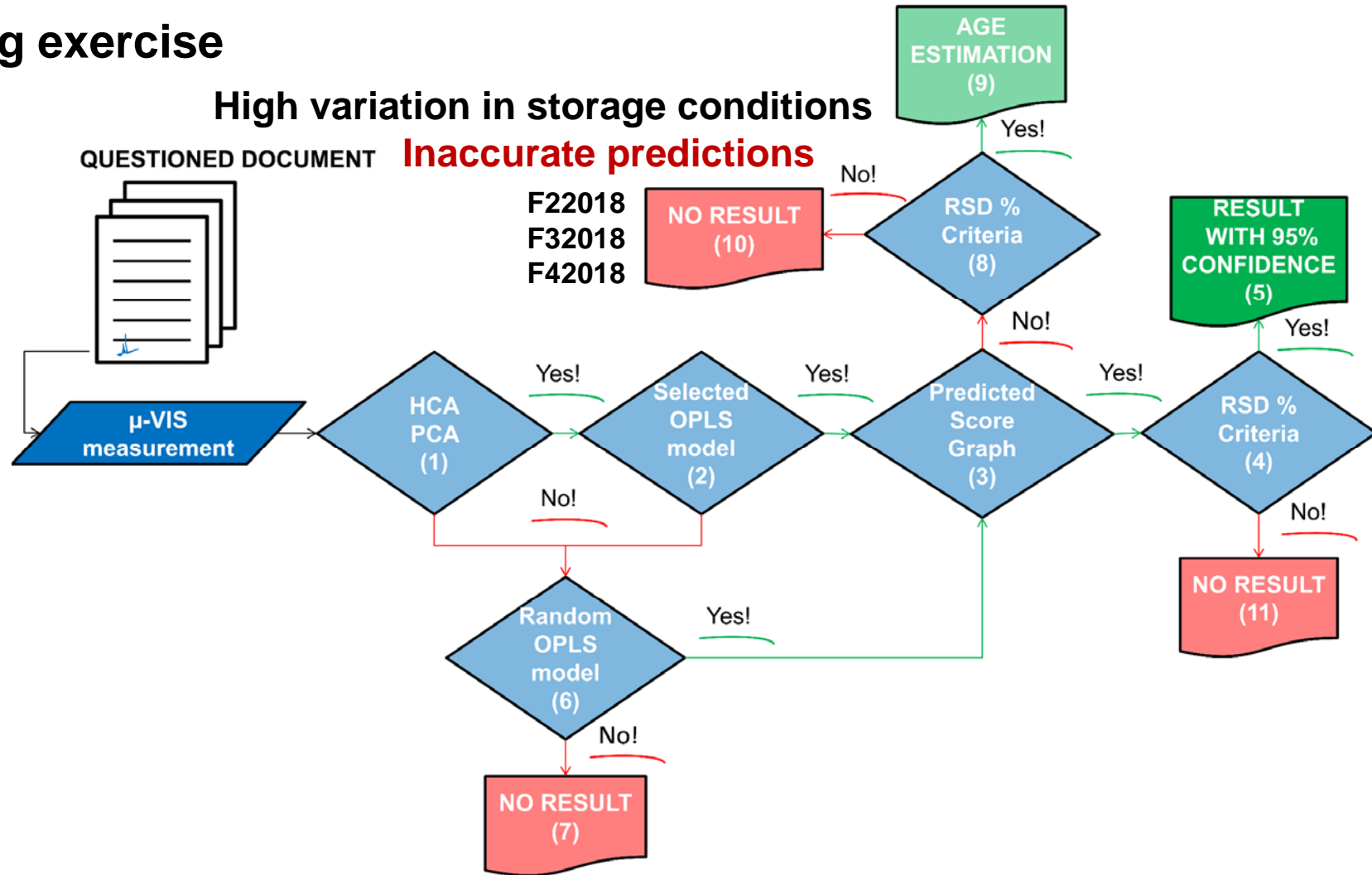


Blind testing exercise

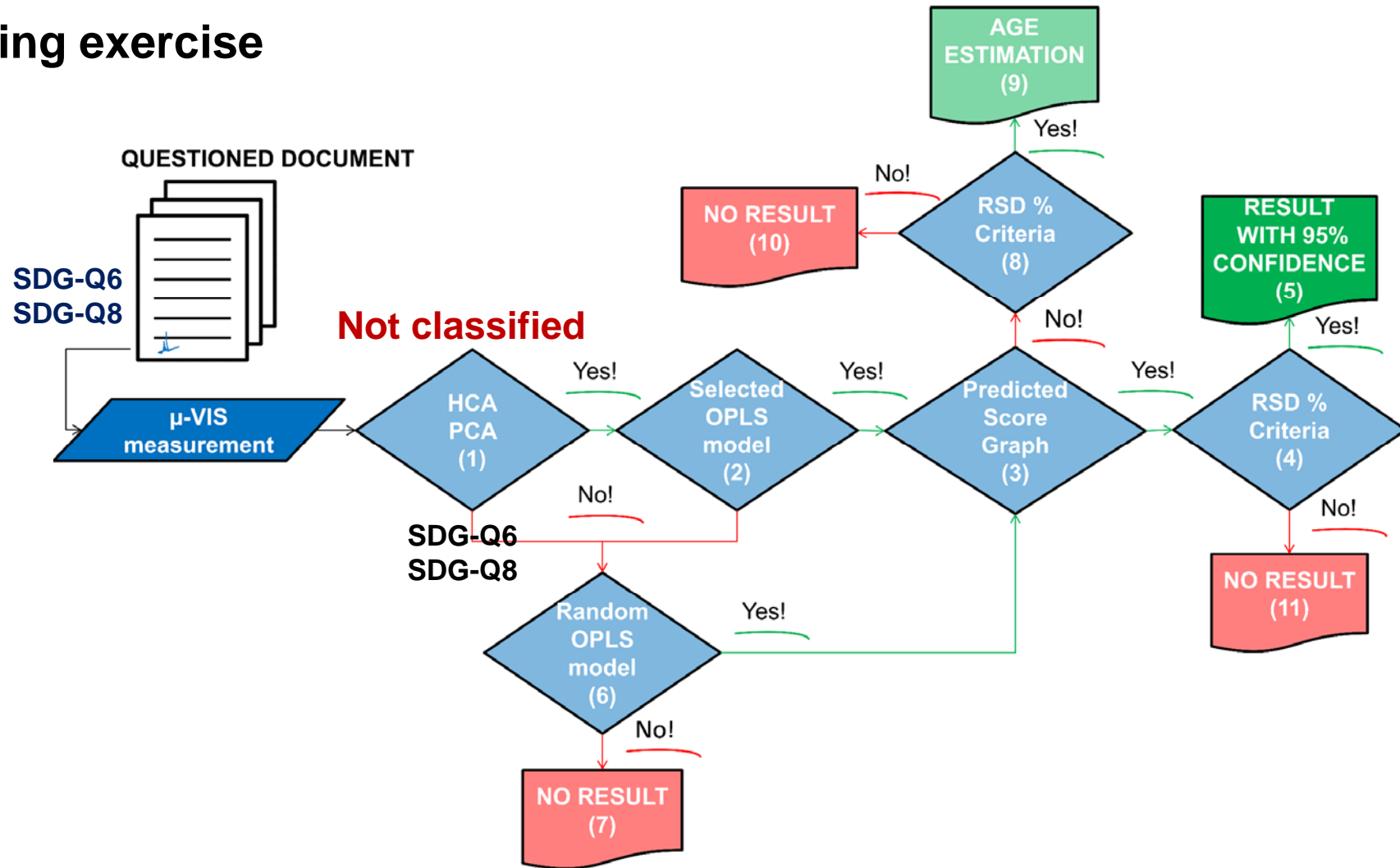


Blind testing exercise

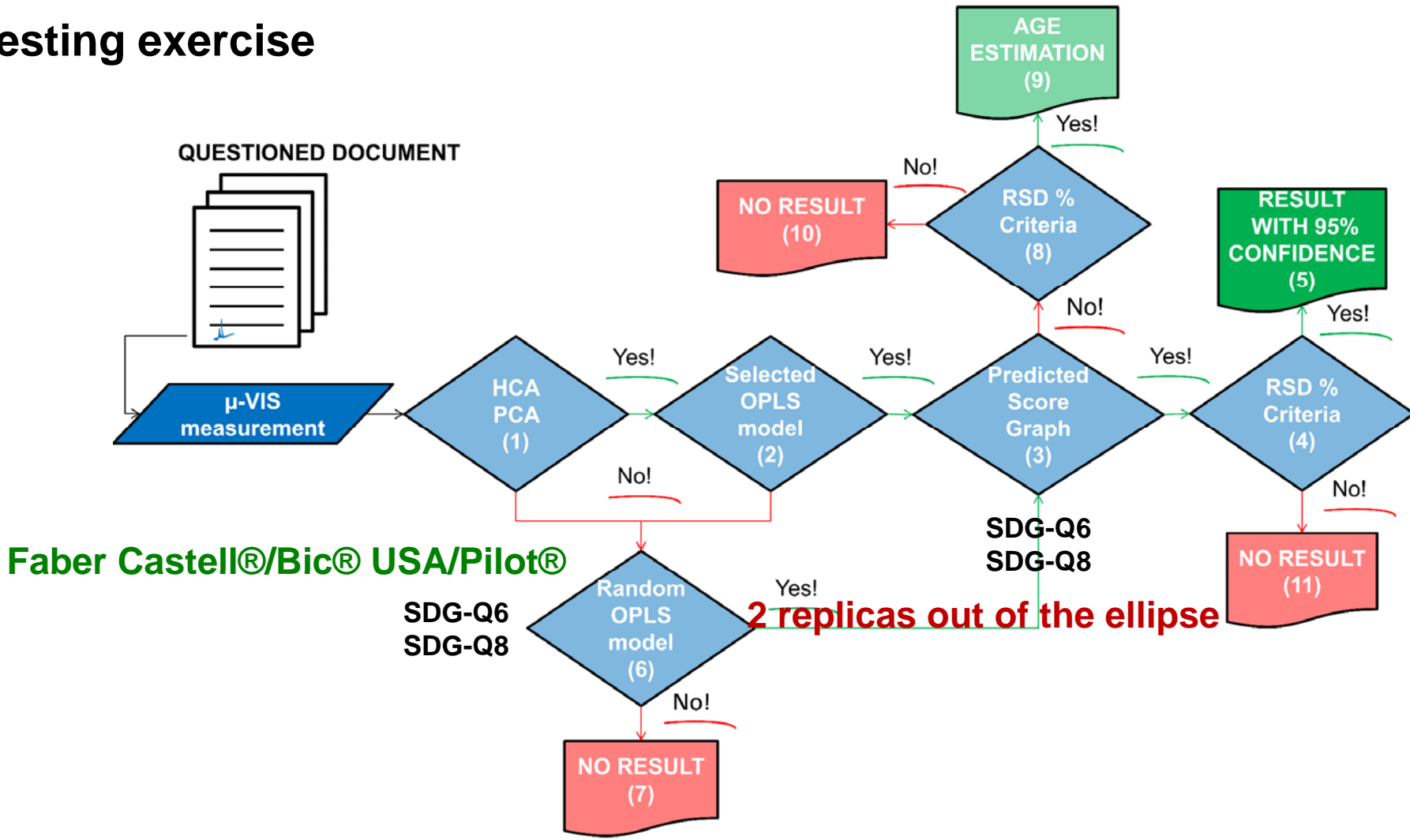
