

# Quality control of choline as a dietary supplement by high performance liquid chromatography coupled to a charged aerosol detector

Authors: Katherine Lovejoy<sup>1</sup> and Paul Gamache<sup>2</sup>

<sup>1</sup>Thermo Fisher Scientific, Germering, Germany

<sup>2</sup>Thermo Fisher Scientific, Chelmsford, MA, USA

Keywords: Charged aerosol detection, CAD, bitartrate, choline, chloride, citrate, hydrophilic interaction chromatography, HILIC, impurities, United States Pharmacopeia, USP, Vanquish Flex

## Goal

To develop a single method with direct detection of choline salts and impurities without the need for derivatization.

## Application benefits

- Direct and simultaneous detection of choline, its counterions, and impurities
- Suitable for assay of choline content and impurities and identification of counterions in commercially available dietary supplements
- Method compatible for both charged aerosol detection and mass spectrometry for determination of impurities



## Introduction

Choline is an essential nutrient involved in cell membrane integrity, lipid transport and metabolism, neurotransmission, and many other functions.<sup>1</sup> The main food sources of choline are animal-based products, beans, and nuts, which contain various forms including free choline, phosphocholine, and phosphatidylcholine.<sup>2</sup> Choline deficiency can lead to muscle damage, liver damage, and nonalcoholic fatty liver disease.<sup>2</sup> To ensure nutritional adequacy, the use of dietary supplements containing choline (e.g., choline bitartrate, choline chloride, choline citrate, phosphatidylcholine, and lecithin) is common.<sup>3</sup> Since the quality of dietary supplements can vary greatly, laboratory testing of raw materials and final products is often needed to ensure safety and effectiveness.

Traditionally, high performance liquid chromatography (HPLC) with ultraviolet (UV) detection is used for the analysis of choline salts. As choline salts respond poorly to this approach, precolumn derivatization with reagents containing a strong chromophore is required. However, this approach is limited: derivatization may interfere with the counter ions and may miss impurities that do not react.<sup>4</sup>

In this study, HPLC with a charged aerosol detector (CAD) was used for measurement of content and purity of choline bitartrate, choline citrate, and choline chloride (see Figure 1 for structures). A mixed-mode column operated in HILIC mode with an ammonium acetate buffered mobile phase and acetonitrile-water gradient was used. This approach enabled the separation of choline, its counterions, and several impurities. CAD provided direct detection of these analytes, thus avoiding the need for derivatization used in older methods. Several impurities were identified based on retention time comparison to individual standards or by mass spectrometry and include Na<sup>+</sup>, K<sup>+</sup>, Cl<sup>-</sup> and O-(2-hydroxyethyl)choline (CAS 41830-55-1) (see Figure 1 for structure). Choline content was measured over the range of 0.08 to 0.12 mg/mL with precision  $\leq 3.0\%$  RSD, recovery from 96.7 to 100.8%, and correlation coefficient ( $r^2$ ) for linear regression of  $> 0.997$ .

This HPLC-CAD method provides advantages over derivatization techniques by allowing more comprehensive measurement of sample composition using a simplified approach. This method is recommended for the analysis of choline citrate in dietary supplements.<sup>4</sup>

