

Determination of gentamicin and related impurities in gentamicin sulfate

Jingli Hu, Manali Aggrawal, and Jeffrey Rohrer, Thermo Fisher Scientific, 1214 Oakmead Parkway, Sunnyvale, CA, USA, 94085

ABSTRACT

Purpose: To demonstrate that the USP Gentamicin Sulfate monograph Content of Gentamicins method and the USP in-process revision Gentamicin Sulfate monograph method for organic impurities method can be successfully executed with a Thermo Scientific™ Dionex™ IonPac™ AmG-3µm C18 column on a Thermo Scientific™ Dionex™ ICS-5000* HPIC™ system.

Methods: The analysis of gentamicin sulfate based on an ion-pairing HPLC-pulsed amperometric detection (PAD) method is described in the U.S. and European Pharmacopoeias (USP and EP). The eluent of the USP and EP monograph methods contains trifluoroacetic acid, pentafluoropropionic acid, and acetonitrile, and its pH is adjusted to 2.6 with sodium hydroxide to avoid silica-bonded phase hydrolysis when exposed to lower pH conditions. Dionex IonPac AmG-3µm C18 columns are specifically designed for superior resistance towards acidic conditions. Therefore, an aqueous TFA solution can be used as the eluent without adjusting its pH to a higher value. In addition to using the USP/EP method, we modified the eluent in the USP/EP monograph in two ways, with each modification used to make a new method. Method A is a simple eluent method that uses a 100 mM TFA eluent. Method B is a fast method that includes the addition of 2% acetonitrile to the eluent (final eluent concentration 98 mM TFA, 2% acetonitrile) to make the separation 2.5 times faster without compromising resolution and column performance.

Results: System suitability was evaluated as described in the USP Gentamicin Sulfate monograph using chromatograms of a system suitability standard and 10 µg/mL sisomicin sulfate. The five congeners (C1, C1a, C2, C2a, and C2b) and sisomicin were well separated. Sisomicin is detected with good sensitivity. The separation, linearity, reproducibility, and sensitivity were found to meet or exceed the current USP/EP Gentamicin Sulfate monograph performance requirements. The separation and sensitivity of both modified methods were found to meet or exceed the current USP/EP Gentamicin Sulfate monograph performance requirements.

INTRODUCTION

Gentamicin is a broad-spectrum, water-soluble antibiotic belonging to the group of aminoglycoside antibiotics. The ratio of major components of gentamicin C and related substances needs to be routinely investigated and controlled in commercial products. The analysis of gentamicin sulfate based on an ion-pairing HPLC-PAD method is described in the USP and EP. The Dionex IonPac AmG-3µm column contains a silica-based packing material used for reversed-phase chromatography but is packed in a PEEK column body rather than stainless steel. This column was specifically designed for ion-pairing reversed-phase separations of aminoglycoside antibiotics and we used it to execute the methods in the USP and EP Gentamicin Sulfate monographs. Here, we apply a four-potential waveform to detect gentamicin components, rather than the three-potential waveform reported in the USP Gentamicin Sulfate monograph for the Content of Gentamicins test. Compared to the three-potential waveform, the four-potential waveform minimizes electrode wear and dramatically improves long-term peak area reproducibility. We also did the analysis with the three-potential waveform.

MATERIALS AND METHODS

Sample Preparation

Gentamicin Sulfate reference standard and sample were purchased from Sigma-Aldrich.

Sample solution (a), 1 mg/mL. Dissolve 25 mg of sample in 25 mL of eluent. Use this sample preparation for impurity analysis.

Sample solution (b), 0.2 mg/mL. Dilute 5 mL of sample solution (a) to 25 mL with eluent. Use this sample preparation for the Content of Gentamicins analysis.

Instrument

A Dionex ICS-5000* HPIC dual system was used. Figure 1 shows a flow diagram of the system configured for carbohydrate detection using electrochemical detection.

In PAD, using the four-potential waveform, the conventional working electrode is pulsed through the different potentials at set times, completing two cycles within one second (Figure 2). This waveform is optimized to provide a clean, stable gold layer in preparation for detection of the next eluting peak.

Data Analysis

Thermo Scientific™ Chromeleon™ Chromatography Data System (CDS), version 7.2.