High variation in storage conditions



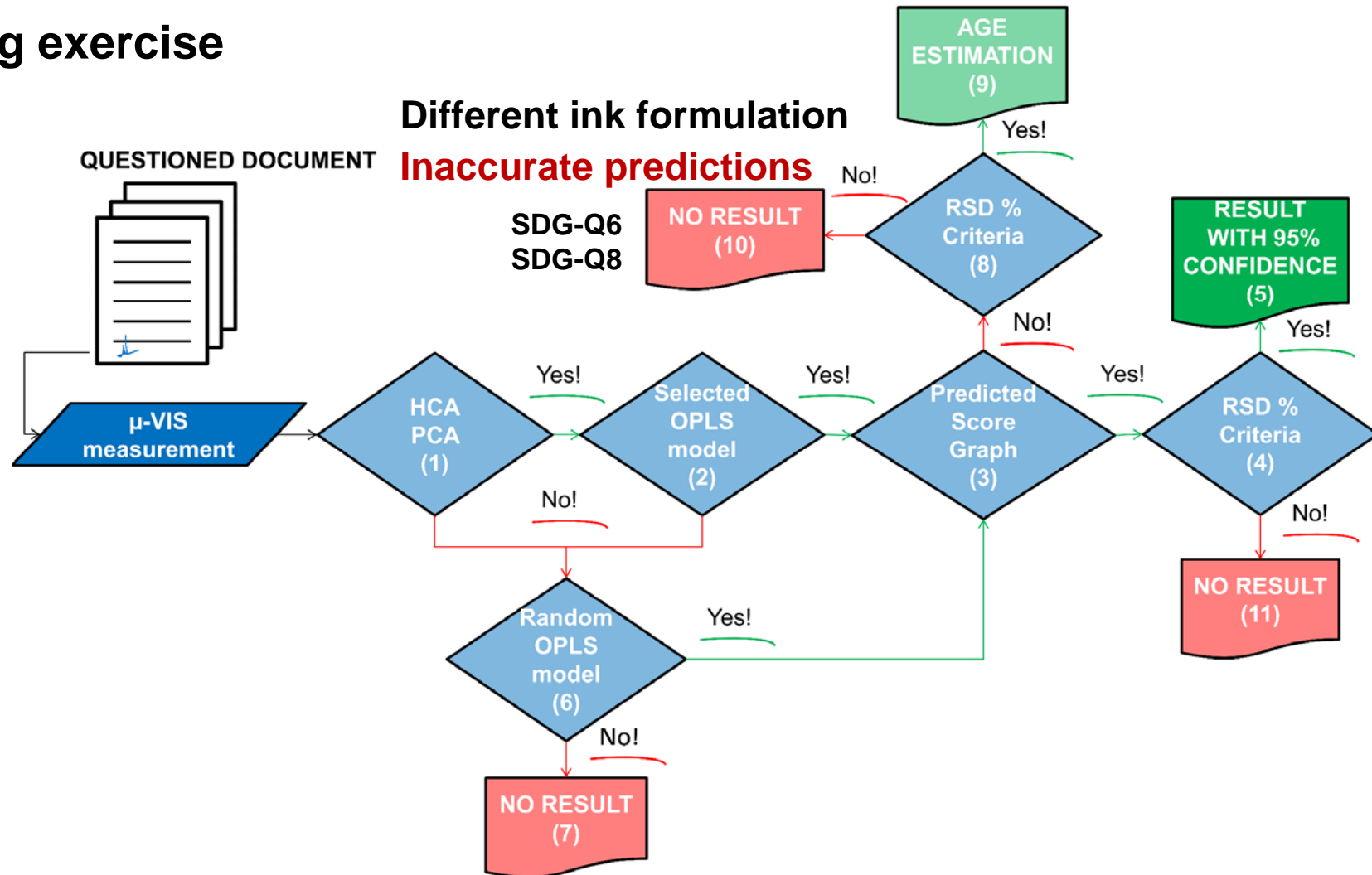
Blind testing exercise



Blind testing exercise



Blind testing exercise



Remarks

