

# HGA Graphite Furnace AA – Automatic Matrix Modification for Improved Analytical Quality and Sample Throughput

## Atomic Absorption



The matrix modification technique is a very important feature in the concept of interference-free trace metal determinations. Applying this technique, the chemical forms, and thereby the physical properties, of the element under study and/or the matrix can be changed by adding a suitable reagent in excess to the sample and standard reference solutions.

### Matrix modification

- Decreases the volatility of the analyte element and prevents its loss during thermal pretreatment. This also allows application of higher pretreatment temperatures for better matrix removal.
- Increases the volatility of matrix components and promotes their removal before atomization. PerkinElmer furnace autosamplers feature automatic matrix modification with the ability to apply one or more matrix modifiers sequentially or simultaneously.

The following tables provide detailed information on how to prepare the most common matrix modification solutions.

**Table 1. HGA Graphite Furnace: Matrix Modifiers for Routine Applications.**

Modifier	Major Application	Absolute Mass Required	Stock Reagents Required	Working Solutions (for a 5- $\mu$ L modifier addition)	
				Preparation	Concentration
Pd + Mg(NO <sub>3</sub> ) <sub>2</sub>	Ag, As, Au, Bi,	15 $\mu$ g Pd	1% (10 g/L) Pd*	Dilute 3 mL of Pd stock solution and 0.3 mL of Mg stock solution to 10 mL with 18 M $\Omega$ deionized water.	0.3% (3 g/L) Pd
	Cd, Cu, Ga, Ge, Hg, In, Sb, Se, Sn, Te, Tl <sup>a</sup>	+ 10 $\mu$ g Mg(NO <sub>3</sub> ) <sub>2</sub>	+ 1% (10 g/L) Mg **		+ 0.2% (2 g/L) Mg(NO <sub>3</sub> ) <sub>2</sub>
Mg(NO <sub>3</sub> ) <sub>2</sub>	Al, Be, Co, Cr Fe, Mn, V, Zn	50 $\mu$ g Mg(NO <sub>3</sub> ) <sub>2</sub>	1% (10 g/L) Mg**	Dilute 1.7 mL of Mg stock solution to 10 mL with 18 M $\Omega$ deionized water.	1.0% (10 g/L) Mg(NO <sub>3</sub> ) <sub>2</sub>

**Table 1. HGA Graphite Furnace: Matrix Modifiers for Routine Applications, cont.**

Modifier	Major Application	Absolute Mass Required	Stock Reagents Required	Working Solutions (for a 5- $\mu$ L modifier addition)	
				Preparation	Concentration
NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>	Pb for D <sub>2</sub> background correction	200 $\mu$ g NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>	Use either: <b>A:</b> 10% (100 g/L) NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> liquid***  or <b>B:</b> solid ultrapure NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>  <b>Caution:</b> do not use (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	<b>A:</b> Dilute 4.0 mL of NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> stock solution and 10 mL with 18 M $\Omega$ deionized water:  or <b>B:</b> Dissolve 0.4 g NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> in 18 M $\Omega$ deionized water and make up to 10 mL.	4.0% (40 g/L) NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub>
NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> + Mg(NO <sub>3</sub> ) <sub>2</sub>	Pb, Cd for Zeeman background correction	For Zeeman technique, use 200 $\mu$ g NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> together with 10 $\mu$ g Mg(NO <sub>3</sub> ) <sub>2</sub>	Use either: <b>A:</b> 10% (100 g/L) NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> liquid*** + 1% (10 g/L) Mg**  or <b>B:</b> solid ultrapure NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> + 1% (10 g/L) Mg**  <b>Caution:</b> do not use (NH <sub>4</sub> ) <sub>2</sub> HPO <sub>4</sub>	<b>A:</b> Dilute 4.0 mL of NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> stock solution and 0.3 mL of Mg stock solution to 10 mL with 18 M $\Omega$ deionized water.  or <b>B:</b> Dissolve 0.4 g NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> in 18 M $\Omega$ deionized water and add 0.3 mL of Mg stock solution and make up to 10 mL.	4.0% (40 g/L) NH <sub>4</sub> H <sub>2</sub> PO <sub>4</sub> + 0.2% (2 g/L) Mg(NO <sub>3</sub> ) <sub>2</sub>

<sup>a</sup>TI = with Zeeman background correction only.

**Table 2. HGA Graphite Furnace: Matrix Modifiers for Special Applications.**

Modifier	Major Application	Absolute Mass Required	Stock Reagents Required	Working Solutions (for a 5- $\mu$ L modifier addition)	
				Preparation	Concentration
Pd + Ca(NO <sub>3</sub> ) <sub>2</sub>	P	20 $\mu$ g Pd + 5 $\mu$ g Ca(NO <sub>3</sub> ) <sub>2</sub>	1% (10 g/L) Pd* + Ca (1 g/L)****	Dilute 4 mL of Pd stock solution and 2 mL Ca(NO <sub>3</sub> ) <sub>2</sub> to 10 mL with 18 M $\Omega$ deionized water	0.4% (4 g/L) Pd + 0.1% (1 g/L) Ca(NO <sub>3</sub> ) <sub>2</sub>
Use either: <b>A:</b> 95% Ar + 5% H <sub>2</sub> matrices or <b>B:</b> HNO <sub>3</sub>	Halide (e.g., NaCl)	Used as an alternate internal gas during dry and pyrolysis steps  NA	None  Ultrapure conc. HNO <sub>3</sub>	To 80 mL 18 M $\Omega$ deionized water, add 20 mL conc. HNO <sub>3</sub>  For a 10- $\mu$ L modifier addition	20% (v/v) HNO <sub>3</sub>
CH <sub>3</sub> OH HF HNO <sub>3</sub>	Boric acid matrix	10 $\mu$ L CH <sub>3</sub> OH + 0.2 $\mu$ L conc. HF + 0.2 $\mu$ L conc. HNO <sub>3</sub>	CH <sub>3</sub> OH + conc. HF + conc. HNO <sub>3</sub>	10 mL CH <sub>3</sub> OH + 200 $\mu$ L conc. HF + 200 $\mu$ L conc. HNO <sub>3</sub> and mix all	

\* Part No. B0190635

\*\* Part No. B0190634

Note: 1% Mg corresponds to 6% Mg(NO<sub>3</sub>)<sub>2</sub> or 10.5% Mg(NO<sub>3</sub>)<sub>2</sub> • 6H<sub>2</sub>O

1  $\mu$ g Mg(NO<sub>3</sub>)<sub>2</sub> corresponds to 0.17  $\mu$ g Mg or 1.75  $\mu$ g Mg(NO<sub>3</sub>)<sub>2</sub> • 6H<sub>2</sub>O

\*\*\* Part No. N9303445

\*\*\*\* Part No. N9300108

Note: 1% Ca corresponds to 4.09% Ca(NO<sub>3</sub>)<sub>2</sub> or 5.89% Ca(NO<sub>3</sub>)<sub>2</sub> • 4H<sub>2</sub>O

1  $\mu$ g Ca(NO<sub>3</sub>)<sub>2</sub> corresponds to 0.24  $\mu$ g Ca or 1.44  $\mu$ g Ca(NO<sub>3</sub>)<sub>2</sub> • 4H<sub>2</sub>O