Figure 1. The Dionex ICS-5000* HPIC system flow diagram configured for ED detection.

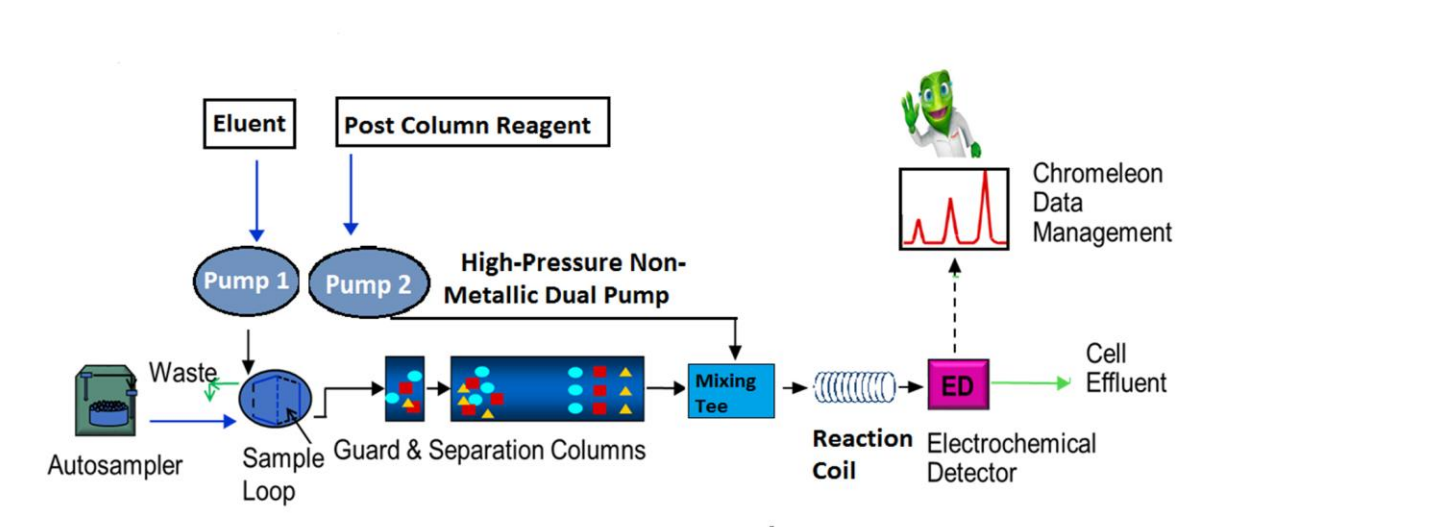
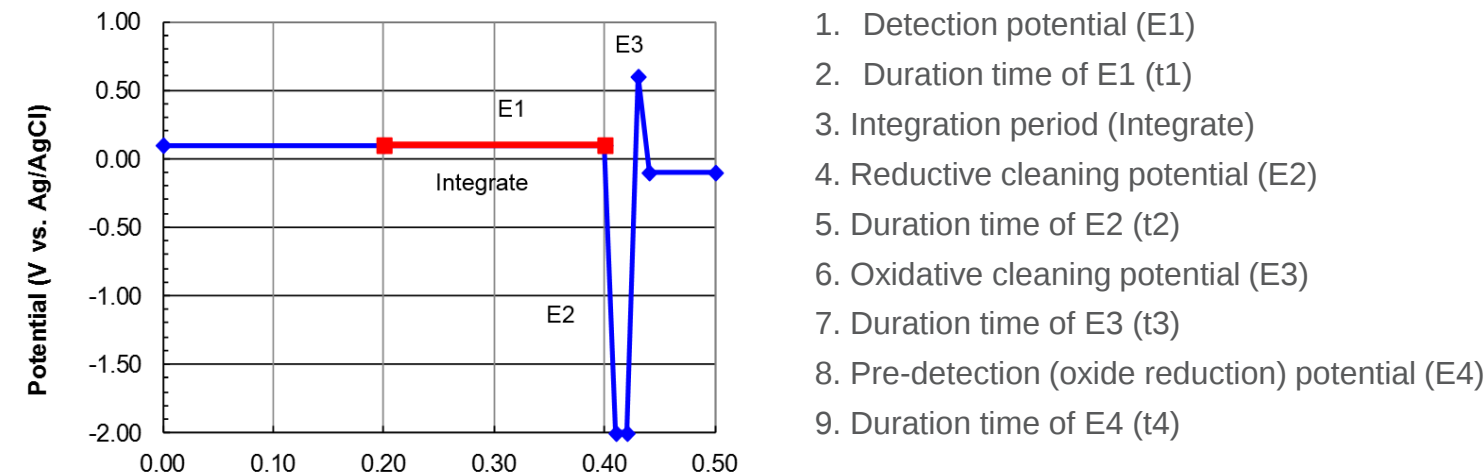


