

# Nitrogen/Protein determination in flours by the Thermo Scientific FlashSmart Elemental Analyzer using argon as carrier gas

## Authors

Dr. Liliana Krotz, Walter Galotta,  
and Dr. Guido Giuzzi  
Thermo Fisher Scientific, Milan,  
Italy

## Keywords

Argon, Combustion, Flour, Food  
Quality, Labeling, Nitrogen,  
Protein

## Goal

This application note presents data on Nitrogen/Protein determination in flour reference materials with different nitrogen concentrations to demonstrate the performance of the system using argon for food quality and labeling purposes, while showing compliance to international standards requirements.

## Introduction

Flour is a fine powder made by grinding cereal grains or other suitable edible plant matter which contains starch as polysaccharides and protein. It is most commonly made from wheat but also rye, barley, rice and corn (similar to “polenta”). Ground legumes and nuts, such as soy, peanuts, almonds, and other tree nuts are also called flours. It is the key ingredient of bread; a staple food in many countries, and its availability has often been a major economic and political issue.

Regulations in many countries require specific control of flour quality. One of the most important parameters is protein content, which can be monitored through the precise and accurate determination of the levels of nitrogen in the product to determine nutritional quality.

Helium supply has been beset with problems over the last few years, with world-wide shortages and subsequent price increases. It has therefore been necessary to test an alternative gas, argon, which is readily available.

The Thermo Scientific™ FlashSmart™ EA (Figure 1) copes effortlessly with a wide array of laboratory requirements such as accuracy, day by day reproducibility and high sample throughput. It operates according to the dynamic flash combustion of the sample and typically uses helium as the carrier gas.



Figure 1. FlashSmart Elemental Analyzer.

## Methods

Samples are weighed in tin containers and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler together with the proper amount of oxygen. After combustion, the resultant gases are carried by an argon flow to a second reactor filled with copper, then swept through CO<sub>2</sub> and H<sub>2</sub>O traps, a GC column and are finally detected by a Thermal Conductivity Detector (TCD). The analytical configuration as well as the TCD Detector are as the same as those used with helium as the carrier gas (see Figure 2).

A complete report is automatically generated by the Thermo Scientific™ EagerSmart™ Data Handling Software and displayed at the end of the analysis. The EagerSmart Data Handling Software provides the option AGO (Argon Gas Option) which modifies the argon carrier flow during the run to optimize the analysis.

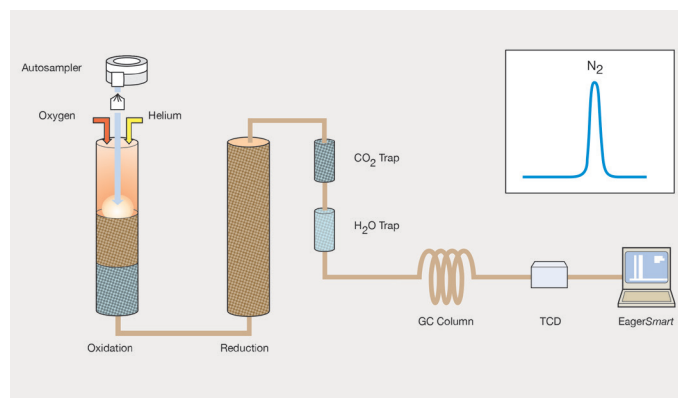


Figure 2. Nitrogen configuration.

## Analytical Conditions

Combustion Furnace Temperature	950 °C
Reduction Furnace Temperature	840 °C
Oven Temperature	50 °C (GC column inside the oven)
Argon Carrier Flow	60 ml/min
Argon Reference Flow	60 ml/min
Oxygen Flow	300 ml/min
Oxygen Injection End	30 sec for standard, Category B (OxyTune) for flours and 30 sec for soy bean meal
Sample Delay	10 sec
Run Time	10 mins

## Official methods requirements for protein determination by combustion

AOAC (Method 992.23) and AACC (Method 46-30, 1999) indicate that a suitable fineness of grind must be determined (for each different material analyzed) to achieve precision that gives RSD of ≤ 2% for 10 successive determinations of nitrogen.