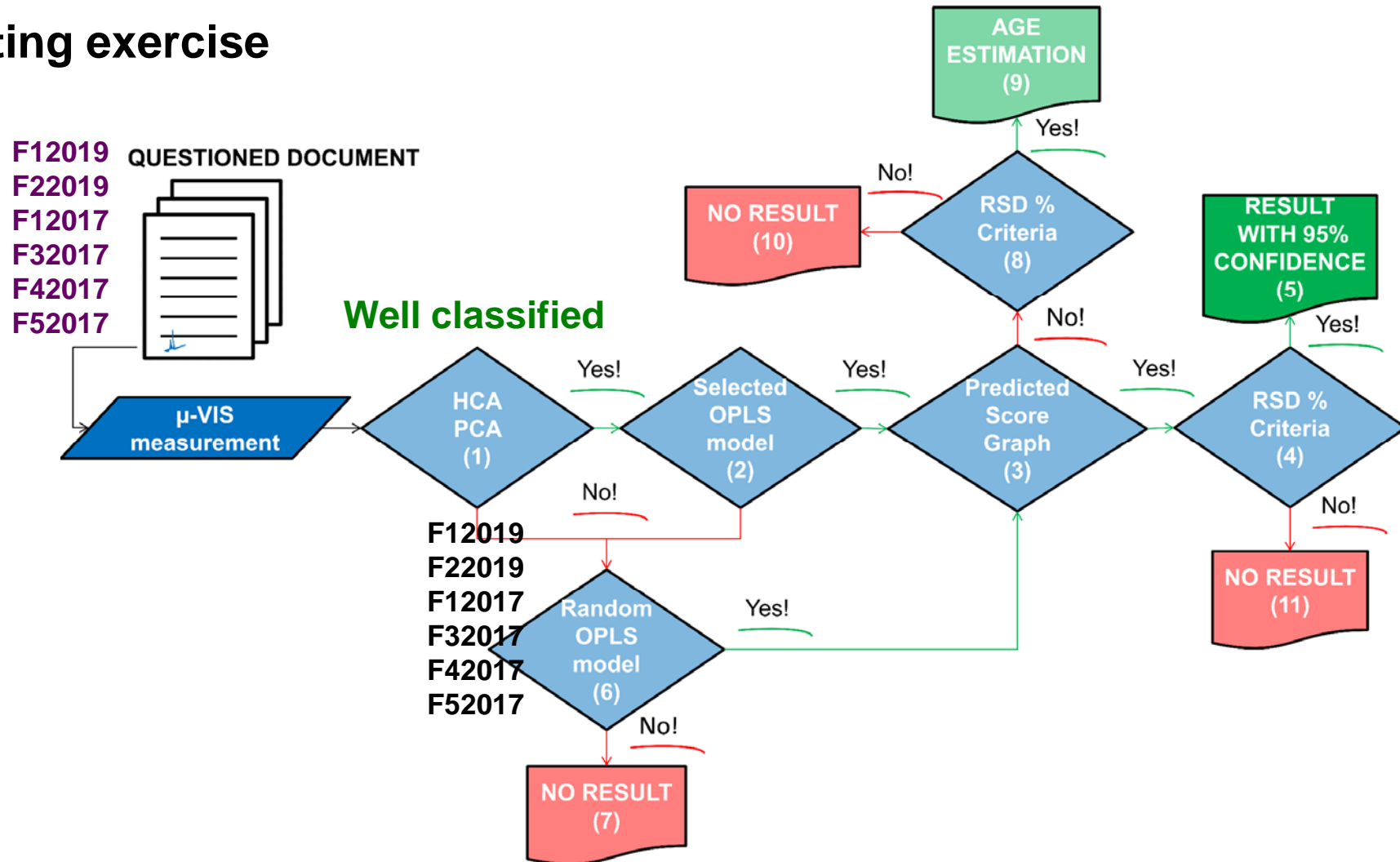
Exact date prediction with **95% confidence** whenever:

- I. **Both replicas** lie **within the ellipse** of the predicted score plot.
- II. The **age** of the ink is **within the time application range** of the corresponding OPLS model.
- III. The ink is **well classified/clustered** into one of the classes/groups of pen brands studied.