Figure 2. Four-potential carbohydrate waveform.



RESULTS

System suitability

In the USP monograph for gentamicin sulfate, resolution >1.5 between gentamicin C2 and gentamicin C2b is specified in the system suitability requirements. The EP gentamicin sulfate monograph includes two additional requirements: Signal-to-noise ratio >20 for 10 µg/mL sisomicin and resolution >1.2 between sisomicin and gentamicin C1a. The system suitability was evaluated using the chromatograms of a system suitability standard and 10 µg/mL sisomicin sulfate. Figure 3 shows this separation using a Dionex IonPac AmG-3µm C18 column set. The five congeners (C1, C1a, C2, C2a, and C2b) and sisomicin were well separated. Figure 4 shows the chromatogram of sisomicin sulfate. Sisomicin is sensitively detected. The system suitability requirements are met for all parameters (Table 1). Peak resolution between C2 and C2b is 4.2, exceeding the USP and EP requirement of 1.5. Peak resolution between sisomicin and C1a is 3.75, exceeding the EP requirement of 1.2. The signal-to-noise ratio of 10 µg/mL sisomicin sulfate is 233, easily exceeding the EP requirement of 20.

Figure 3. Separation of a system suitability standard (gentamicin 100 µg/mL + sisomicin 20 µg/mL) using a Dionex IonPac AmG-3µm C18 column.

