

# LSC in Practice

## Solubilization of Insect Larvae

### Problem

A researcher was using PerkinElmer's Soluene®-350 (PerkinElmer part number 6003038) to solubilize insect larvae. The larvae were about 1 cm long and 2 mm wide. The larvae were cut up into three or four pieces before they were solubilized. The researcher observed that the cuticle of the larvae would become clear, but would not solubilize. The researcher was concerned that the remaining artifact of the cuticle could cause physical quenching. The nuclide of interest was  $^{14}\text{C}$ .

### Discussion

The cuticle is the exoskeleton which supports the body of the larvae, and will not be solubilized by Soluene-350. To solubilize the cuticle, an acid solubilization should be performed as described in LSC *Sample Preparation by Solubilization* Application Note (007902\_31).

The researcher could try one of the acid methods recommended in the Application Note – probably the perchloric/nitric mixed acid system – to solubilize the entire larvae. However, this could lead to a potential loss of  $^{14}\text{C}$  as  $^{14}\text{CO}_2$  during such a solubilization procedure.

Fortunately, since this researcher is looking at  $^{14}\text{C}$ , and the cuticle is transparent after solubilization, there should not be any physical quench problem when counting. This is especially true since  $^{14}\text{C}$  is a high energy beta emitter.

### Recommendation

The primary concern is the potential loss of signal due to physical quenching. To verify proper counting conditions, the researcher should utilize one of the following recommendations:

1. Stay with the present procedure and confirm recovery by arranging to have a larva checked by combustion in an Oxidizer.
2. After solubilization with Soluene-350 and counting, remove the cuticle and carefully wash it. Add the cuticle to another vial, then add Hionic-Fluor™ and count to see if there is any activity originating from the cuticle.

It could be that the cuticle has not assimilated any  $^{14}\text{C}$  in this experiment.

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)

©2007 PerkinElmer, Inc. All rights reserved. The PerkinElmer logo and design are registered trademarks of PerkinElmer, Inc. Hionic-Fluor is a trademark and Soluene 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.