

# LSC in Practice

## Counting Potassium Hydroxide May Produce High Background Due to $^{40}\text{K}$

### Problem

A researcher called one of our offices to report a suspected contamination problem with one batch of PerkinElmer's Hionic-Fluor™ (PerkinElmer part number 6013319).

The researcher was using Hionic-Fluor to count  $^{14}\text{CO}_2$  that had been trapped in 2 M potassium hydroxide, which acted as an absorber in this sample collection system. After collection, by bubbling the gas through the absorber solution, 2 mL of sample were added to 10 mL of Hionic-Fluor. When counting the prepared homogeneous sample, the researcher was observing backgrounds that were over 100 CPM (0-156 keV) and stable for longer than 12 hours.

The stability of the high backgrounds indicated contamination, rather than chemiluminescence, which prompted the researcher to report the problem to a PerkinElmer Application Laboratory.

### Discussion

Upon receipt of the report, we immediately tested our retained samples of the suspected lot against an alternate lot.

Testing was conducted under the following conditions:

1. Instrument	PerkinElmer Tri-Carb® Model 1900 operated at 20 °C
2. Windows	0-156 keV and 4-156 keV
3. Vials	High performance glass vials (PerkinElmer part number 6001015)
4. Cocktails	Hionic-Fluor
5. Cocktail volume	10.0 mL
6. Sample	2.0 mL of 2 M KOH

Under the conditions listed below, we obtained backgrounds of 70 to 100 CPM for both lots of Hionic-Fluor. These backgrounds did not decay over a 12 hour period and the spectrum strongly suggested an isotopic contamination.

We repeated our comparison testing but substituted 2 M NaOH as the sample and only observed normal background levels for both lots of cocktail.

Consequently, we determined that the background problem was associated with the potassium hydroxide, due to high levels of naturally occurring radioactive  $^{40}\text{K}$ .

### Recommendation

To avoid this unexpected source of contamination, we recommended that the  $\text{CO}_2$  absorber solution be changed from potassium hydroxide to sodium hydroxide. Our prior research indicates that the same  $\text{CO}_2$  capacity will be achieved without the potential for unwanted backgrounds.

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



For a complete listing of our global offices, visit [www.perkinelmer.com/lasoffices](http://www.perkinelmer.com/lasoffices)

©2008 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. Hionic-Fluor is a trademark and Tri-Carb 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.