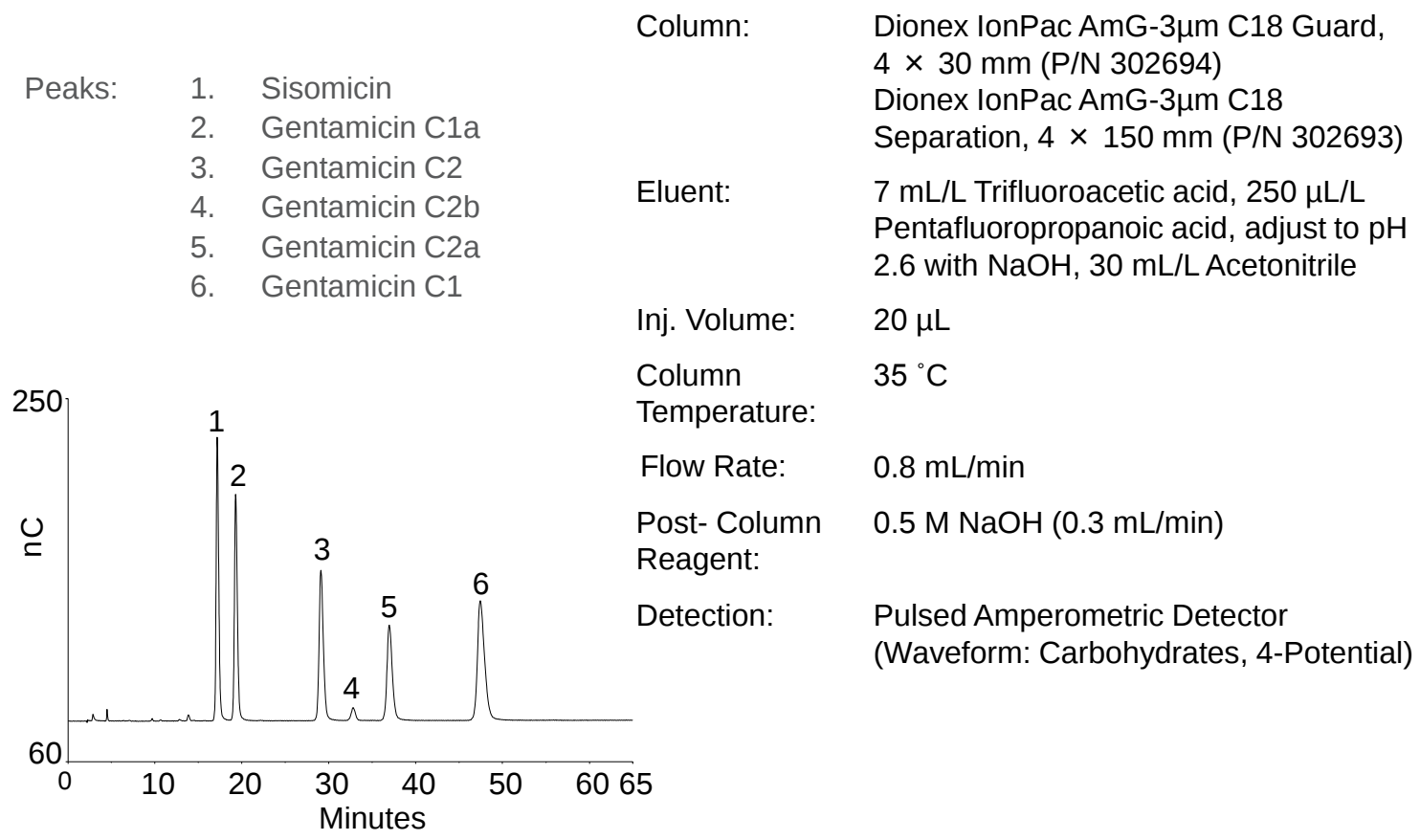


Figure 4. Sisomicin USP standard (10 µg/mL).

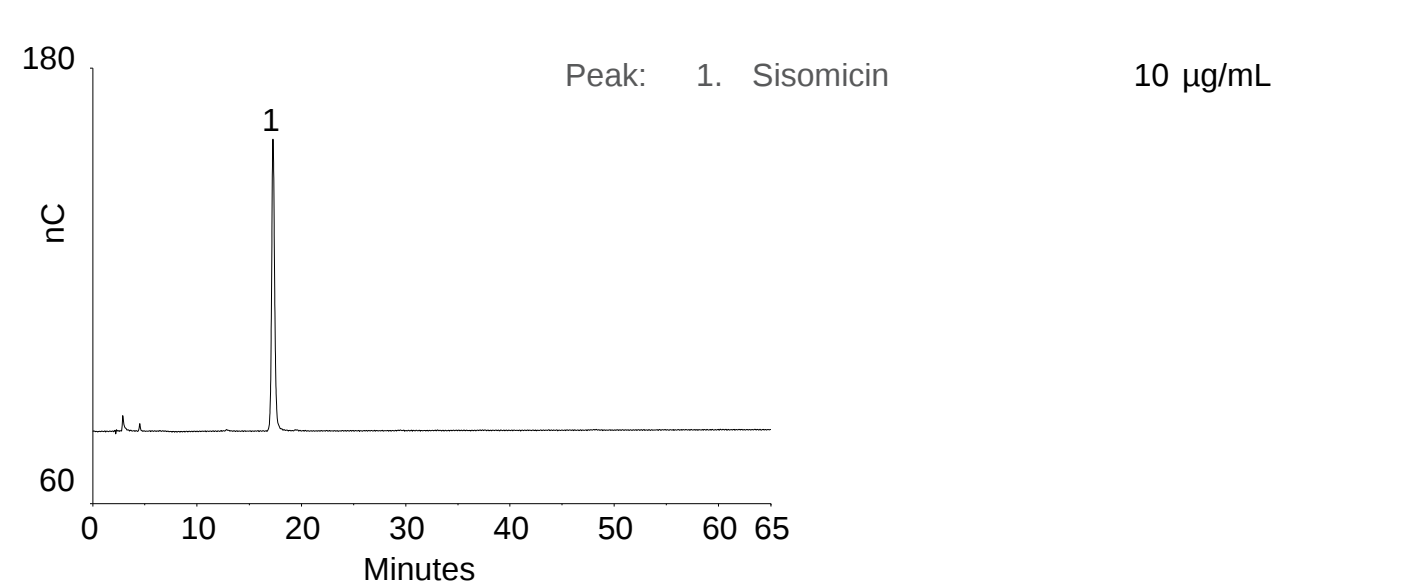


Table 1. System suitability using the four-potential waveform.