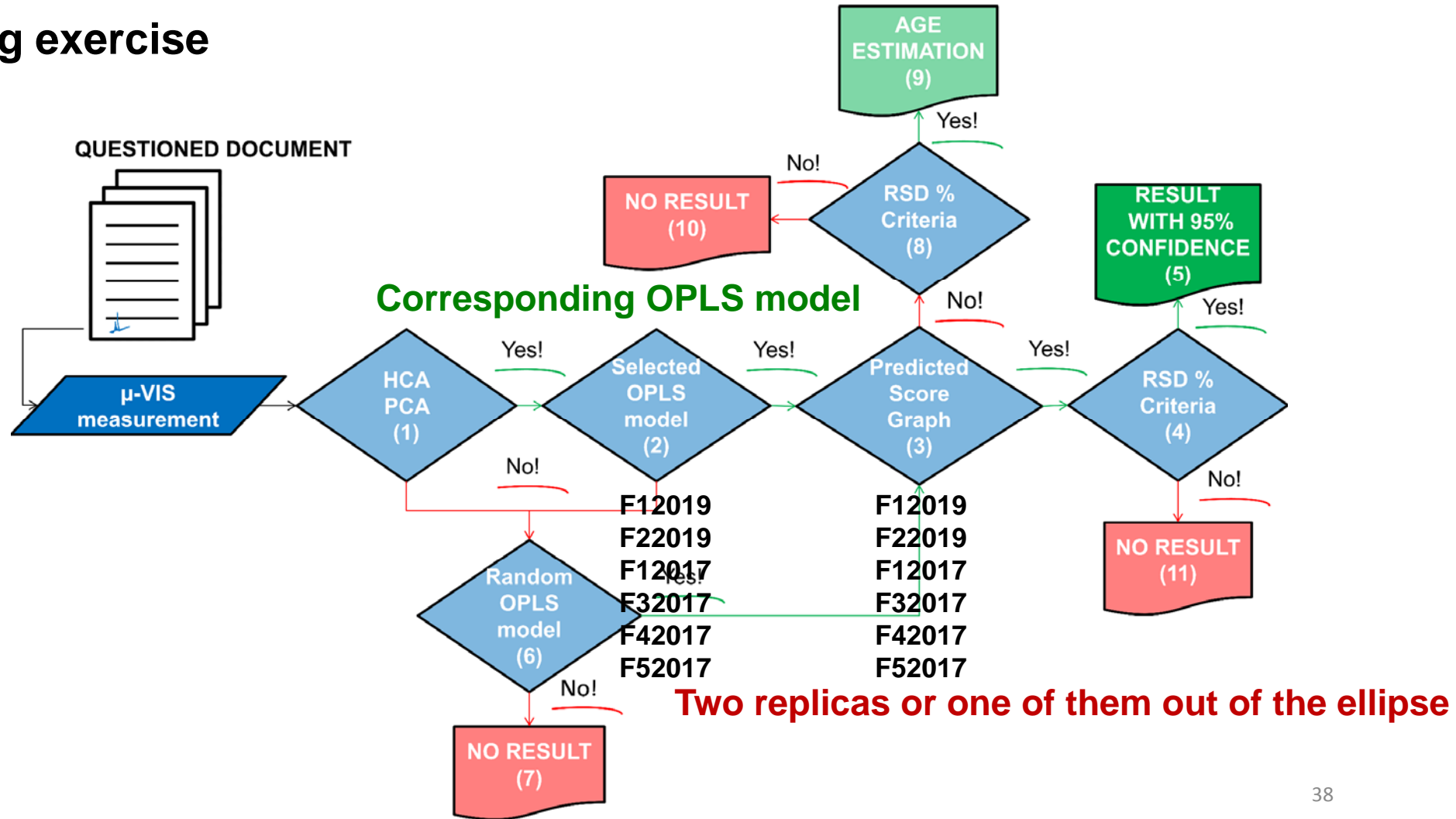
\approx



Blind testing exercise

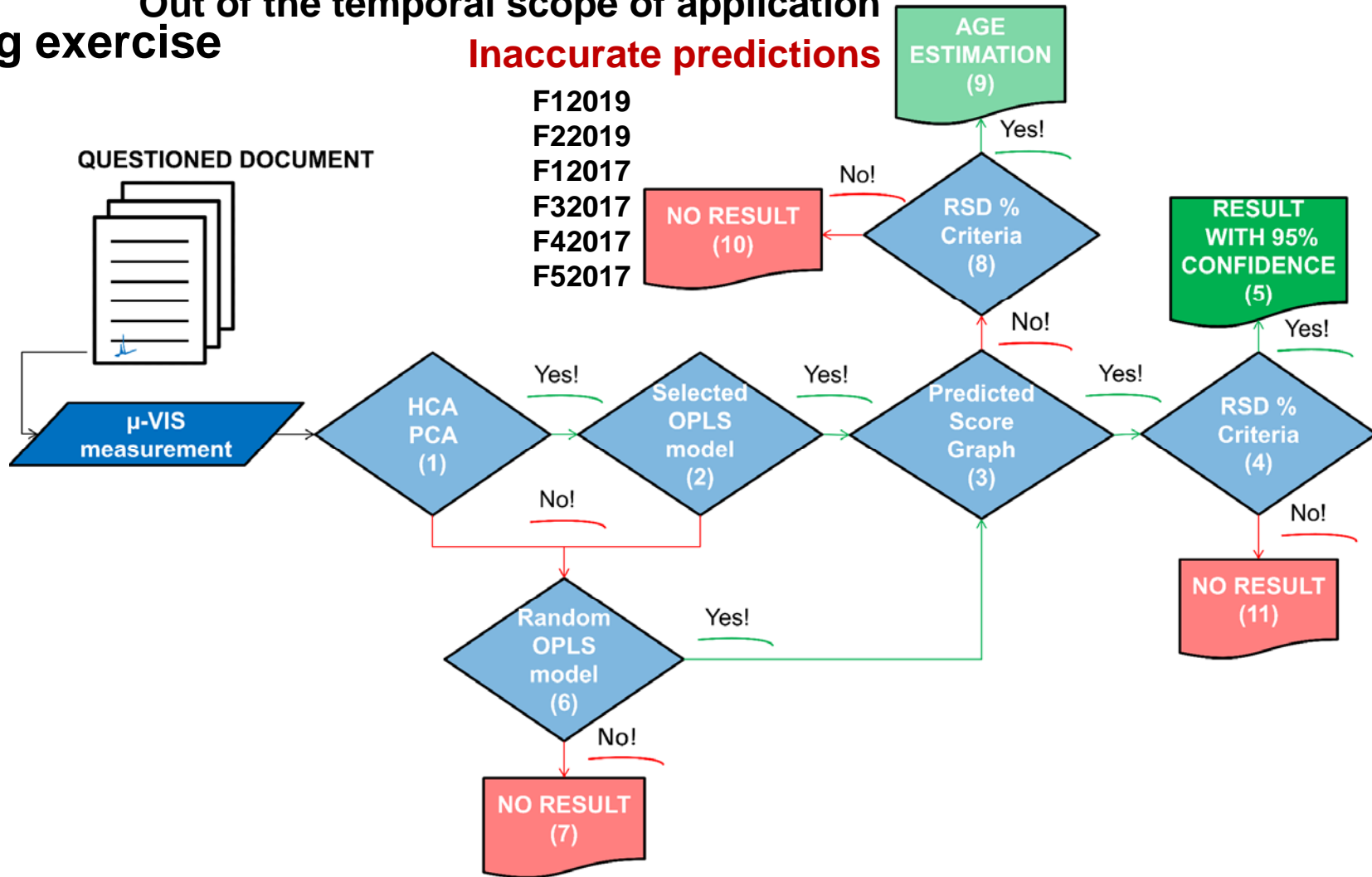


Blind testing exercise



Blind testing exercise

Out of the temporal scope of application
Inaccurate predictions



Remarks

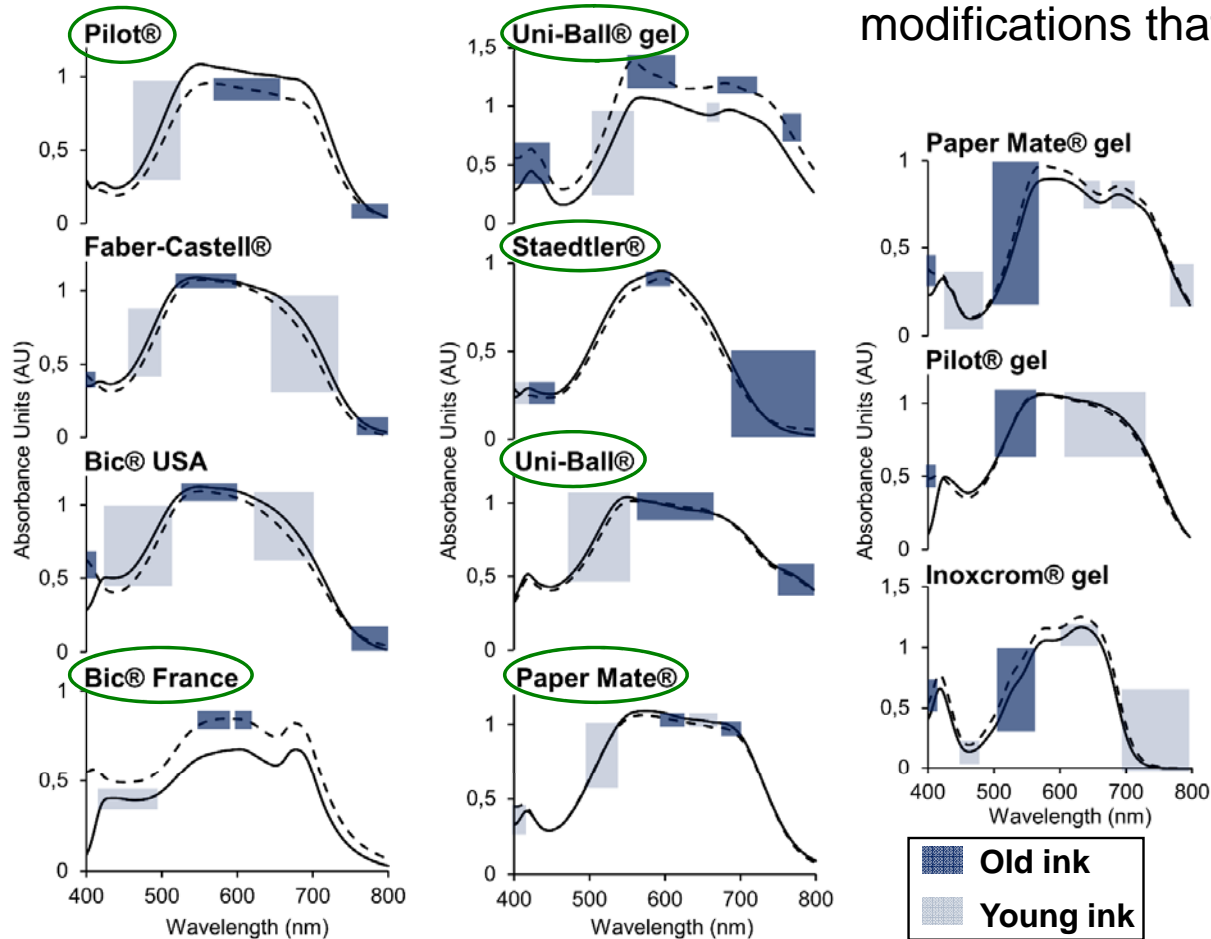
Ability to detect with 95% confidence inks out of their temporal application range

whenever:

- I. The ink is **correctly clustered/classified** in the PCA/HCA model with one of the pen brands studied.
- II. The two **ink replicas** or one of them falls **out of the ellipse** of the predicted score plot.

