

Application Note 179

LC/MS CHROMASOLV® Solvents

In traditional liquid chromatography using ultraviolet (LC/UV) detection, certain criteria are placed on the mobile phase solvents and modifiers. For example the U.S. Pharmacopoeia (USP) defines "HPLC-Grade" water by setting maximum absorbance criteria at various wavelengths as well as limiting residue on evaporation.

In LC/MS the criteria for selection of solvents and modifier reagents are significantly different than in LC/UV. Phosphate buffers that are commonplace in LC/UV exhibit poor volatility and suppress ionization in LC/MS and thus are avoided. In much the same way, metal content of solvents, which do not significantly affect most LC/UV analyses, must be examined closely for use in LC/MS.

This report discusses the positive effect of controlling sodium and potassium metals in LC/MS solvents.

Investigation of Cluster Formation in Presence of Alkali Metal Ions

In LC/MS the formation of pseudomolecular ions $[M+H]^+$ is highly desired for ease in structural interpretation of the data. When significant concentrations of sodium and potassium ions are present in the mobile phase solvents, adducts are often formed. The formation of adducts can complicate data interpretation and adversely affect sensitivity. Human gastrin (MW=2097) is used as a model peptide for this series of LC/MS studies. The normal ionization product using electrospray ionization (ESI) is the protonated, doubly-charged molecular ion with $m/z = 1049.8 [M+2H]^{2+}$.

When human gastrin is dissolved in water (0.2% formic acid) with a very low amount of sodium and potassium ions (<0.1ppm), few metal ion clusters appear, and the

Table 1. Specifications for LC/MS CHROMASOLV Solvents

LC/MS Solvent:	Water	Acetonitrile	Methanol	2-Propanol	Ethyl Acetate
Specifications					
Assay (GC) (min)	-	99.9%	99.9%	99.9%	99.7%
Fluorescence at 254nm (max)	1 ppb	0.5 ppb	1 ppb	1 ppb	-
Fluorescence at 365nm (max)	1 ppb	0.5 ppb	1 ppb	1 ppb	-
Chloride (Cl) (max)	0.000001%	-	-	-	-
Fluoride (F) (max)	0.000001%	-	-	-	-
Nitrate (NO ₃) (max)	0.00001%	-	-	-	-
Sulfate (SO ₄) (max)	0.00001%	-	-	-	-
Free acid (max)	-	0.001%	0.001%	0.001%	-
Free alkali (as NH ₃) (max)	-	0.0002%	0.0005%	0.0005%	0.0005%
Non-volatile matter (max)	0.001 %	0.0002%	0.0005%	0.0005%	0.0005%
Water (Karl Fischer) (max)	-	0.01%	0.02%	0.05%	0.03%
Transmittance at 200nm (min)	95%	95%	-	-	-
Transmittance at 230nm (min)	99%	99%	75%	75%	-
Transmittance at 260nm (min)	-	-	98%	98%	50%
HPLC gradient (254nm) (max)	1 mAU	0.2mAU	2 mAU	2 mAU	-
Silver (Ag) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	-
Aluminum (Al) (max)	0.5 ppm	0.5 ppm	0.5 ppm	0.5 ppm	-
Barium (Ba) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	-
Calcium (Ca) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm
Cadmium (Cd) (max)	0.05 ppm	0.05 ppm	0.05 ppm	0.05 ppm	-
Cobalt (Co) (max)	0.02 ppm	0.02 ppm	0.02 ppm	0.02 ppm	-
Chromium (Cr) (max)	0.02 ppm	0.02 ppm	0.02 ppm	0.02 ppm	-
Copper (Cu) (max)	0.02 ppm	0.02 ppm	0.01 ppm	0.02 ppm	-
Iron (Fe) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	-
Potassium (K) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm
Magnesium (Mg) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm
Manganese (Mn) (max)	0.02 ppm	0.02 ppm	0.01 ppm	0.02 ppm	-
Sodium (Na) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm
Nickel (Ni) (max)	0.02 ppm	0.02 ppm	0.02 ppm	0.02 ppm	-
Lead (Pb) (max)	0.1 ppm	0.1 ppm	0.02 ppm	0.1 ppm	-
Tin (Sn) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	-
Zinc (Zn) (max)	0.1 ppm	0.1 ppm	0.1 ppm	0.1 ppm	-
Particle test	+	+	+	-	-
LC/MS suitability test	+	+	+	+	+

pseudomolecular ion mass response can easily be determined (Figure A). The observed clusters $[M+H+Na]^{2+}$ and $[M+H+K]^{2+}$ are in relatively low abundance.

When the content of sodium and potassium is higher (10ppm), the adducts formed with these metal ions become dominant (Figure B). The result is a complex set of responses due to the adduct formations possible. In this case responses at m/z 1049.9, 1060.9, 1068.9 (not labeled), 1072.1 and 1080.1 correspond to $[M+2H]^{2+}$, $[M+H+Na]^{2+}$, $[M+H+K]^{2+}$, $[M+2Na]^{2+}$ and $[M+Na+K]^{2+}$ adducts, respectively.