## Results

To evaluate the performance of the system for flour analysis, samples with different nitrogen content were chosen. Instrument calibration was performed with approximately 50–70 mg of aspartic acid standard (10.52 N%), using K factor as calibration method. To evaluate the calibration, Thermo Scientific Pasta Reference Material was analyzed. The protein factor 6.25 was used to calculate the protein content.

Table 1 shows the Nitrogen/Protein data obtained of the analysis of Pasta Reference Material. The sample weight was 150–160 mg. The certified N% is 2.227 and the uncertainty declared by the supplier is 0.097.

**Table 1. Nitrogen/Protein data of pasta reference material.**

N%	RSD%	Protein %	RSD%
2.21	0.542	13.81	0.535
2.23		13.94	
2.19		13.69	
2.20		13.75	
2.22		13.87	
2.21		13.81	
2.20		13.75	
2.22		13.87	
2.21		13.81	
2.20		13.75	

Five Flour Reference Materials from 1.36 to 7.55 N% were chosen to correlate the experimental results to the expected values. Table 2 shows the certified N% and the uncertainty declared by the supplier.

**Table 2. Expected Nitrogen/Protein data of flour reference materials.**

Sample description	Supplier specification	
	N%	Uncertainty ( $\pm$ )
Wheat flour reference material	1.36	0.25
Rice flour reference material	1.38	0.05
Barley flour reference material	1.9	0.04
Oatmeal reference material	1.9	0.10
Soy bean meal reference material	7.5	0.05

Table 3 shows the experimental N/Protein data obtained which were very satisfactory. The sample weight was 140–150 mg.

In all cases the RSD% was less than 2% as requested in the Official Methods.

**Table 3. Experimental Nitrogen/Protein data of flours reference material.**

Sample	Wheat flour		Rice flour		Barley flour		Oatmeal	
	N%	Protein %	N%	Protein %	N%	Protein %	N%	Protein %
%	1.37	8.57	1.36	8.49	1.89	11.80	1.88	11.76
	1.36	8.51	1.37	8.59	1.91	11.94	1.88	11.73
	1.34	8.36	1.40	8.73	1.91	11.91	1.86	11.61
	1.35	8.41	1.40	8.74	1.91	11.95	1.90	11.90
	1.36	8.52	1.40	8.75	1.94	12.10	1.86	11.64
	1.36	8.49	1.38	8.65	1.91	11.91	1.85	11.56
	1.33	8.28	1.38	8.65	1.87	11.66	1.92	11.97
	1.36	8.51	1.38	8.62	1.89	11.80	1.87	11.67
	1.33	8.30	1.41	8.83	1.90	11.89	1.92	12.01
	1.36	8.52	1.39	8.69	1.92	11.97	1.89	11.79
Average %	1.35	8.45	1.39	8.67	1.90	11.89	1.88	11.76
RSD%	1.034	1.210	1.130	1.106	0.997	0.997	1.302	1.307

Table 4. Experimental Nitrogen/Protein data of soy bean meal reference material (130–190 mg).

Weight (mg)	N%	Av. N%	RSD%	Protein %	Av. Protein %	RSD%
130.48 132.67 129.90	7.50 7.51 7.50	7.50	0.077	46.86 46.91 46.84	46.87	0.077
140.57 141.41 140.28	7.49 7.53 7.45	7.49	0.534	46.81 47.05 46.57	46.81	0.513
161.30 160.29 162.44	7.48 7.51 7.58	7.52	0.682	46.73 46.95 47.38	47.02	0.703
170.53 171.02 171.22	7.47 7.53 7.51	7.50	0.407	46.66 47.06 46.91	46.88	0.431
182.89 182.54 180.50 91.57	7.49 7.56 7.55 7.54	7.53	0.413	46.81 47.23 47.22 47.12	47.09	0.417

## Conclusions

Good repeatability, accuracy and precision was obtained with the FlashSmart Elemental Analyzer using argon as the carrier gas. The RSD% obtained was less than 2% according to the AOAC and AACC Official Methods Performance Requirements.

No memory effect was observed when changing the type of sample, indicating complete combustion and detection of the element.

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