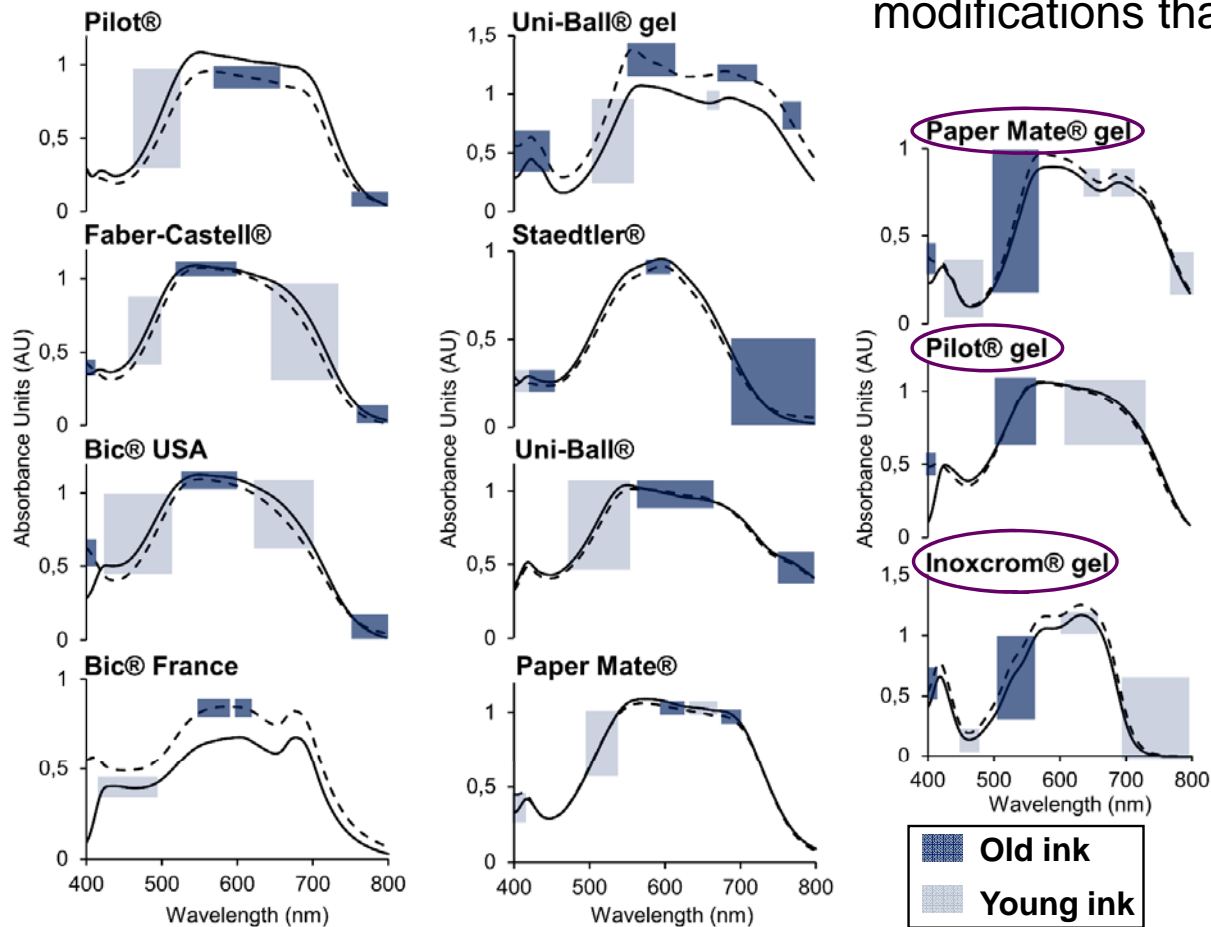
Loadings plot

Ink brands characterized over time by spectral modifications that **shift toward higher wavelengths**



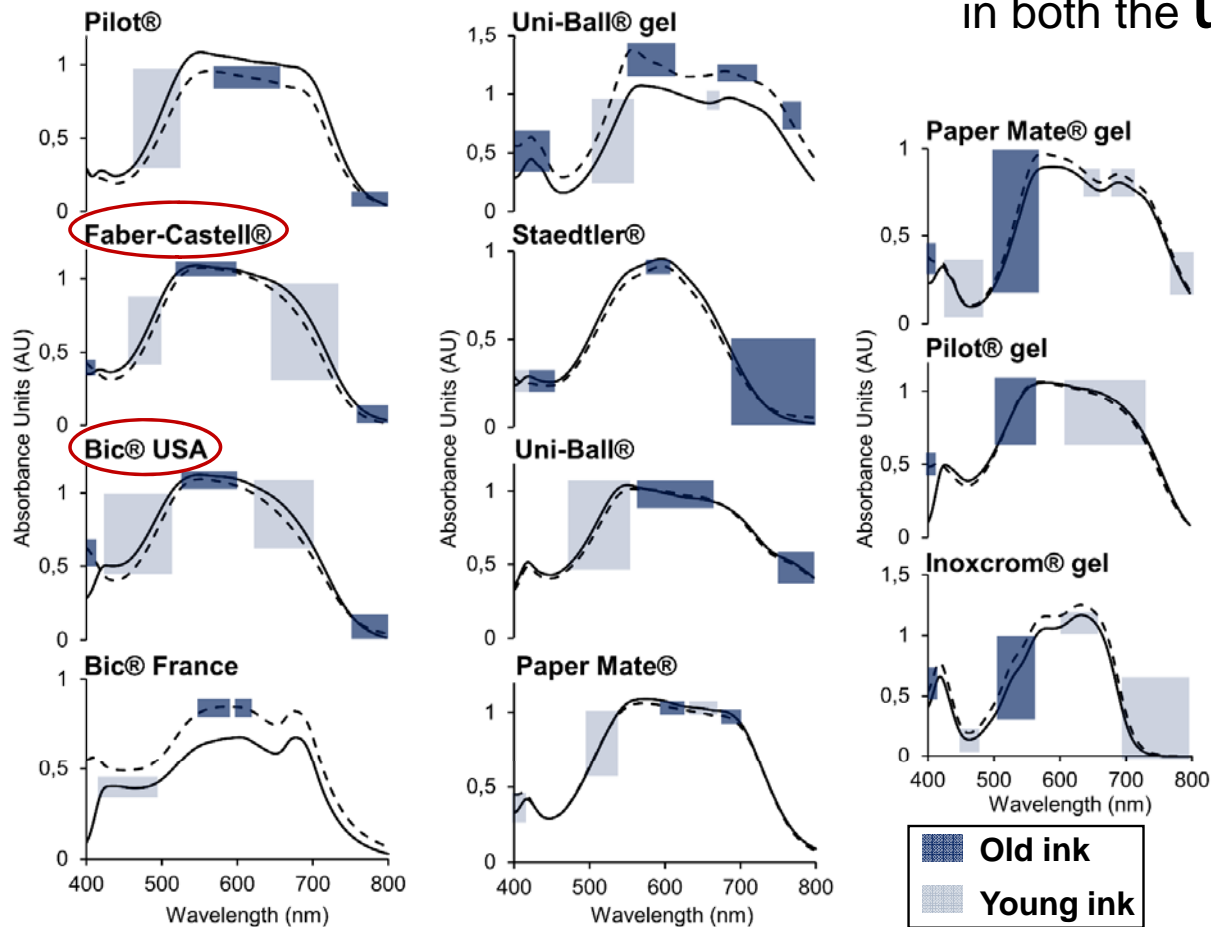
Loadings plot

Ink brands characterized over time by spectral modifications that **shift toward lower wavelengths**



