

LSC in Practice

Safer Cocktail Selection for Use with Potassium Chloride

Problem

One of PerkinElmer's field representatives requested help for a researcher currently using PerkinElmer cocktails that contained toluene and xylene. The researcher hoped to "skip" two generations of cocktails and move up to PerkinElmer's newest, ULTIMA Gold™, line of cocktail. The researcher's primary concern was compatibility of the cocktail and sample.

The researcher's sample consisted of 0.4 M potassium chloride.

Discussion

We evaluated the sample loading capacities for several members of PerkinElmer's ULTIMA Gold family of cocktails using 0.4 M potassium chloride at 20°C.

The results of our testing are tabulated below:

Recommendations

The results from Table 1 show that PerkinElmer's ULTIMA Gold XR and ULTIMA Gold LLT are both highly suited for this researcher's application.

If the researcher is using an 8 mL vial, we recommend using 1.5 mL of 0.4 M KCl with 5.0 mL of ULTIMA Gold XR or ULTIMA Gold LLT. If a 6 mL vial is being used, then 1.5 mL of 0.4 M KCl may be added to 4.0 mL of cocktail.

Cocktail	Maximum Capacity (mL/10.0 mL cocktail)	Maximum Capacity (mL/5.0 mL cocktail)
ULTIMA Gold (part number 6013329)	1.5 mL 0.4 M KCl	0.75 mL 0.4 M KCl
ULTIMA Gold MV (part number 6013159)	1.5 mL 0.4 M KCl	0.75 mL 0.4 M KCl
ULTIMA Gold XR (part number 6013119)	6.0 mL 0.4 M KCl	3.0 mL 0.4 M KCl
ULTIMA Gold LLT (part number 6013377)	6.0 mL 0.4 M KCl	3.0 mL 0.4 M KCl

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
Phone: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/lasoffices

©2007 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. ULTIMA Gold is a trademark of PerkinElmer, Inc. or its subsidiaries, in the United States and other countries. All other trademarks not owned by PerkinElmer, Inc. or its subsidiaries that are depicted herein are the property of their respective owners. PerkinElmer reserves the right to change this document at any time without notice and disclaims liability for editorial, pictorial or typographical errors.