Test	EP criteria	Measured
Resolution between sisomicin and C1a	>1.2	3.75
Resolution between C2 and C2b	>1.5	4.20
Signal-to-noise ratio (sisomicin 10 µg/mL)	>20	233

Linearity

The linearity of gentamicin was investigated in the concentration range of 10–200 µg/mL (10, 25, 50, 100, 200 µg/mL). For all gentamicin species, the coefficients of determination were better than 0.997. This reveals that a sample concentration of 200 µg/mL is within the response linear range and can be used for analysis.

Method precision

Method precision was evaluated with five replicate injections of gentamicin sample (0.2 mg/mL). The relative standard deviation (RSD) for five injections of sample #2 ranged between 0.1 to 0.5%.

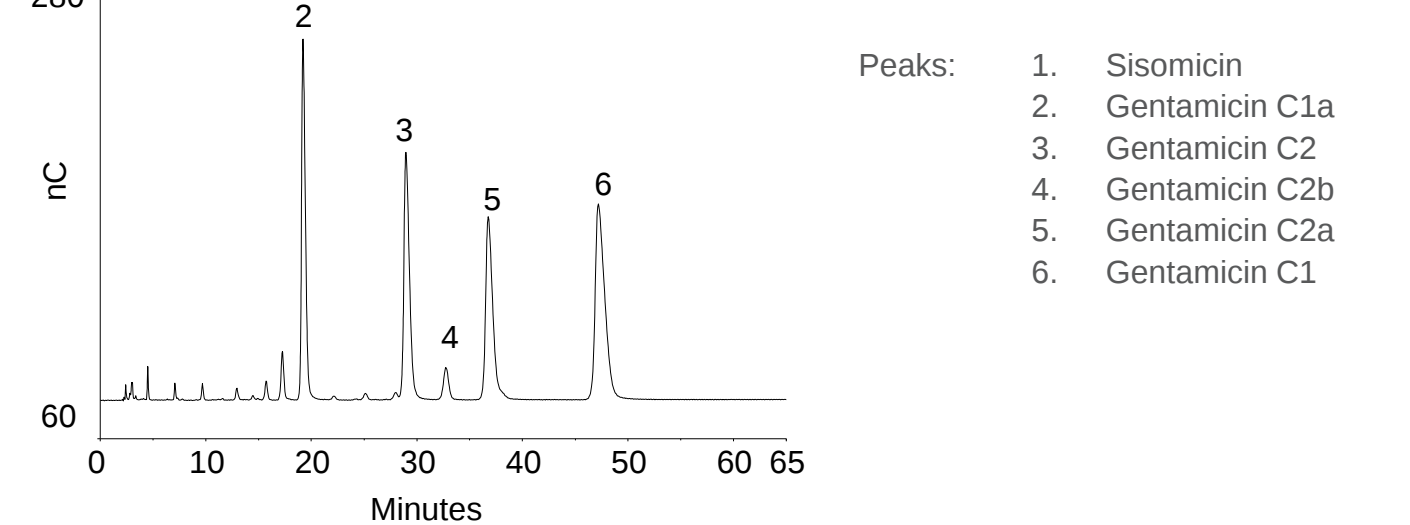
Method LOD and LOQ

The USP method for validation specifies a signal to noise (S/N) ratio of 3 for the determination of the LOD and a signal-to noise (S/N) ratio of 10 for the determination of the LOQ. The LOD and LOQ were then determined by injecting sisomicin sulfate (0.2 µg/mL). The LOD and LOQ in the sample solution were 0.173 µg/mL 0.577 µg/mL, respectively.