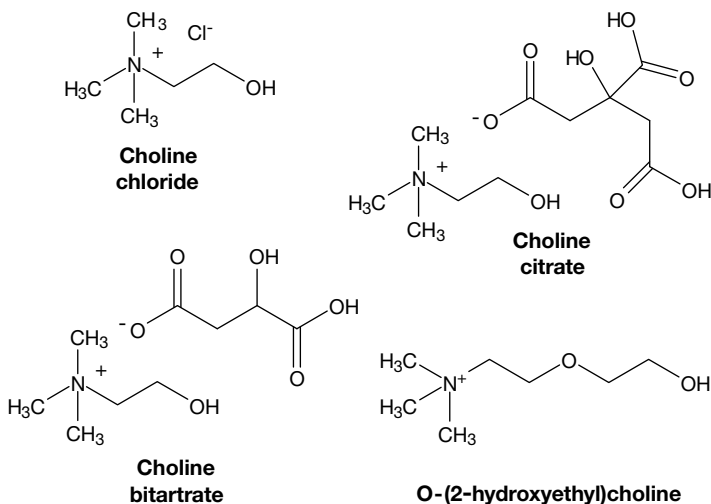


Figure 1. Structures of choline salts and the impurity O-(2-hydroxyethyl)choline

## Experimental

### Reagents and standards

| Chemical name   | Part number |
|---|-------------|
| Deionized water, 18.2 MΩ·cm resistivity or higher—from a Thermo Scientific™ Barnstead™ GenPure™ xCAD Plus Ultrapure Water Purification System | 50136149    |
| Fisher Scientific™ acetonitrile, Optima™ LC/MS grade  | A955-212    |
| Fisher Scientific™ ammonium acetate, LC/MS grade  | A114-50     |
| Fisher Scientific™ acetic acid, LC/MS grade   | A113-50     |
| Fisher Scientific™ sodium chloride (>99.9%)   | S/3165/53   |
| Choline bitartrate, choline chloride, and choline citrate were purchased from reputable vendors   |             |

### Equipment

| Item name  | Part number                      |
|--|----------------------------------|
| Fisher Scientific™ Fisherbrand™ Mini Vortex Mixer  | 14-955-152                       |
| Thermo Scientific™ Orion™ 3 Star pH Benchtop Meter   | 13-644-928                       |
| Thermo Scientific™ Finpipette™ F1 Variable Volume Single-Channel Pipettes: 100-1000 μL, 10-100 μL, 1-10 μL | 4641100N<br>4641070N<br>4641030N |
| Thermo Scientific™ PP Crimp/Snap Top autosampler vials   | C401114                          |
| Thermo Scientific™ 11 mm autosampler vial crimp caps (Chlorobutyl, PTFE)                                   | 11568150                         |

## Mobile phase preparation

Prepare an 80.4 mM ammonium acetate buffer by adding 6.2 g of ammonium acetate to 900 mL deionized water and adjust the pH to 4.7 with glacial acetic acid. Bring the volume to 1.0 L in a volumetric flask. This buffer is used for mobile phase and diluent preparation. Note: pH electrodes can be a significant source of contamination causing higher baseline noise and baseline artifacts. For best results, it is recommended to determine the volume of glacial acetic acid needed to adjust the pH and then use only volumetric and gravimetric techniques to prepare buffers used for diluent and mobile phase. This avoids contamination from the pH electrode and typically provides more reproducible retention times between batches.

For mobile phase A, combine 200 mL of the prepared buffer with 800 mL of acetonitrile and mix thoroughly. For mobile phase B, combine 200 mL of the prepared buffer with 300 mL deionized water and 500 mL acetonitrile and mix thoroughly.

## Standard and sample preparation

The diluent for standard and sample preparation was prepared as 80.4 mM ammonium acetate buffer/acetonitrile (30/70, v/v).

Standard solutions of choline bitartrate, choline chloride, and choline citrate were prepared from commercially available standards in diluent. Standard solutions of 0.08 mg/mL, 0.10 mg/mL, and 0.12 mg/mL were used for assay of content and for identification while a 2.0 µg/mL standard solution was used for impurities.

Samples were prepared at 0.10 mg/mL for assay of choline content and at 2.0 mg/mL for impurities. Assay method validation was done based on choline citrate with solutions of 0.08, 0.10, 0.12 mg/mL for calibration and accuracy/precision determination. Instrument precision was calculated as %RSD from six consecutive injections of 0.10 mg/mL solution.

A solution containing related salt counter ions was prepared as 0.10 mg/mL sodium chloride.

## Instrumentation

Chromatographic separation was performed on a Thermo Scientific™ Vanquish™ Flex Quaternary UHPLC system. Analytes were detected using a charged aerosol detector. An optional single quadrupole mass spectrometer was used to help identify impurities and artifacts.

| Module  | Part Number |
|---|-------------|
| Thermo Scientific™ Vanquish™ System Base Flex               | VF-S01-A    |
| Thermo Scientific™ Vanquish™ Quaternary Pump                | VF-P20-A    |
| Thermo Scientific™ Vanquish™ Split Sampler HT               | VH-A10-A    |
| Thermo Scientific™ Vanquish™ Column Compartment H           | VH-C10-A-03 |
| Thermo Scientific™ Vanquish™ Charged Aerosol Detector F     | VF-D20-A    |
| Thermo Scientific™ ISQ™ EM Single Quadrupole Mass Detector* | ISQEM-ESI   |

\*For some experiments, a post-column tee was used to split the flow 1:1 between the CAD and the single quadrupole mass detector.

