

DD-Fluorescent Protein Reporter Systems Protocol-at-a-Glance (PT4088-2)

This protocol is provided for use with the DD-Fluorescent Protein Reporter Systems using vector-based (Cat. Nos. 631079, 631081, 631083, 631085, 631087, 631089, 632190, 632191 & 632192) or lentiviral delivery (Cat. Nos. 631748, 631751 & 631753). For a general introduction to the theory of the ProteoTuner™ and Lenti-X™ Lentiviral Expression Systems, please refer to their respective User Manuals (PT4039-1 for the ProteoTuner Systems and PT5135-1 for the Lenti-X Systems). Both manuals can be found at www.clontech.com/manuals. For the specifics of the DD-Fluorescent Protein Reporter Systems protocol, refer only to this Protocol-at-a Glance.

In brief, the experimental protocol is as follows. First, transfect your cells with the pDD-Fluorescent Protein Vector or pLVX-DD-Fluorescent Protein Vector containing the precloned NF κ B or CRE promoter, or your promoter of interest. Then, treat the cells under three conditions: without Shield1 or an inducer (the background control), with Shield1 but not the inducer (the negative control), or with Shield1 plus the inducer (the test condition). In the absence of Shield1, the reporter protein is rapidly targeted to and degraded by proteasomes, which minimizes background fluorescence from leaky promoters. When Shield1 is added, the reporter is stabilized; and when the inducer is added together with Shield1, reporter molecules expressed during promoter activation will contribute to the fluorescence signal, providing a high signal-to-noise ratio.

A. Clone your promoter of interest

(Promoterless systems only; Cat. Nos. 631748, 631751, 631753, 632190, 632191 & 632192)

1. Using either standard or In-Fusion™ cloning techniques, insert your promoter of interest into the multiple cloning site (MCS) of the vector included with your system. This promoter will now drive the expression of the fluorescent reporter (DD-AmCyan1, DD-ZsGreen1, or DD-tdTomato).
2. Transfect the resulting construct into your cells of interest by your method of choice.
3. Culture cells for 12–24 hr.

B. Study promoter activity *(all systems)*

1. Split the cells into at least three samples, depending on how many samples you would like to analyze.

NOTE: We recommend performing all experiments in duplicate or triplicate.

- **For adherent cells:** Split your cells into at least three parallel cultures. Allow the cells to adhere.

NOTE: We recommend the use of 6-well plates; however other plate formats can also be used.

- **For cell suspensions:** Distribute your cell suspension evenly into at least three wells.
2. Prepare the following culture media solutions. The total volume required depends on the number of wells/plates in your experiment.
 - **Background control** (for basal promoter activity; without Shield1): Prewarmed (37°C) culture medium.
 - **Negative control** (no promoter activation; with Shield1): Dilute the Shield1 stock solution 1:500 in prewarmed (37°C) culture medium, for a final concentration of 1 μ M.
 - **Test condition** (with promoter activation *and* Shield1): Dilute the Shield1 stock solution 1:500 in prewarmed (37°C) culture medium, for a final concentration of 1 μ M. Add the inducer or other compound of interest to this medium, at the concentration defined by your experimental protocol.



Clontech

United States/Canada
800.662.2566

Asia Pacific
+1.650.919.7300

Europe
+33.(0)1.3904.6880

Japan
+81.(0)77.543.6116

Clontech Laboratories, Inc.
A Takara Bio Company
1290 Terra Bella Ave.
Mountain View, CA 94043
Technical Support (US)
E-mail: tech@clontech.com
www.clontech.com

(PR023487; published May 2010)

3. Remove the culture medium (from Step B.1) and replace with the appropriate culture medium (from Step B.2), to determine the background, the negative control, and the effect of the inducer.
 - **For adherent cells:** after the cells have adhered, remove the media and replace with the appropriate culture medium (from Step B.2).
 - **For cell suspensions:** centrifuge the cells for 5 min at $\leq 1,000$ RPM and remove the media. Resuspend each pellet in the appropriate culture medium (from Step B.2) and transfer each suspension back into the appropriate well/plate/flask.
4. Culture the cells (adherent or suspended) at 37°C for the time required by your experimental protocol.
5. Collect the cells for analysis.
 - **For adherent cells:** collect the cells by trypsinizing; then wash and pellet the cells.
 - **For cell suspensions:** wash and collect the cells by gentle centrifugation for 5 min at $\leq 1,000$ RPM.
6. Resuspend the cell pellet in PBS and analyze the fluorescent signal of the cells using flow cytometry. Use the background control from transfected cells cultured in the absence of Shield1 to determine the overall background. Alternatively, the reporter can be detected using a fluorescence plate reader, especially if the experiment was performed in a 96-well format.
7. Calculate the “fold induction” of promoter activity:

$$\frac{\text{Fluorescence intensity of cells treated with Shield1 \& the inducer of interest}}{\text{Fluorescence intensity of cells treated with Shield1 alone}}$$

