

# EA-IRMS: No Intra-sample Memory for $\delta^{34}\text{S}$ Measurements of Inorganic Sulfur Materials Using the EA IsoLink IRMS System

Christopher Brodie, Oliver Kracht,  
Alexander Hartwig, Andreas Hilker,  
Charles Douthitt  
Thermo Fisher Scientific,  
Bremen, Germany

## Key Words

$\delta^{34}\text{S}$ , Carry-over, EA-IRMS, Intra-sample Memory

## Goal

To demonstrate that the EA IsoLink IRMS System is free from intra-sample memory for  $\delta^{34}\text{S}$  measurements by using inorganic international reference materials.



Figure 1. Thermo Scientific EA IsoLink IRMS System.



## Introduction

Accurate and precise  $\delta^{34}\text{S}$  measurements by EA-IRMS require that the system can process samples without intra-sample effects, generally referred to as “memory effect” or “carry-over”. Any intra-sample memory effect or carry-over will produce inaccurate  $\delta^{34}\text{S}$  values and significantly reduce data reproducibility as well as reducing system productivity and consuming valuable laboratory resources.

This application brief demonstrates that the Thermo Scientific™ EA IsoLink™ IRMS System (Figure 1) does not show an intra-sample memory effect on  $\delta^{34}\text{S}$  measurements and that accurate and precise  $\delta^{34}\text{S}$  values are produced.

**Intra-sample performance for  $\delta^{34}\text{S}$  analysis**

Figure 2 illustrates a series of 30 sequential  $\delta^{34}\text{S}$  measurements of international standard reference materials IAEA S1, IAEA S2, IAEA S3, IAEA SO5 and IAEA SO6. Samples were weighed in amounts of

200–280  $\mu\text{g}$ , which equates to 25–30  $\mu\text{g}$  of sulfur. A 3-point calibration was produced, with an  $r^2 = 0.999$ . The standard deviation of the mean is 0.19‰ across a 56‰ range (Figure 3). The  $\delta^{34}\text{S}$  precision was  $\leq 0.18\text{‰}$  for  $n = 6$  showing excellent accuracy (Table 1).

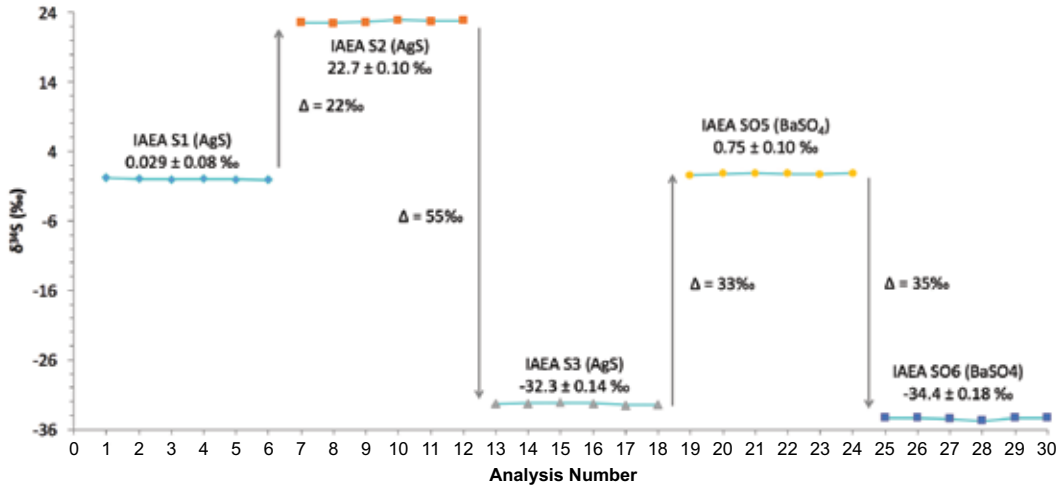


Figure 2. Sequence of 30 sequential measurements of international reference materials.

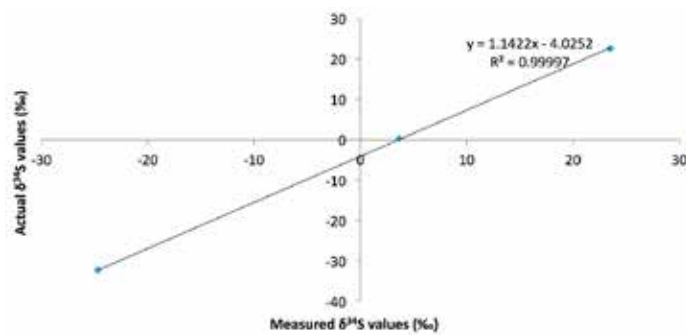


Figure 3.  $\delta^{34}\text{S}$  3-point calibration curve.

**Summary**

The analysis of inorganic sulfur materials shows no intra-sample memory or carry-over effect on  $\delta^{34}\text{S}$  measurements on the EA IsoLink IRMS System, therefore:

- There is no need to discard the first or second sample points from replicate measurements prior to analyzing the data.
- There are no trends in the replicate analysis of individual samples.
- There is no requirement to undertake memory correction.

The data presented in Table 1 are application data and are not warranted because they exceed product specifications. The warranted product specification for  $\delta^{34}\text{S}$  is  $\pm 0.2\text{‰}$  (1 sd) for 50  $\mu\text{g}$  of sulfur measured on Sulfanilamide.

Table 1. Summary of measured  $\delta^{34}\text{S}$  values and actual  $\delta^{34}\text{S}$  values.

Sample	Weight (mg)	n	$\mu\text{g S}$	Measured $\delta^{34}\text{S}_{\text{VCDT}}$ (‰) [Mean $\pm 1\sigma$ ]	Actual $\delta^{34}\text{S}_{\text{VCDT}}$ (‰) [Mean $\pm 1\sigma$ ]
IAEA S1 (AgS)	0.202–0.285	6	30.4	0.03 $\pm$ 0.08	-0.30 $\pm$ 0.20
IAEA S2 (AgS)	0.192–0.219	6	25.8	22.67 $\pm$ 0.10	22.70 $\pm$ 0.20
IAEA S3 (AgS)	0.191–0.234	6	26.0	-32.30 $\pm$ 0.15	-32.30 $\pm$ 0.20
IAEA SO5 (BaSO <sub>4</sub> )	0.196–0.268	6	30.4	0.75 $\pm$ 0.10	0.50 $\pm$ 0.20
IAEA SO6 (BaSO <sub>4</sub> )	0.188–0.272	6	29.2	-34.43 $\pm$ 0.18	-34.1 $\pm$ 0.20

Find out more at [thermofisher.com/EALink](http://thermofisher.com/EALink)