## Data processing and software

Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS), version 7.2.9 was used for data acquisition and processing.

Table 1. Chromatographic conditions

| Parameter               | Setting  |
|-------------------------|--|
| Column                  | MilliporeSigma™ SeQuant™ ZIC™ –pHILIC, 4.6 × 150 mm, 5 µm (P/N 1.50461.0001)   |
| Mobile phase A          | 80.4 mM ammonium acetate buffer pH 4.7 / acetonitrile (20/80, v/v)   |
| Mobile phase B          | 80.4 mM ammonium acetate buffer pH 4.7 / water / acetonitrile (20/30/50, v/v/v)  |
| Injection volume        | 10 µL  |
| Flow rate               | 0.5 mL/min   |
| Gradient                | See Table 2  |
| Column temperature      | 30 °C still air mode, 30 °C active preheater   |
| Autosampler temperature | 10 °C  |
| CAD conditions          | Data collection rate: 10 Hz; Filter: 3.6 s; Evap T: 35 °C; PFV: 1.00   |
| MS conditions           | Vaporizer temp.: 144 °C; Ion transfer tube: 300 °C; Source voltage: 3000 V (positive), -2000 V (negative); Sheath gas: 32.3 psig; Aux gas: 3.6 psig; Sweep gas: 0.5 psig |
| MS full scan, positive  | Time: 0–35 min; Spectrum type: centroid; Mass range: 50–400 <i>m/z</i> ; Dwell time: 0.1 s; Source CID voltage: 15 V   |
| MS SIM, positive        | Time: 0–35 min; SIM mass: 148.2 <i>m/z</i> ; SIM width: 0.1 amu; Dwell time: 0.1 s; Source CID voltage: 15 V   |
| MS full scan, negative  | Time: 0–35 min; Spectrum type: centroid; Mass range: 50–400 <i>m/z</i> ; Dwell time: 0.1 s; Source CID voltage: 15 V   |

**Table 2. Gradient conditions**

| Time [min] | %A | %B  |
|------------|----|-----|
| 0          | 90 | 10  |
| 3          | 90 | 10  |
| 20         | 0  | 100 |
| 22         | 0  | 100 |
| 26         | 85 | 15  |
| 28         | 90 | 10  |
| 35         | 90 | 10  |

## Results and discussion

Choline, its counterions, and impurities were separated using a mixed-mode column containing a zwitterionic stationary phase using gradient elution (Table 2). Using this approach analytes are resolved by both HILIC and ion exchange mechanisms enabling the simultaneous separation of cations, anions, and neutral analytes. Click [here](#) for further information about mixed-mode columns.

The charged aerosol detector is a universal detector and, under isocratic conditions, shows uniform response for all non-volatile analytes, independent of their chemical structure. For ionic solutes like choline and its counterions, CAD response is based on the salt formed between the analyte and oppositely charged mobile phase additives. In this method, choline is measured as its acetate salt while citrate, bitartrate, and chloride are measured as their ammonium salts.