Sample analysis

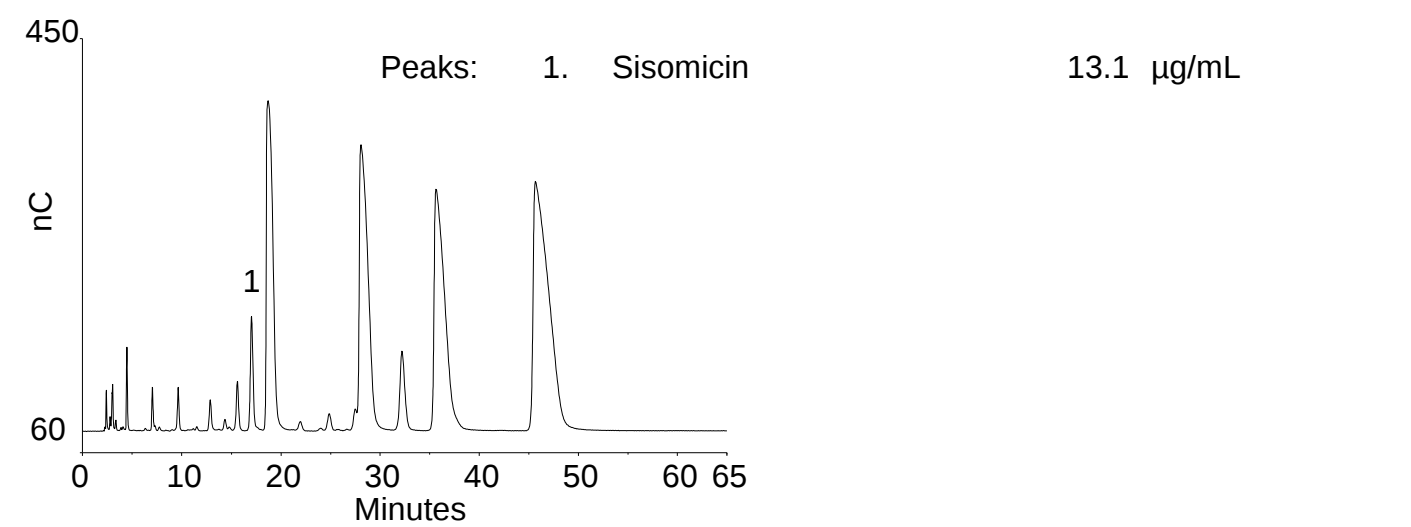
Sample solution (b) was used for content of gentamicins analysis. Figure 5 shows the separation of gentamicin sample (0.2 mg/mL). A few impurities were detected and they were separated from the five gentamicin constituents.

Figure 5. Separation of gentamicin sample (0.2 mg/mL) using a Dionex IonPac AmG-3µm C18 column.



Sample solutions (a) were used for impurities analysis. Figure 6 shows the chromatograms of samples. The five times greater concentration of these samples compared to the samples used for the Content of Gentamicins analysis allows the impurity peaks to be more easily observed.

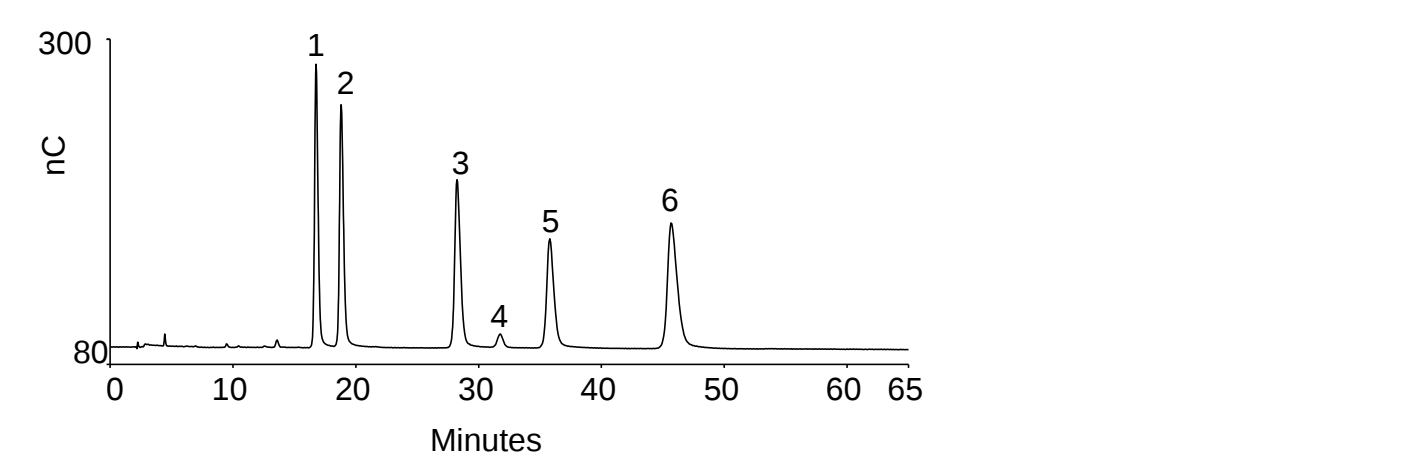
Figure 6. Separation of gentamicin sample (1 mg/mL) using a Dionex IonPac AmG-3µm C18 column.



