

Research Use Only. Not for use in diagnostic procedures.

**Bioware® Brite Cell Line 4T1-Red-FLuc**

**Product No.: BW124087**

**We recommend using this line in Nude and SCID mouse models.  
An immune response may occur in Balb/c mice.**

**Material Provided**

**Cells:** 2 x 1 mL frozen aliquots (BW124087V)

**Format:** 1.0 x 10<sup>6</sup> cells / mL in 95% FBS, 5% DMSO

DESIGNATION	4T1-Red-FLuc
Tissue	Mouse: mammary gland; adenocarcinoma
Source of Parental Line	ATCC (CRL-2539)
Gene Transfer Vehicle	Red-FLuc-Puro 3d generation lentivirus
Bioluminescence In Vitro	At least <b>15,000</b> photons/cell/sec. Exact number will vary depending on imaging and culturing conditions.
Recommended Media and FBS	RPMI 1640 ATCC Cat. No. 30-2001. Supplement the above with 10% Hyclone Fetal Bovine Serum (FBS) GE HealthCare Cat. No. SH300071
Recommended Storage Conditions	Remove frozen cells from dry ice packaging and immediately place cells at a temperature below -130° C, preferably in liquid nitrogen vapor, until ready to use.
Average Doubling Time	14 hours
Other Recommendations	When initially thawing, use T25 flask or 10cm plate. Cells should be ready to expand within 1-4 days. Antibiotics can be used in the media if desired after the initial thaw. (puromycin at 5ug/mL).  <b>Refer to Cell Culture Guidelines for more detailed instructions.</b>

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 products, visit [www.perkinelmer.com](http://www.perkinelmer.com).

Copyright ©2015, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

## The Features

### Perkin Elmer Bioware® Brite cell line models offer researchers the ability to:

- Monitor early tumor development
- Monitor tumor growth and metastases *in vivo*
- Quantify tumor burden in the whole animal
- Follow responses to therapeutic treatments non-invasively in longitudinal studies using the same cohorts of mice

### Murine Pathogen Free

All Perkin Elmer cell lines are confirmed to be pathogen free by the IMPACT Profile I (PCR) at the University of Missouri Research Animal Diagnostic and Investigative Laboratory.

### Cell Line Stability

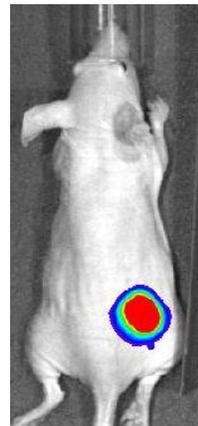
Cell may undergo genotypic changes resulting in reduced responsiveness over time in normal cell culture conditions. Genetic instability is a biological phenomenon that occurs in all stably transfected cells. Therefore, it is recommended to prepare an adequate number of frozen stock at early passages.

### Product Warranty

PerkinElmer warrants that cells will be viable upon shipment from PerkinElmer for a period of thirty days, provided they have been properly stored and handled during this period.

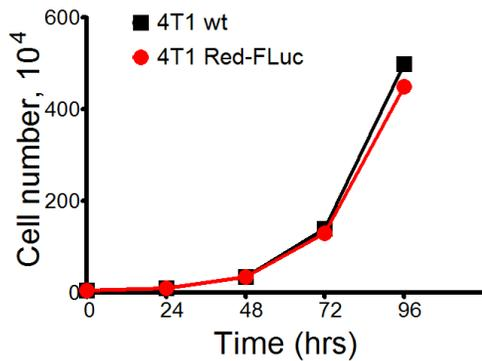
### Mouse Mammary Gland Cancer Cell Line: 4T1-Red-FLuc

4T1-Red-FLuc is a luciferase expressing cell line which was stably transfected with firefly luciferase gene from *Luciola Italica* (Red-FLuc). The cell line was established by transducing lentivirus containing Red-FLuc luciferase gene under the control of human ubiquitin C promoter. These cells will serve as a new tool to detect drug efficacy *in vitro* and *in vivo* with high sensitivity in nude and SCID mouse models.



