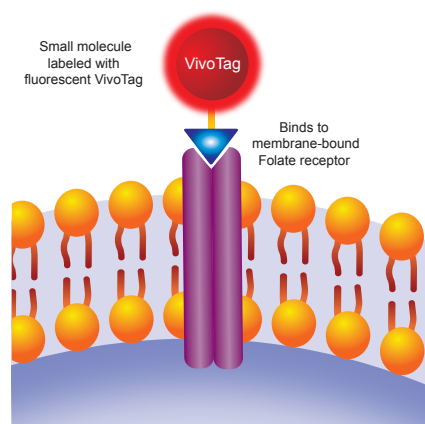


QUANTITATE TUMOR METABOLISM

FolateRSense™ 680 Targeted Fluorescent Pre-clinical Imaging Agent

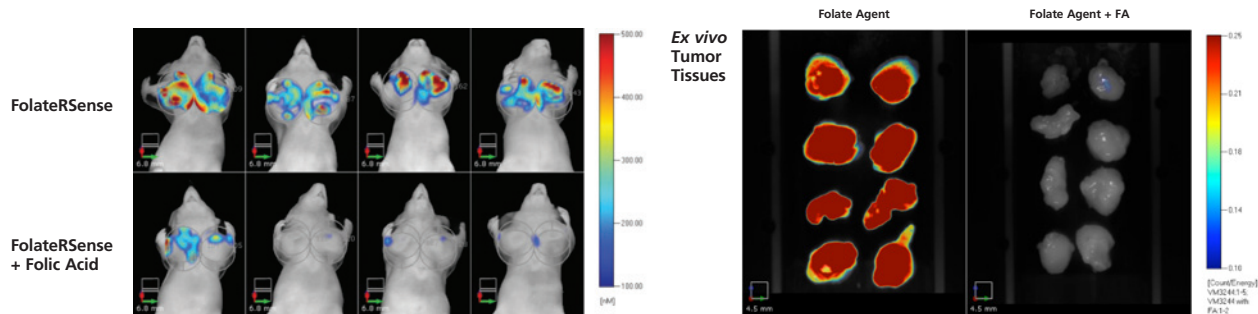
Overexpression of Folate Receptor (FR) protein is associated with tumor growth, and uptake of folic acid is a necessary part of tumor metabolism. The FolateRSense imaging agent is highly specific and sensitive in its detection of FR and can be used to closely monitor and quantitate tumor growth and metabolism.

- **Highly specific** to Folate Receptor protein, which is expressed by many growing tumors.
- Equivalent sensitivity and specificity to anti-folate receptor antibody, and **optimized for use *in vivo***.
- Quantitates folate receptor protein **even in low-expressing tumor models**.
- Rapid clearance—five minute half life—**enables longitudinal studies of tumor metabolism**.
- Excitation wavelength compatible with fluorescence microscopy: **use the same probe *in vitro*, *in vivo*, and *ex vivo***.



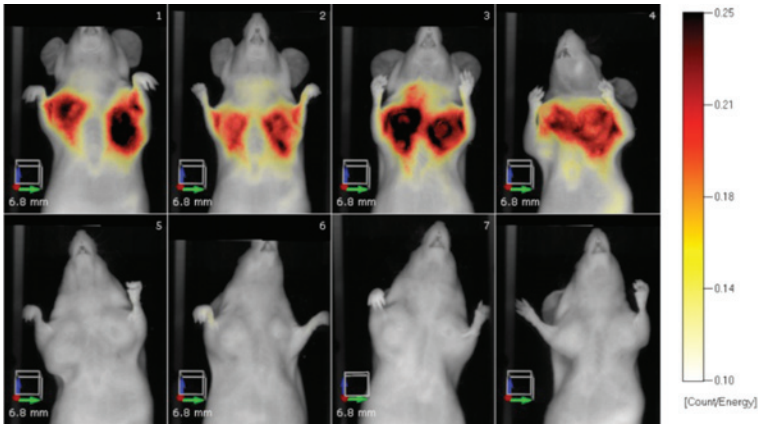
FolateRSense agent mechanism of action

Highly Specific Quantitation



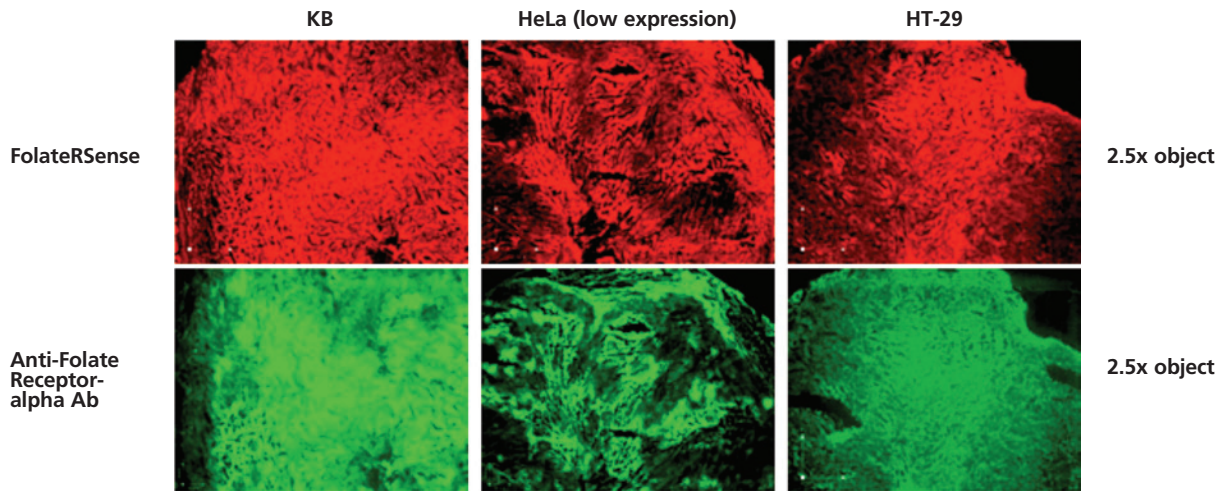
KB Tumor Xenografts: 2 mM (100x) free folic acid competition shows specific binding of FolateRSense agent, *in vivo*, in four mice.

Sensitivity for Low FR Expressing Tumors



Mice were implanted with low FR-Expressing HT-29 cells subcutaneously. Even low FRA-expressing tumors are effectively and specifically imaged with the FolateRSense. Bottom panel shows mice that were injected with Folic acid resulting in blockage of fluorescent signal.

Equivalent to Antibody Performance



Ex vivo tissue study shows FolateRSense as sensitive and specific as anti-FR antibody, even in low FR expressing HeLa cells.



The FolateRSense agent is optimized for use on PerkinElmer *in vivo* imaging Systems.

Learn more at www.perkinelmer.com/invivo

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



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

Copyright © 2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

010184_01

Printed in USA

March 2012