Waveform comparison

The analysis of gentamicin was evaluated using the three-potential carbohydrate waveform that is in the USP and EP Gentamicin Sulfate monographs. Figure 7 shows the separation of a system suitability standard using the three-potential waveform. The five congeners (C1, C1a, C2, C2a, and C2b) and sisomicin were well separated. The system suitability requirements are met for all parameters.

Figure 7. Separation of a system suitability standard (gentamicin 100 µg/mL + sisomicin 20 µg/mL) using a Dionex IonPac AmG-3µm C18 column with the three-potential waveform.



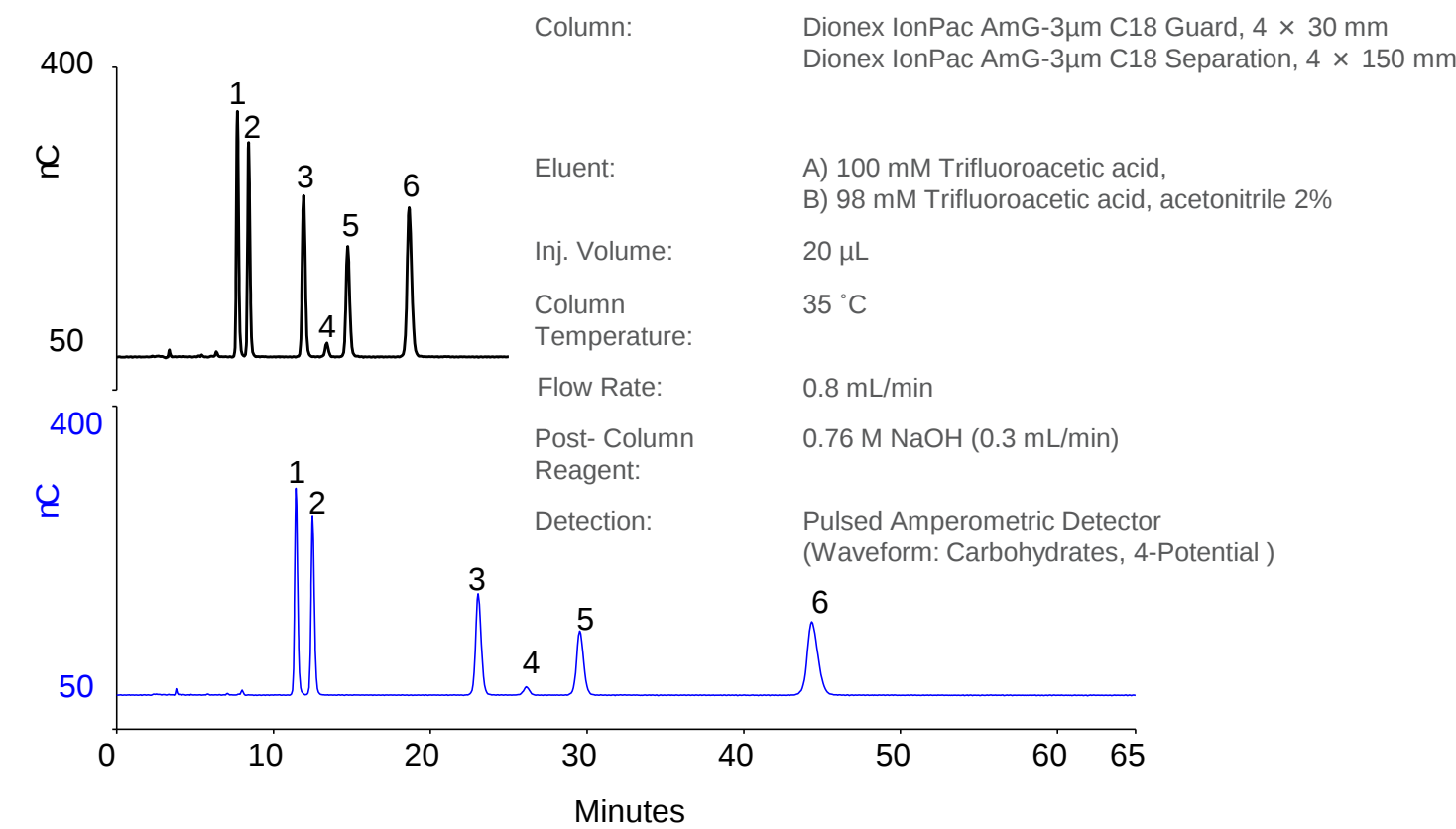
Modified methods

We modified the eluent in the USP/EP monographs in two ways, with each modification used to make a new method. Method A is a simple eluent method that uses a 100 mM TFA eluent. Method B is a fast method that involves the addition of 2% acetonitrile to the eluent (final eluent concentration 98 mM TFA, 2% acetonitrile) to make the separation 2.5 times faster without compromising resolution and column performance. The system suitability was evaluated using the chromatograms of the system suitability standard and 10 µg/mL sisomicin sulfate. Figure 8 shows this separation with a Dionex IonPac AmG-3µm C18 column set using the two methods. The five congeners (C1, C1a, C2, C2a, and C2b) and sisomicin were well separated using both methods. TFA acts as the ion-pairing agent and plays an important role in the gentamicin separation. Gentamicin separation is normally completed in 60 min when using 100 mM TFA as the eluent (Figure 8A). To accelerate the separation, 2% acetonitrile was added to the 100 mM TFA eluent (final eluent concentration 98 mM TFA, 2% acetonitrile), and this resulted in a separation that was less than 25 min (Figure 8B). Using either method the system suitability requirements are met for all parameters (Table 2).

Table 2. System suitability using modified methods.

Test	EP criteria	Measured (Method A)	Measured (Method B)