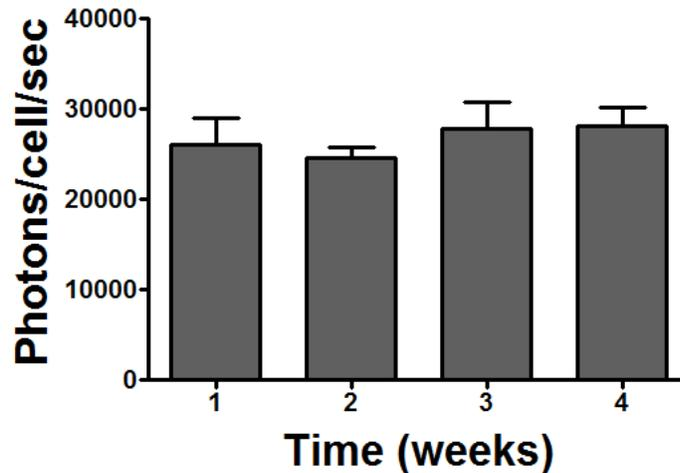
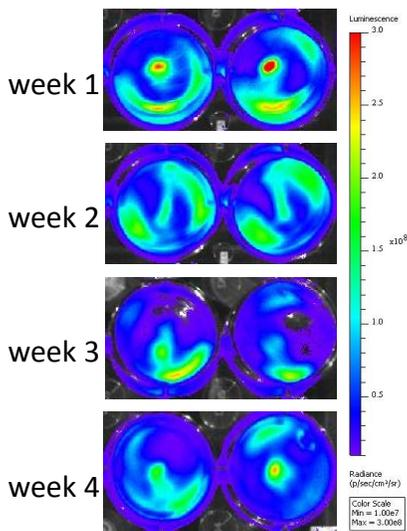
Bioluminescence image  
of 4T1 Red-FLuc  
subcutaneous tumor

## Growth Curve of 4T1-Red-FLuc Cells



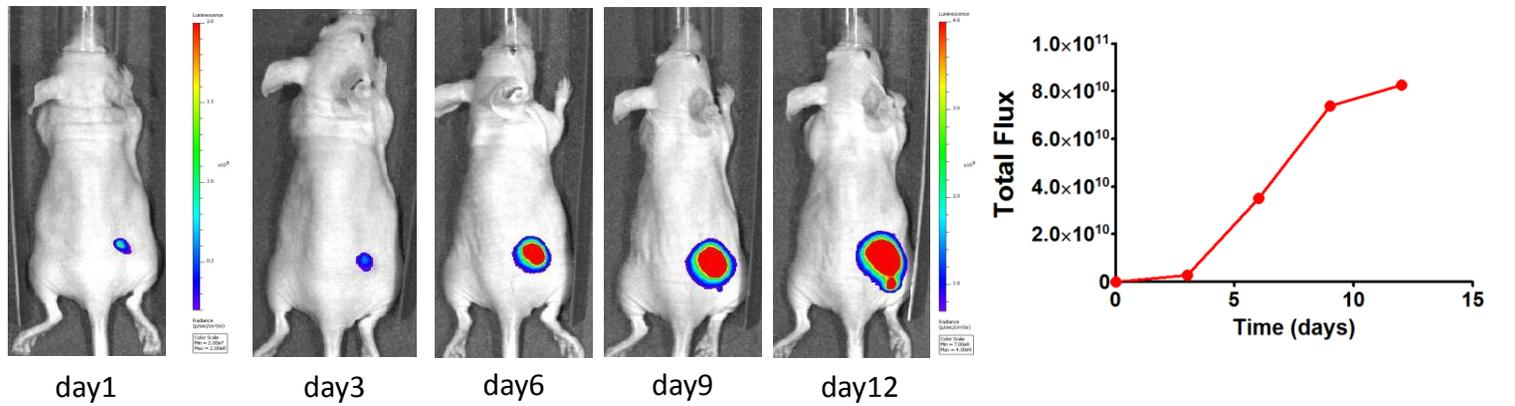
**Figure 1.**  $5 \times 10^4$  cells were plated on a 6cm plate. The total numbers of cells were counted every 24 h using a Nexcelom automatic cell counter.

