

## Injection of Tumor Cells

1. Tumor cells, at a prescribed cell concentration, can be injected subcutaneously (20-100 uL), intraperitoneally (200-750 uL), intravenously (50-100 uL), by an intracardiac route (50-100uL), or orthotopically (20-50 uL).

Recommendations: Use a 25g 5/8" needle for s.c. and i.p. injections of adult mice, 26 g 1/2" needle for i.v. injections, 26-27g 1/2" needle for i.c. injections, and a 30g 1/2" needle for some orthotopic injections.

2. Dilute cells in DPBS and mix well before each injection. Draw the cells into the syringe without a needle to prevent shearing. Attach the needle to the syringe, and before injecting flick or invert the syringe to ensure the cells are in suspension.

### Notes:

- a. Do not pass cells through needle more than once since this will shear cells and reduce total cell number.
- b. Be sure to resuspend cells well in the syringe barrel before each injection.
- c. It is recommended that no more than three mice be injected per syringe.
- d. Concentrating cells in a small volume helps to localize and increase the initial signal *in vivo*

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (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/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.