Resolution between Sisomicin and C1a	>1.2	2.85	2.63
Resolution between C2 and C2b	>1.5	4.53	3.97
Signal to noise ratio (sisomicin 10 µg/mL)	>20	248	242

Figure 8. Separation of a system suitability standard using method A and B.



CONCLUSIONS

This work demonstrated that the USP Gentamicin Sulfate monograph Content of Gentamicins method and the USP in-process revision Gentamicin Sulfate monograph method for organic impurities method could be successfully executed with a Dionex IonPac AmG-3µm C18 column using either the four- or three-potential carbohydrate waveform on a Dionex ICS-5000* HPIC system.

The separation, linearity, reproducibility, and sensitivity were found to meet or exceed the current USP/EP Gentamicin Sulfate monograph performance requirements. This method is reliable and can be used for the routine monitoring of gentamicin.

Gentamicin sulfate and related impurities can be separated with a Dionex IonPac AmG-3µm C18 column using two modified methods. Method A is a simple eluent method (100 mM TFA). Method B is a fast method that involves the addition of 2% acetonitrile to the eluent (final eluent concentration 98 mM TFA, 2% acetonitrile) to make the separation 2.5 times faster without compromising resolution and column performance. The separation and sensitivity of both methods were found to meet or exceed the current USP/EP Gentamicin Sulfate monograph performance requirements.

REFERENCES

- Thermo Scientific Application Note 72647: Determination of gentamicin and related impurities in gentamicin sulfate
- Thermo Scientific Application Update 72648: Determination of gentamicin and related impurities in gentamicin sulfate using simple eluents
- Thermo Scientific Technical Note 21: Optimal Settings for Pulsed Amperometric Detection of Carbohydrates Using the Dionex ED40 Electrochemical Detector

TRADEMARKS/LICENSING

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is not intended to encourage use of these products in any manner that might infringe the intellectual property rights of others.