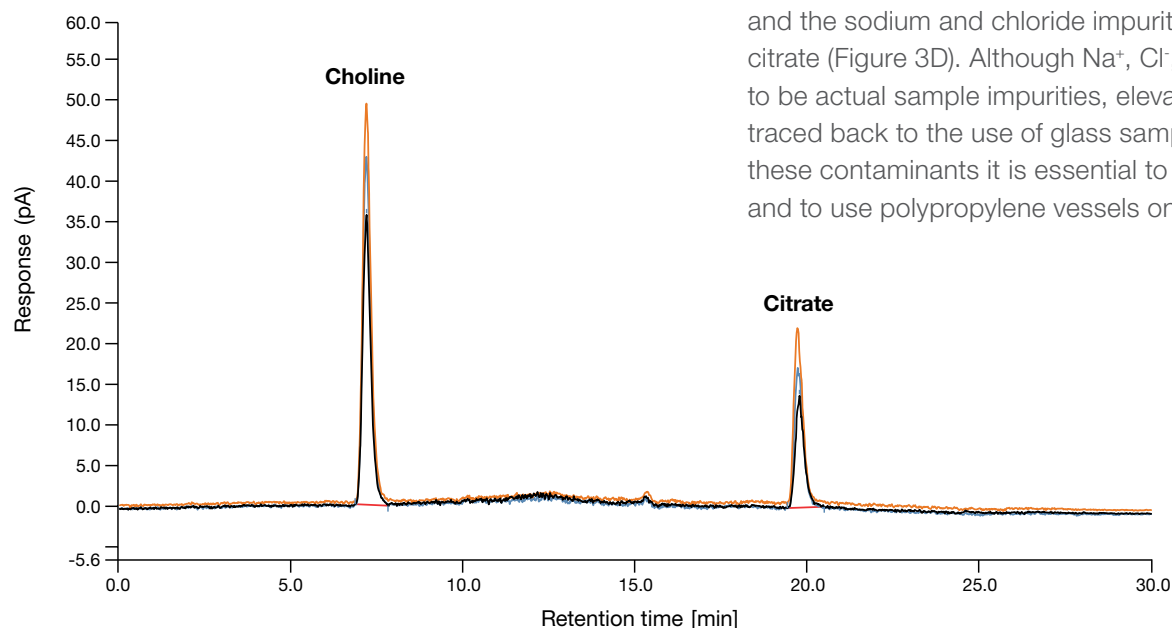


Figure 2. Chromatogram of choline citrate (black = 0.08 mg/mL; blue = 0.10 mg/mL; orange = 0.12 mg/mL)

## Choline citrate method

Figure 2 shows the separation of choline and its counterion citrate at three different concentration levels (0.08, 0.10, and 0.12 mg/mL). Choline was detected as its acetate salt, while citrate was detected as its ammonium salt. Both peaks were well separated and easily quantified.

The validation data for choline citrate analysis are presented in Table 3. The precision was <2% on two separate days. Linearity (0.08, 0.10, 0.12 mg/mL) and recovery (at 80, 100, and 120% of the content specification) were based on linear least squares regression.

Table 3. Validation data for choline citrate on precision, linearity, and recovery

| Parameter                           | Choline citrate |         |         |         |
|-------------------------------------|-----------------|---------|---------|---------|
|                                     | Choline         |         | Citrate |         |
|                                     | Day 1           | Day 2   | Day 1   | Day 2   |
| Precision 0.10 mg/mL (% RSD, n = 6) | 1.83            | 1.23    | 1.33    | 1.76    |
| Linearity ( $r^2$ )                 | 0.99997         | 0.99784 | 0.99778 | 0.99918 |
| Recovery (%) at level               | Day 1           | Day 2   | Day 1   | Day 2   |
| 80                                  | 98.3            | 100.2   | 94.2    | 105.5   |
| 100                                 | 100.8           | 99.4    | 98.1    | 102.2   |
| 120                                 | 99.1            | 96.7    | 99.8    | 98.7    |

## Measurement of other choline salts and impurities

The method can be used to evaluate other choline salts (e.g., choline chloride and choline bitartrate) and impurities (Figure 3), for example, the sodium impurity found in choline tartrate (Figure 3B) and choline chloride (Figure 3C) and the sodium and chloride impurities found in choline citrate (Figure 3D). Although  $\text{Na}^+$ ,  $\text{Cl}^-$ , and possibly  $\text{K}^+$ , seem to be actual sample impurities, elevated levels can also be traced back to the use of glass sample vials. To minimize these contaminants it is essential to avoid the use of glass and to use polypropylene vessels only.<sup>4</sup>

An impurity found in choline citrate samples (RT ~6.0 min) was further studied using a single-quadrupole mass spectrometer (Figure 4). Based on its mass spectrum, the

impurity was determined to be O-(2-hydroxyethyl)choline, a known by-product formed during the production of choline hydroxide.