Notice to Purchaser

Clontech products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Clontech products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without written approval of Clontech Laboratories, Inc.

cPPT Element:

This product and its use are the subject of U.S. Pat. No. 6,682,907. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer cannot disclose information, sell or otherwise transfer this product, its components or materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for any commercial purposes. If the buyer is not willing to accept the limitations of this limited use statement, Clontech is willing to accept return of the product with a full refund. For information on purchasing a license to the DNA-Flap technology for purposes other than research, contact the Transfer of Technology Office, Institut Pasteur, 28 rue du Docteur Roux, 75 724 Paris Cedex 15 (www.pasteur.fr).

Fruit Fluorescent Proteins and DsRed-Monomer:

DsRed-Monomer and the Fruit Fluorescent Proteins are covered by one or more of the following U.S. Patents: 7,157,566, 7,393,923, 7,005,511, and 7,250,298.

Living Colors® Fluorescent Protein Products:

Not-For-Profit Entities: Orders may be placed in the normal manner by contacting your local representative or Clontech Customer Service at 650.919.7300. At its discretion, Clontech grants Not-For-Profit Entities a non-exclusive, personal, limited license to use this product for non-commercial life science research use only. Such license specifically excludes the right to sell or otherwise transfer this product, its components or derivatives thereof to third parties. No modifications to the protein coding sequence may be made without express written permission from Clontech. Any other use of this product requires a license from Clontech. For license information, please contact a licensing representative by phone at 650.919.7320 or by e-mail at licensing@clontech.com.

For-Profit Entities wishing to use this product are required to obtain a license from Clontech. For license information, please contact a licensing representative by phone at 650.919.7320 or by e-mail at licensing@clontech.com.

ProteoTuner™ Protein Stabilization/Destabilization Products:**Reef Coral Fluorescent Proteins (RCFPs):**

The RCFPs (including DsRed-Express and DsRed-Express2) are covered by one or more of the following U.S. Patents: 7,338,784, 7,338,783, 7,442,522, 7,157,565, 7,217,789, 7,537,915, 6,969,579, 7,150,979, and 7,166,444.

WPRE:

Clontech has a license to sell products containing WPRE, under the terms described below. Any use of WPRE outside of Clontech's product or the product's intended use, requires a license as detailed below. Before using the product containing WPRE, please read the following license agreement. If you do not agree to be bound by its terms, contact Clontech within 10 days for authorization to return the unused product containing WPRE and to receive a full credit.

Patents: The WPRE technology is covered by one or more of the following U.S. Patents and corresponding patent claims outside the U.S.: 6,136,597; 6,284,469; 6,312,912; 6,287,814, issued to The Salk Institute for Biological Studies

Individual License Agreement: Clontech grants you a non-exclusive license to use the enclosed product containing WPRE in its entirety for its intended use. The product is being transferred to you in furtherance of, and reliance on, such license. Any use of WPRE outside of Clontech's product or the product's intended use, requires a license from the Salk Institute for Biological Studies.

Termination of License: This license agreement is effective until terminated. You may terminate it at any time by destroying all products containing WPRE in your control. It will also terminate automatically if you fail to comply with the terms and conditions of the license agreement. You shall, upon termination of the license agreement, destroy all products containing WPRE in your control, and so notify Clontech in writing. This License shall be governed in its interpretation and enforcement by the laws of the State of California.

Contact for WPRE Licensing: The Salk Institute for Biological Studies, 10010 North Torrey Pines Road, La Jolla, CA 92037, Attn.: Office of Technology Management Phone: 858.453.4100 ext. 1275 Fax: 858.546.8093

Clontech, the Clontech logo and all other trademarks are the property of Clontech Laboratories, Inc., unless noted otherwise. Clontech is a Takara Bio Company. ©2010 Clontech Laboratories, Inc.