

LSC in Practice

LSC Cocktails — Elemental Composition

Problem

A researcher wanted to know the percentages of the elements present in various commonly used cocktails to satisfy regulations concerning the disposal of cocktail waste by incineration.

Discussion

This information was calculated mathematically from the cocktail formulation and the elemental composition of the components. The data, expressed as percentages, are presented in the following table:

PerkinElmer Cocktail Part Number	Carbon C Weight	Hydrogen H	Nitrogen N	Oxygen O	Phosphorus P	Sulphur S	Sodium Na	Average Molecular
ULTIMA Gold™ 6013329	78.9%	9.6%	0.2%	9.5%	1.4%	0.2%	0.2%	254.7
ULTIMA Gold XR 6013119	73.2%	10.0%	0.2%	15.2%	0.9%	0.3%	0.2%	297.5
ULTIMA Gold AB 6013309	76.3%	9.7%	0.05%	13.8%	0.1%	N/A	N/A	293.42
ULTIMA Gold MV 6013159	78.0%	10.1%	0.2%	10.2%	1.1%	0.2%	0.2%	261.68
ULTIMA Gold LLT 6013377	76.4%	9.7%	0.05%	13.85%	0.01%	N/A	N/A	293.28
ULTIMA Gold F 6013179	90.46%	9.39%	0.07%	0.08%	N/A	N/A	N/A	211.97
Insta-Gel® 6013399	70.5%	9.8%	0.02%	19.7%	N/A	N/A	N/A	315.6

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 and Insta-Gel is a registered 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.