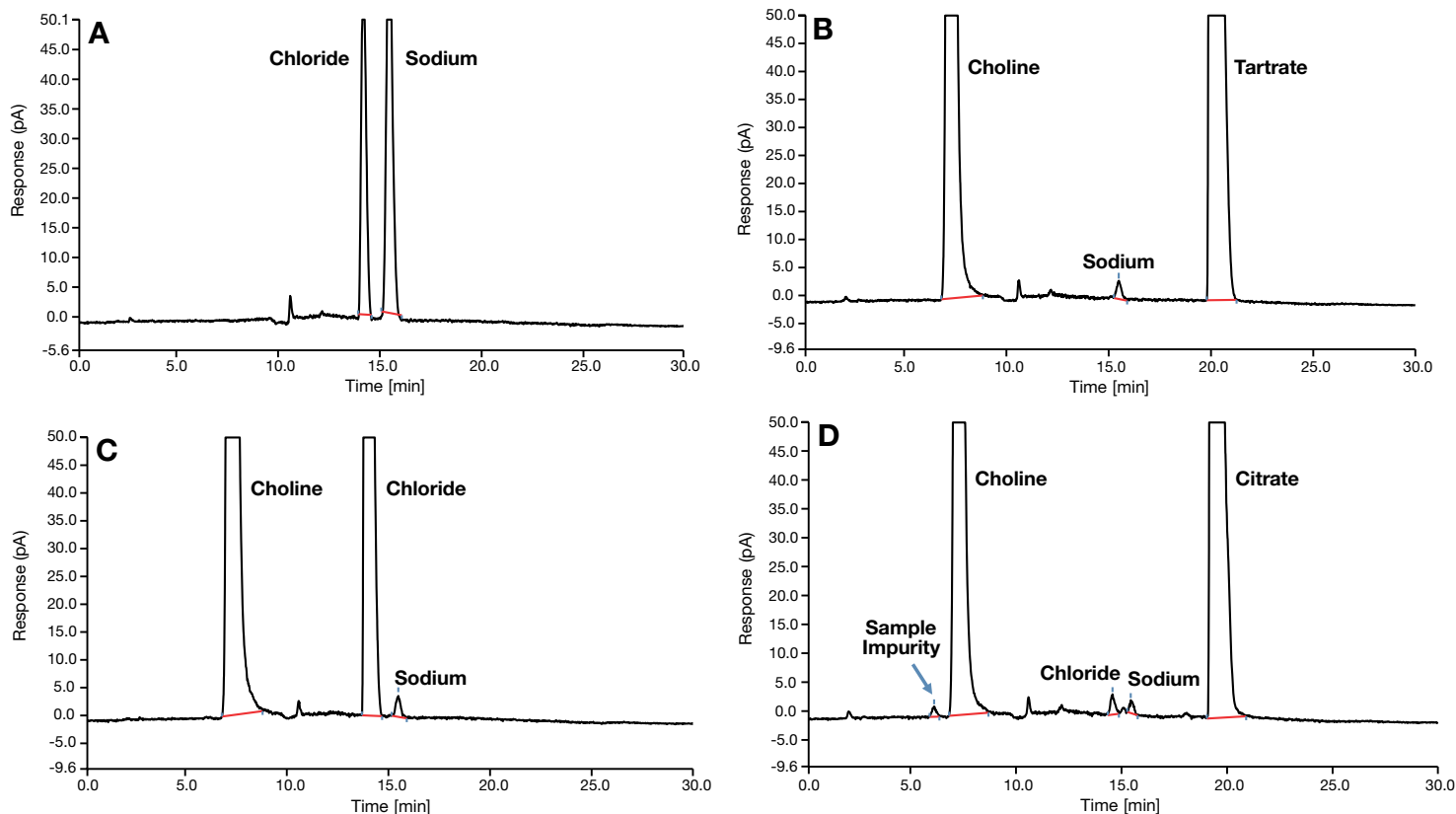


Figure 3. (A) Chromatogram of sodium chloride (0.1 mg/mL)—sodium detected as its acetate salt, chloride as its ammonium salt; (B) analysis of choline bitartrate (2.0 mg/mL) showing sodium impurity; (C) analysis of choline chloride (2.0 mg/mL) showing sodium impurity; (D) analysis of choline citrate (2.0 mg/mL) showing sodium, chloride, and a choline-related sample impurity (eluting at 6.0 min)

