

# Elemental Analysis: Nitrogen/Protein determination in starch by flash combustion in alternative to Kjeldahl method

Author: Liliana Krotz, Thermo Fisher Scientific, Milan, Italy

Keywords: CHNS, combustion, nitrogen, protein, starch, slurry

## Goal

To assess the performance of the elemental analyzer for N/Protein determination for low and trace nitrogen content in starch and CHNS determination.

## Introduction

As starch is the most common carbohydrate in the human diet (being contained in large amounts in staple food), and since it is largely used for diverse industrial uses, its elemental content must be monitored and tested for quality control.

Pure starch is a white, tasteless and odorless powder. Starch is largely used in the food industry as food additive, modified starches and sugars. For the starch refining process and from a food safety point of view, its protein content is used as an important parameter. As additive for food processing, food starches are typically used as thickeners and stabilizers in foods, such as puddings, custards, soups, sauces, gravies, pie fillings, and salad dressings, and are exceptional binders in processed meats.



For the production of ethanol in the manufacturing of beer, whisky and biofuel, starch is converted into sugars. Many of the sugars used in processed foods are produced by using starch. Mixing starches in warm water produces a paste, which can be used as a thickening, stiffening or gluing agent.

Starch is also used in industries other than food: for example as an adhesive in the papermaking process, as additive before ironing, as an excipient, as tablet disintegrant, and as binder for pharmaceutical uses.

In the production process, the protein content of starch, calculated through the nitrogen determination, is periodically monitored, and tested for quality control. Considering that starch is used also in the preparation of animal feed, the determination of N/Protein is critical. Therefore, it is very important to have an automated method allowing the analysis of N/Protein with an excellent reproducibility.

The Dumas method (combustion method) allows faster, safer, and more reliable determination of nitrogen than the traditional Kjeldahl method and has been approved by different associations.

The Thermo Scientific™ FlashSmart™ Elemental Analyzer (Figure 1), based on the dynamic flash combustion of the sample, meets a wide array of requirements of laboratories such as stability, accuracy, day by day reproducibility and high sample throughput.



Figure 1. Thermo Scientific FlashSmart Elemental Analyzer.

The FlashSmart Elemental Analyzer provides rapid and automated nitrogen determination without the use of hazardous chemicals, offering advantages over traditional methods. The FlashSmart EA allows analyses of high and low nitrogen levels with no configurational change and without matrix effect. The protein content is calculated automatically using the dedicated conversion factor in the Thermo Scientific™ EagerSmart™ Data Handling Software.

The FlashSmart EA also allows the simultaneous determination of nitrogen, carbon, hydrogen and sulfur by combustion method using the same system with a specific single combustion-reduction reactor for the complete characterization of all type of matrices.

## Method

The Elemental Analyzer operates according to the dynamic flash combustion of the sample. The sample is weighed in tin containers and introduced into the combustion reactor via the Thermo Scientific™ MAS Plus Autosampler with oxygen.

For nitrogen determination, after combustion, the produced gases are carried by a helium 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 finally detected by a Thermal Conductivity Detector (TCD) (Figure 2). A complete report is automatically generated by the EagerSmart Data Handling Software and displayed at the end of the analysis. The EagerSmart Data Handling Software controls all analytical parameters of the instrument including the oxygen flow and the timing of oxygen injection. It automatically calculates the amount of oxygen, relative to the sample matrix and sample weight, through the dedicated Thermo Scientific™ OxyTune Function (patented) ensuring the complete combustion of the sample. Through this optimization also decreases the cost per analysis by not wasting oxygen or consuming the copper unnecessarily. Figure 3 shows the OxyTune Categories.

