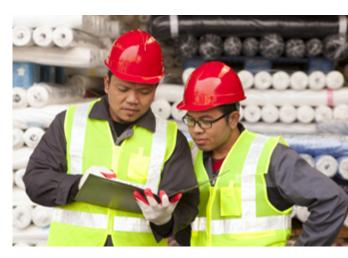
Determination of Trace Anions in Concentrated Caustics Using Single-Pass Neutralization Pretreatment and IC

Aaron Rose¹ and Paul Voelker²

¹Thermo Fisher Scientific, Bannockburn, IL, USA

²Thermo Fisher Scientific, Sunnyvale, CA, USA



Analytical Conditions

Column:	Dionex IonPac AS19 Analytical
	$(2 \times 250 \text{ mm})$
Eluent:	10 mM potassium hydroxide (KOH)
	0-12 min; 10-45 mM KOH
	12–25 min
Eluent Source:	Thermo Scientific Dionex
	EGC 500 KOH Eluent Generator
	Cartridge
Flow Rate:	0.5 mL/min
Temperature:	30 °C
Detection:	Suppressed conductivity, Dionex
	AERS 500 suppressor, 2 mm
	(P/N 082541), AutoSuppression,
	recycle mode



Introduction

Caustic soda (sodium hydroxide) formed from the electrolysis of brine is used as feedstock in pulp and paper, plastics, detergents, textiles, and other industries. Monitoring the formation of contaminants and byproducts during salt electrolysis enables purity objectives to be achieved in a safe and efficient manner.

Ion chromatography (IC) with suppressed conductivity detection is the preferred method for determining ionic species in aqueous media, especially at ppm and ppb concentrations. Yet, analysis of anions by IC in strong bases (or acids) is a challenge. Direct injections of concentrated base samples overload the column, resulting in poor chromatography and quantification, while dilution sacrifices sensitivity for the anion of interest.

Autoneutralization, introduced by Dionex (now part of the Thermo Scientific brand) in the mid '90s, eliminated the need to dilute the concentrated base, but involved a tedious two-step process requiring either a double-pass or park-and-neutralize step. With the introduction of the new Thermo Scientific[™] Dionex[™] AERS[™] 500 Anion Electrolytically Regenerated Suppressor, base neutralization can now be performed with a single pass.

Experimental

Sample Preparation Conditions

Sample:	Diaphragm caustic 1:500 dilution
	with DI water
Eluent:	DI water with Thermo Scientific [™] Dionex [™] IonPac [™]
	ATC-HC 500 Anion Trap Column $(9 \times 75 \text{ mm})$
Flow Rate:	1 mL/min
Injection Volume:	25 μL
Temperature:	30 °C
Neutralization:	Dionex AERS 500 suppressor, 4 mm
	(P/N 082540)
Concentrator	

Column: Dionex IonPac UTAC-XLP1

Results

Figure 1 depicts the two-valve configuration for single-pass neutralization with either the Thermo Scientific[™] Dionex[™] ICS-5000⁺ HPIC[™] system, capable of high-pressure IC, or the Dionex ICS-2100 IC system.

• Valve 1 employs external water regeneration for the Dionex AERS 500 suppressor by pump (Thermo Scientific Dionex AXP Auxiliary Pump for a Dionex ICS-2100 IC system or Thermo Scientific Dionex DP1 Dual Pump for a Dionex ICS-5000⁺ HPIC system) with a Thermo Scientific Dionex ATC-HC trap column and valve switch that pushes a 25 μ L loop of caustic sample through a 4 mm Dionex AERS 500 suppressor (the neutralizer mode) and then on to Valve 2.

• Valve 2 employs a KOH eluent and valve switch to elute the neutralized sample off the Dionex IonPac UTAC-XLP1 column onto a Dionex IonPac AS19 column and through a 2 mm Dionex AERS suppressor run in recycle mode.

