



Paraben Standard for Gradient Module Testing Instruction Manual

Catalog Numbers

125-0083

125-0124

125-0131

343-0012

BIO-RAD

For Initial System Test

The Paraben Standard, for gradient module testing, is a lyophilized mixture of methyl-, ethyl-, and propyl parabens which absorbs light at 280 nm and separates well on a reversed phase guard cartridge. The standard is used to test the capabilities of Bio-Rad's gradient HPLC systems. It is supplied as a set of six vials containing lyophilized standard, with a total paraben content of approximately 3 mg per vial.

Instructions

Rehydrate the standard in one of the vials by adding 1.0 ml of water and 1.0 ml of 70% isopropanol. Swirl gently to dissolve the parabens, and allow the vial to stand 2-3 minutes. Swirl again; filter through a micro Prep-Disc® filter (catalog number 343-0012) to remove fine particulates before injection.

Elution Conditions

Use the Bio-Sil® ODS-5S guard cartridge in a Standard Cartridge Holder. These are included with each new gradient module.

Form a gradient by mixing two solvents: Solvent A is water and solvent B is 70% isopropanol in water. Use a 7 minute linear gradient from 20 to 50% B at a 1.5 ml/min flow rate. Inject 20 µl of standard, and detect at 280 nm. A chromatogram similar to that in Figure 1 should be obtained.

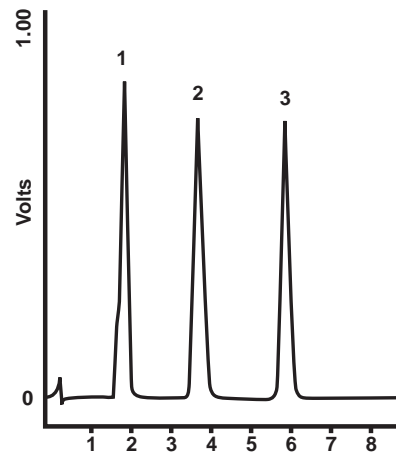


Fig. 1. Example of a test chromatogram. 1 = methyl paraben, 2 = ethyl paraben, 3 = propyl paraben.

Ordering Information

Catalog Number	Product Description
125-0083	Paraben Standard
125-0124	Bio-Sil ODS-5S Cartridge, 2
125-0131	Standard Cartridge Holder
343-0012	Micro Prep-Disc Filters, 0.45 mm, 50