Conclusions

In this study we have shown that the alkali metal content of solvents used in LC/MS has a significant impact on the results. High levels of sodium and potassium, commonly found in traditional LC/UV solvents can both complicate mass spectral interpretation and decrease sensitivity. Such artifacts make spectral interpretation significantly more challenging. In addition, the relative response of the pseudomolecular ion is decreased 10-fold when the alkali metal concentrations are high.

Ordering Information

Other solvents and sizes are available. Please call or visit our website.

Description	Size	Cat.No.
Water, LC/MS	1L	39253
Acetonitrile, LC/MS	1L and 2.5L	34967
Methanol, LC/MS	1L and 2.5L	34966
2-Propanol, LC/MS	1L and 2.5L	34965
Ethyl Acetate, LC/MS	1L and 2.5L	34972
Gastrin 1 Human	250µg, 0.5mg, and 1mg	G 9020

Trademark

CHROMASOLV – Sigma-Aldrich Co.

For expert answers to your questions, contact our Technical Service Department:

Phone **800-359-3041, 814-359-3041**

Fax **800-359-3044, 814-359-5468**

E-mail **techservice@sial.com**

To download Supelco's free technical literature visit us at

sigma-aldrich.com/supelco-literature

Figure A. Mass Spectrum of Human Gastrin at Low ppm Alkali Metal Concentrations

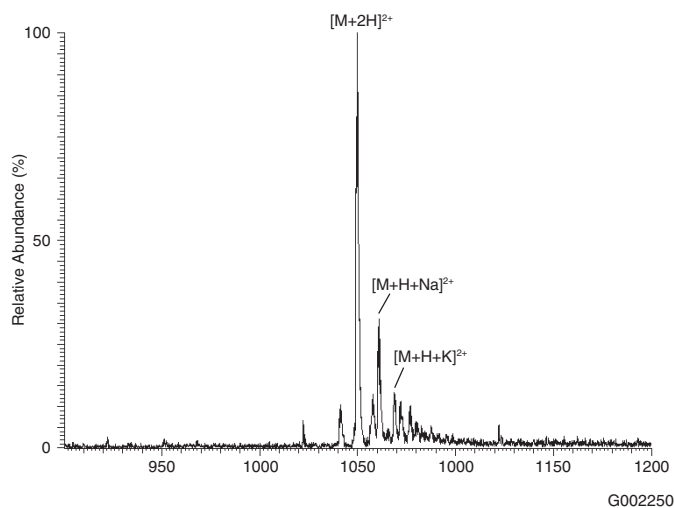
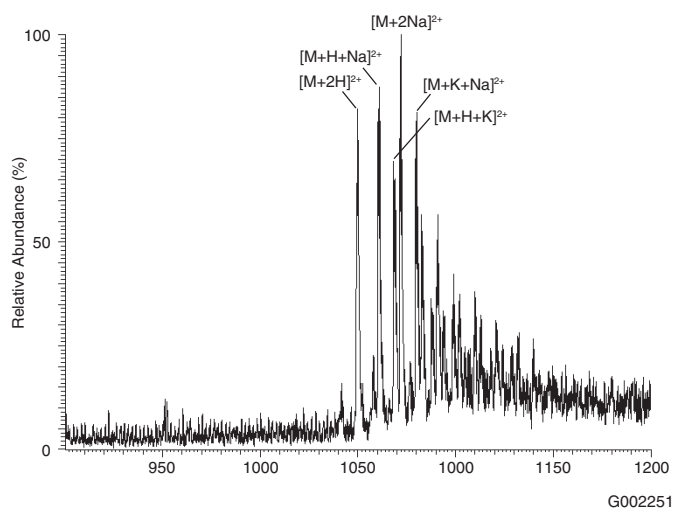


Figure B. Mass Spectrum of Human Gastrin at High ppm Alkali Metal Concentrations



Riedel-deHaën®

sigma-aldrich.com/supelco

Order/Customer Service 800-247-6628, 800-325-3010 • Fax 800-325-5052 • E-mail supelco@sial.com

Technical Service 800-359-3041, 814-359-3041 • Fax 800-359-3044, 814-359-5468 • E-mail techservice@sial.com

SUPELCO • 595 North Harrison Road, Bellefonte, PA 16823-0048 • 814-359-3441

We are committed to the success of our Customers, Employees and Shareholders through leadership in Life Science, High Technology and Service.

The SIGMA-ALDRICH Family  SIGMA  ALDRICH  Fluka  Riedel-deHaën  SUPELCO

© 2004 Sigma-Aldrich Co. Printed in USA. Supelco brand products are sold through Sigma-Aldrich, Inc. Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.



T304179
GND

ISO 9001
REGISTERED