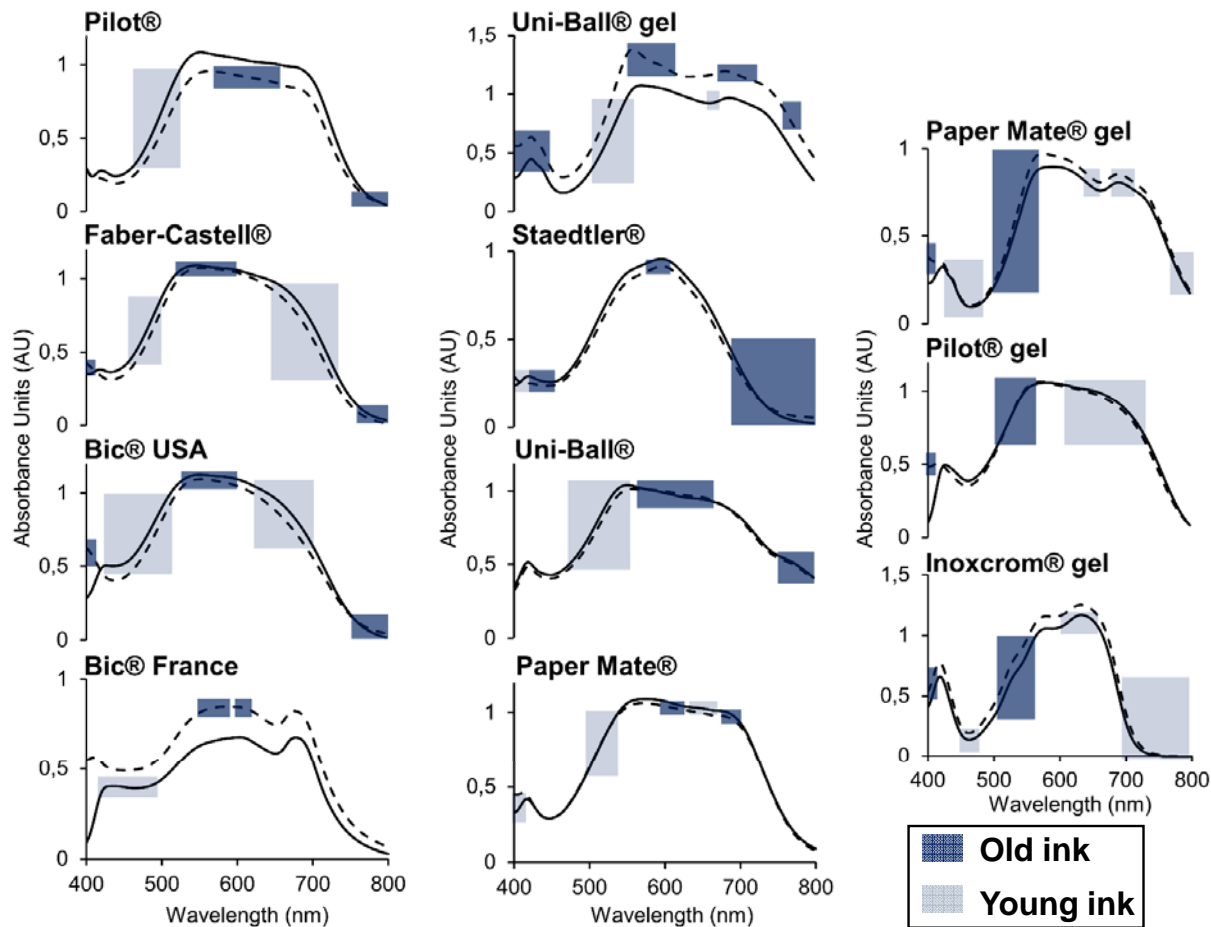
Loadings plot

Ink brands characterized over time by spectral modifications in both the **upper and lower wavelengths**



Loadings plot

Inability to obtain universal OPLS models



- **Multivariate regression methods** demonstrate great **versatile** of:
 - Application to **different forensic fields**.
 - Easily coupling with the most commonly used **analytical techniques**.

- The preparation of **custom-made** synthetic laboratory **samples** aged under **accelerated conditions**:
 - Overcome the **unavailability of samples**.
 - Enable **broad temporal scope** methodologies in a rapid manner.

- Multivariate regression **models** are easily **updated** as new samples are added.

- **New** chemometric-based dating **methodologies**:
 - Respond to **complex real-life claims** that conventional methodologies fail to cope with.
 - Improve and overcome the **disadvantages** or **objectives not reached** by their predecessors.

- The **statistical basis** of chemometrics provides **reliable** and **objective** age estimates.

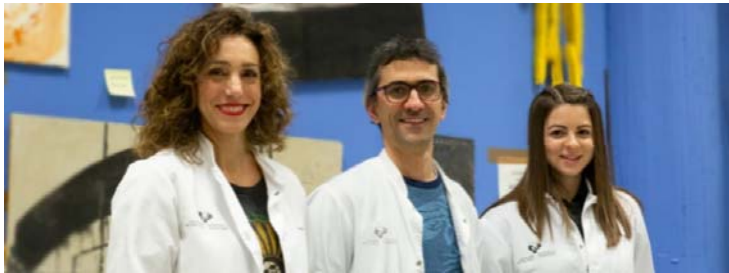
- The **guideline** developed could be a valuable tool for use by the **forensic community**.

- **Testing** for the real **applicability** of the methodologies should involve the **participation** of experts.

Team and collaborating academic and law enforcement institutions



FACULTY OF SCIENCE AND TECHNOLOGY
UNIVERSITY OF THE BASQUE COUNTRY



Contact information:

luis.bartolome@ehu.eus laura.ortiz@ehu.eus