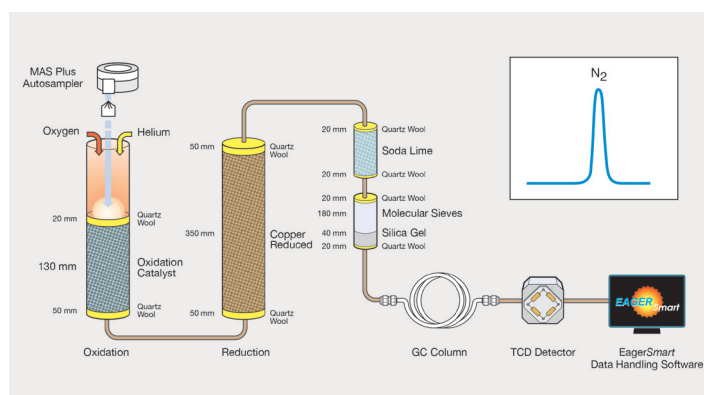


Figure 2. Nitrogen configuration.

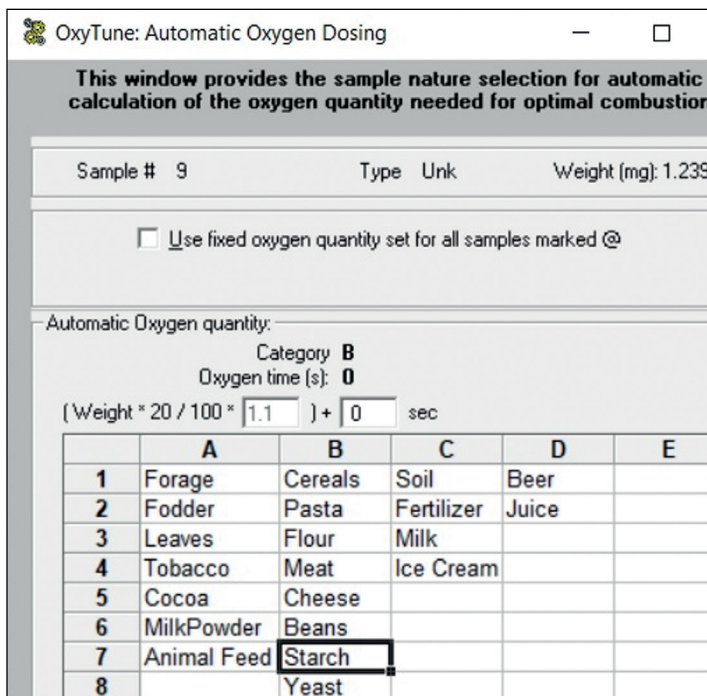


Figure 3. OxyTune EagerSmart Data Handling Software window.

For CHNS determination, after combustion, the produced gases are carried in a helium carrier gas to a layer filled with copper. The analyte then enters the GC column, which separates the produced gases before detection by a Thermal Conductivity Detector (TCD) (Figure 4). For weight percent determination a complete report is automatically generated by the EagerSmart Data Handling Software and displayed at the end of the analysis.

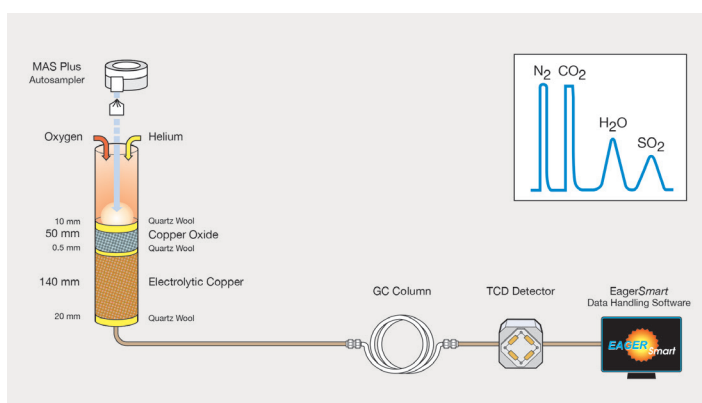


Figure 4. CHNS configuration.

## Results

The starch samples analyzed were chosen on basis of the different nitrogen content. The calibration of the Analyzer was performed with 10–12 mg of atropine using K factor as calibration method. Due to the low nitrogen content present in starch samples, the blank was evaluated by the analysis of an empty tin container. The starch sample weight was about 150–250 mg.