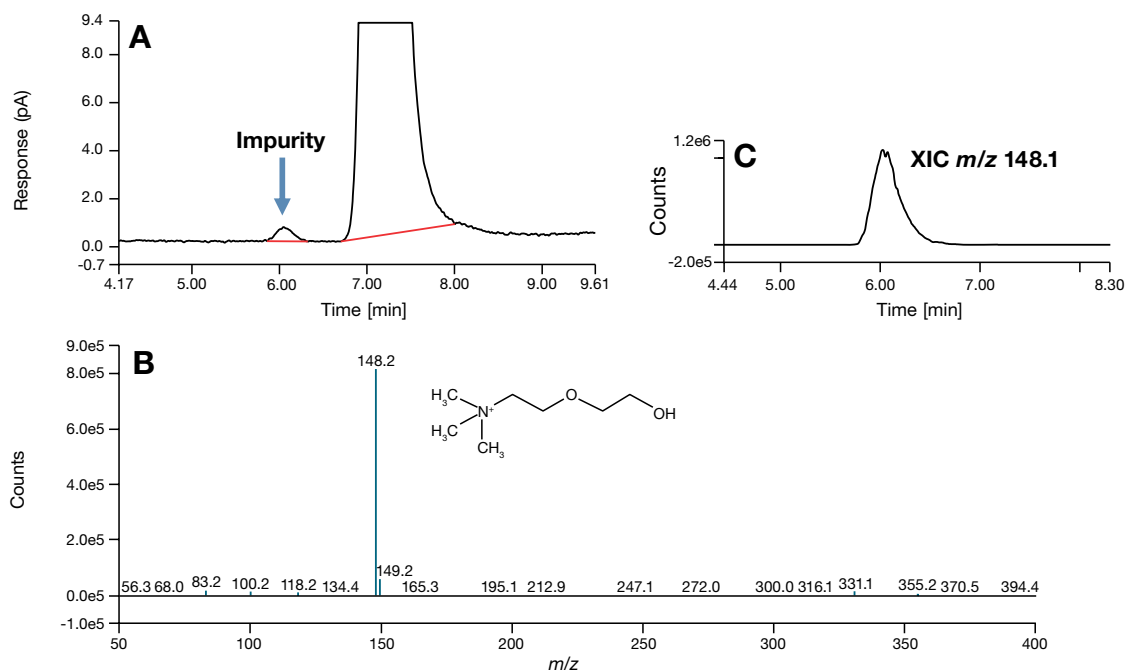


Figure 4. (A) CAD chromatogram of choline citrate (2.0 mg/mL) showing an impurity at RT ~6 min; (B) mass spectrum at the peak apex of the impurity peak at RT ~6 min; (C) extracted ion chromatogram (XIC) of  $m/z$  148.1

## Conclusion

- Mixed-mode chromatography with charged aerosol detection enables the simultaneous analysis of anions and cations, and was used to evaluate choline bitartrate, choline chloride and choline citrate samples
- The low ng sensitivity of CAD allows control of the O-(2-hydroxyethyl)choline impurity and individual unspecified impurities within acceptance criteria of not more than (NMT) 0.1%.<sup>4</sup>
- Use of MS in parallel with CAD facilitated the structural identification of choline-related impurity O-(2-hydroxyethyl)choline in commercial samples.

## References

1. Zeisel, S. H. Dietary Choline: Biochemistry, Physiology, And Pharmacology. *Ann. Rev. Nutr.* **1981**, *1*, 95–121.
2. Zeisel, S. H.; da Costa, K-A. Choline: An Essential Nutrient for Public Health. *Nutr. Rev.* **2009**, *67*, 615–623.
3. Choline—Linus Pauling Institute, Micronutrient Information Center.
4. Choline USP monograph. This document is in the PF stage (USP's journal of public notice and comment). It is not official and may never become official.

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