## In Vitro BLI Signal Stability



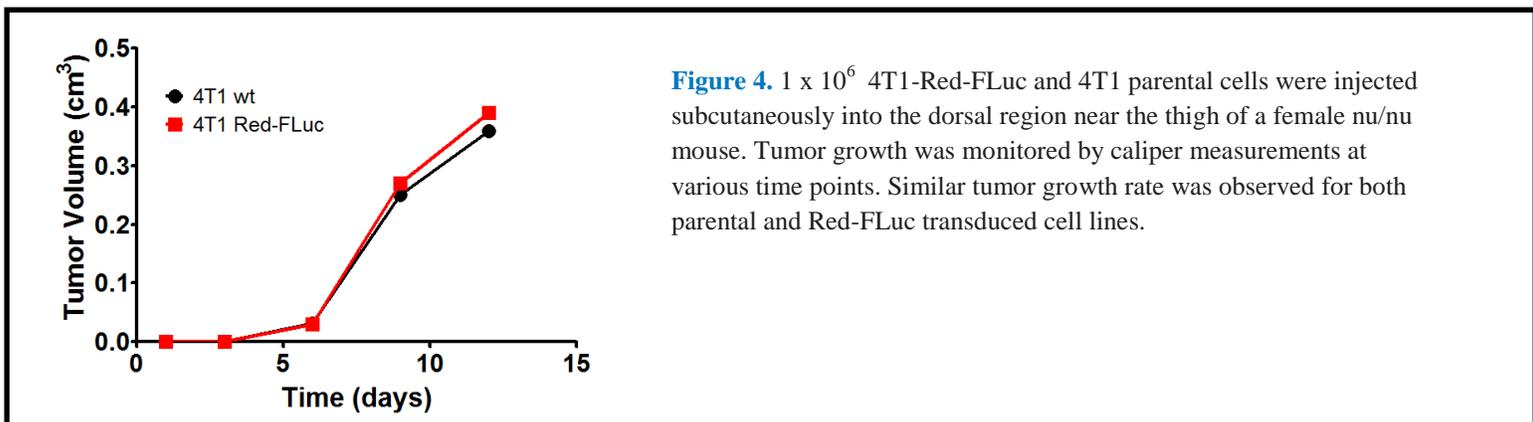
**Figure 2.**  $5 \times 10^4$  cells were plated per well in 24-well plates. Cells were incubated at 37 °C for recovery overnight and luciferase assay was performed using the PerkinElmer IVIS<sup>®</sup> SpectrumCT. Each experiment was done in quadruplicates. The cells were maintained in continuous culture over four weeks and weekly luciferase assay was performed. Bioluminescence data was analyzed using the Living Image 4.0 software.

## Subcutaneous Tumor Growth in a Nu/nu Mouse



**Figure 3.**  $1 \times 10^6$  4T1-Red-FLuc cells were injected subcutaneously into the dorsal region near the thigh of female nu/nu mouse. Tumor growth was monitored for luciferase expression using the PerkinElmer IVIS<sup>®</sup> Spectrum at various time points. Mice were imaged 10 minutes post i.p. injection of luciferin at 150mg/kg at various time points. The image above shows tumor growth from a representative mouse.

## Tumor Growth Comparison Between Wild Type and Red-FLuc Cells



**Figure 4.**  $1 \times 10^6$  4T1-Red-FLuc and 4T1 parental cells were injected subcutaneously into the dorsal region near the thigh of a female nu/nu mouse. Tumor growth was monitored by caliper measurements at various time points. Similar tumor growth rate was observed for both parental and Red-FLuc transduced cell lines.

For more information on our *in vivo* imaging agents, please visit our website: [www.perkinelmer.com/bioware](http://www.perkinelmer.com/bioware).

**This product is sold for *in vivo* animal research use only and is not intended for any diagnostic use or procedures. Excluding purchases by authorized PerkinElmer distributors, this product is sold for use by the original purchaser and is not for resale. Unless otherwise agreed to in writing by PerkinElmer pursuant to a separate written agreement, no commercial use of this product is allowed. “Commercial use” means any and all uses of this product and/or its derivatives by a party for money or other consideration and may include without limitation: (1) product manufacture; (2) providing services, information or data to another party for remuneration; and/or (3) resale of the product or its derivatives, whether or not such product or derivatives are resold for use in research. Commercial use does not include the original purchaser providing the product to its contractor solely for use on the original purchaser’s research; provided that all product materials are returned to the original purchaser and/or destroyed by the contractor upon completion of such project.**