Table 1 shows the nitrogen results obtained of starch samples at 0.25–0.35 N% analyzed in triplicate. While Table 2 shows the N/Protein data obtained of starch and slurry starch samples at trace level (100–500 ppm N) analyzed in triplicate. Slurry starch is a combination of starch (usually corn starch, flour, potato starch or arrowroot) and cold water which is mixed together and used to thicken a soup or sauce. If the starch is solely added directly to a hot liquid, the starch granules cannot disperse easily and clumps form. Once mixed with water, the slurry can be added directly to the hot liquid.

Table 1. Nitrogen data of starch samples.

Sample	N%	RSD%
1	0.2531	0.65
	0.2558	
	0.2561	
2	0.3963	0.64
	0.3991	
	0.3940	
3	0.3589	0.80
	0.3563	
	0.3532	
4	0.3029	0.61
	0.3012	
	0.3049	

**Table 2. N/Protein data of starch and slurry starch samples at trace level.**

Sample	N ppm	Protein %	RSD%
Starch 1	124	0.0775	1.38
	126	0.0789	
	123	0.0768	
Starch 2	162	0.1015	1.54
	167	0.1046	
	164	0.1025	
Starch 3	177	0.1107	0.59
	179	0.1120	
	178	0.1115	
Starch 4	351	0.2191	0.39
	350	0.2185	
	352	0.2202	
Starch 5	371	0.2322	0.83
	366	0.2285	
	370	0.2312	
Starch 6	380	0.2378	0.89
	382	0.2387	
	387	0.2419	
Starch 7	404	0.2528	0.18
	404	0.2525	
	403	0.2519	
Starch 8	476	0.2975	0.25
	474	0.2962	
	474	0.2962	
Starch 9	484	0.3024	0.15
	483	0.3016	
	484	0.3024	
Starch 10	546	0.3409	0.53
	551	0.3445	
	548	0.3426	
Slurry starch 1	146	0.0912	1.35
	150	0.0937	
	148	0.0925	
Slurry starch 2	166	0.1038	1.29
	165	0.1029	
	162	0.1012	

At last the data of nitrogen, carbon, hydrogen and sulfur obtained simultaneously by combustion method is shown in Table 3. The calibration was performed with 2–3 mg BBOT standard (2,5-Bis (5-ter-butyl-benzoxazol-2-yl) thiophene) using K factor as calibration method. The calibration was evaluated by the analysis of BBOT and aspartic acid as unknown. Table 3 shows the CHNS data of the samples analyzed five times (sample weight 3–4 mg).

**Table 3. CHNS data of a starch sample.**

N%	RSD%	C%	RSD%	H%	RSD%	S%	RSD%
2.53	0.33	31.01	0.20	5.40	1.00	0.399	0.90
2.52		30.85		5.45		0.396	
2.54		31.00		5.42		0.391	
2.52		30.96		5.37		0.392	
2.53		30.97		5.31		0.398	

### Conclusions

For nitrogen determination of starches, the Thermo Scientific FlashSmart Elemental Analyzer, based on the combustion method (Dumas), offers advantages over the Kjeldahl Method in terms of automation, ease of use and cost per sample. Differently from the traditional methods, the FlashSmart Elemental Analyzer doesn't require the use of toxic chemicals.

For N/Protein and CHNS determination, good repeatability was obtained, and no memory effect was observed, indicating complete combustion and detection of the element independent of the element content.

With the FlashSmart Analyzer laboratories can analyze nitrogen in a wide range from high to trace nitrogen content.

The Dumas Combustion method has been approved and adopted by several Official Organizations such as ASBC, AOAC, AACC, AOCS, IDF, IFFO and ISO.

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

**For Research Use Only. Not for use in diagnostic procedures.** © 2021 Thermo Fisher Scientific Inc. All rights reserved. AOAC is a trademark of The Association of Official Analytical Chemists; AOCS is a trademark of The American Oil Chemists' Society; AACC is a trademark of The American Association of Cereal Chemists; ASBC is a trademark of The American Society of Brewing Chemists; IFFO is a trademark of the International Fishmeal and Fish Oil Organization; ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representatives for details. Not all products are available in all countries. Please consult your local sales representatives for details. **AN42514-EN 0121C**

**ThermoFisher**  
SCIENTIFIC