New Optimized Single Pass Neutralization

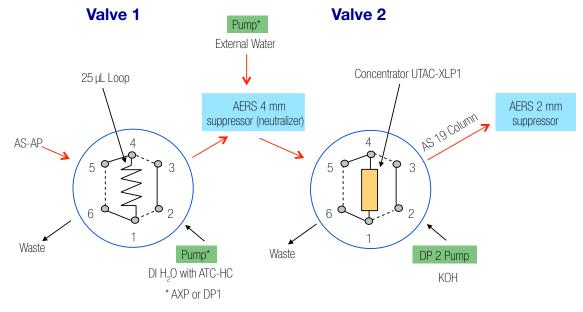


Figure 1. Single-pass pretreatment and IC analysis configuration.

Traditional Double Pass Neutralization

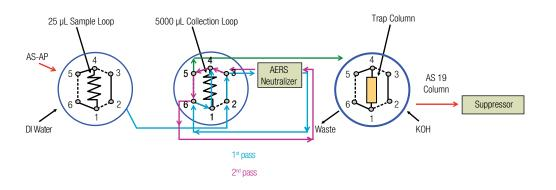


Figure 2. Double-pass neutralization and IC analysis configuration.

Figure 2 depicts the now outdated configuration for double-pass neutralization using either the Dionex ICS-5000⁺ HPIC system or the Dionex ICS-2100 IC system.

- Here the caustic sample is first loaded into a 25 μL sample loop. The sample is then switched in line and the concentrated base from the sample loop is directed to the Thermo Scientific[™] Dionex[™] ASRN[™] Anion Self-Regenerating Neutralizer. There the sample is partially neutralized, then transferred to the 5000 μL loop.
- The recycle valve is then actuated to direct the sample through the Dionex ASRN neutralizer a second time to complete the neutralization. Following the doublepass neutralization process, the sample anions can be focused onto a concentrator column before eluting them onto an analytical column.

Conclusion

The single-pass neutralization pretreatment with the new Dionex AERS 500 suppressor (operated in neutralizer mode) provides a more efficient alternative to the double-pass neutralization method for the analysis of anions in concentrated caustics. Results from the single-pass method with IC satisfy purity objectives in chlor-alkali processes by measuring trace levels of chlorate, chloride, and sulfate in a faster, more effective manner (Figure 3).

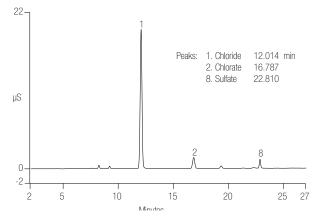


Figure 3. Single-pass neutralization and IC analysis.

www.thermoscientific.com

©2014 Thermo Fisher Scientific Inc. All rights reserved. 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 representative for details.

Africa +43 1 333 50 34 0 Australia +61 3 9757 4300 Austria +43 810 282 206 Belgium +32 53 73 42 41 Canada +1 800 530 8447 China 800 810 5118 (tree call domestic) 400 650 5118 FL90024-EN 0115S $\begin{array}{l} \textbf{Denmark} & +45\ 70\ 23\ 62\ 60\\ \textbf{Europe-Other} & +43\ 1\ 333\ 50\ 34\ 0\\ \textbf{Finland} & +358\ 9\ 3291\ 0200\\ \textbf{France} & +33\ 1\ 60\ 92\ 48\ 00\\ \textbf{Germany} & +49\ 6103\ 408\ 1014\\ \textbf{India} & +91\ 22\ 6742\ 9494\\ \textbf{Italy} & +39\ 02\ 950\ 591\\ \end{array}$

Japan +81 45 453 9100 Korea +82 2 3420 8600 Latin America +1 561 688 8700 Middle East +43 1 333 50 34 0 Netherlands +31 76 579 55 55 New Zealand +64 9 980 6700 Norway +46 8 556 468 00



Russia/CIS +43 1 333 50 34 0 Singapore +65 6289 1190 Spain +34 914 845 965 Sweden +46 8 556 468 00 Switzerland +41 61 716 77 00 UK +44 1442 233555 USA +1 800 532 4752



A Thermo Fisher Scientific Brand