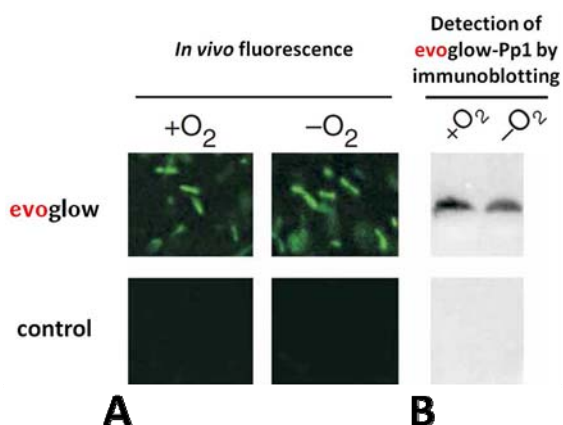


## evoglow<sup>®</sup> Antibodies, whole antiserum

Cat. No.	Amount
FP-21051	70 blots
FP-21052	70 blots

### Product Information

evoglow<sup>®</sup> antibodies are produced in rabbits using PAGE-purified evoglow<sup>®</sup> fluorescent protein as immunogen. They are approved for specific detection of the fluorescence reporters by Western Blot in bacteria<sup>1</sup> as well as yeast<sup>2</sup> (e.g. see Fig.1).



**Fig. 1:** (A) *In vivo*-fluorescence of bacteria cells expressing evoglow<sup>®</sup> under aerobic and anaerobic conditions. (B) Detection of the fluorescence reporter in the cultures shown in (A) by immunoblotting. Cross-reactivity was excluded by cultures expressing harboring the corresponding empty-vector.

### Preparation Instructions

For detection of evoglow<sup>®</sup> protein by Western Blot analysis, the supplied antibody solution is recommended to be diluted 1:10.000, for example with TBS

(Tris-buffered saline). Titers for other applications may vary.

### Storage / Stability

For continuous use, store at 2 - 8 °C for up to one month. For extended storage, the solution may be frozen in working aliquots. Repeated freeze/thaw cycles are not recommended. Storage in "frost-free" freezers is not recommended. If slight turbidity occurs upon prolonged storage, clarify the solution by centrifugation before use.

### Precautions and Disclaimer

For laboratory use only. Not for drug, household or other uses. The antibodies are not intended for the use in human medical or veterinarian medical diagnostic purposes, or in the animal or human body; any liability of the company Jena Bioscience GmbH is excluded.

### References

- 1: Drepper, T. *et al.* (2007) *Nat. Biotechnol.* 25: 443-445.
- 2: Tielker, D. *et al.* (2009) *Eukaryot Cell.* 8(6):913-915.

## evoglow<sup>®</sup> Antibodies, whole antiserum

### Example protocol for immunodetection of immobilized evoglow<sup>®</sup> on a blotting membrane

Note: For a detailed description of the western blot process itself, please use the recommendations provided by the manufacturer of your blotting device

### Required material / working solutions:

- blotting membrane with immobilized evoglow<sup>®</sup> fluorescent protein
- evoglow<sup>®</sup> antibodies
- secondary antibodies of choice (appropriate type: anti-rabbit specificity). For further information please refer to your antibody supplier.
- TBST\* (working solution):  
50 mM Tris-HCl, pH 6.8  
150 mM NaCl  
1 mM MgCl<sub>2</sub>  
0.2 % (v/v) Tween 20
- blocking solution: TBST supplemented with 5 % (w/v) skim milk

### Detailed procedure for immunodetection of evoglow<sup>®</sup> fluorescent proteins

General handling hints: Perform all incubation and washing steps with slight agitation at 30 °C (e.g. on a rocking platform). Handling of membranes should be carried out with gloves and clean trays to avoid contamination.

1. Incubate the membrane with the immobilized evoglow protein for 1 h (alternatively incubate at 4 °C overnight) with blocking solution. This step will saturate the membrane to prevent direct binding of the antibody and unspecific background signals.
2. Wash the membrane 2 times for 5 min with TBST to remove the skim milk.
3. During incubation (step 2), prepare a 1:10.000 dilution of the evoglow<sup>®</sup> antibody solution in TBST.
4. Incubate the membrane with the evoglow<sup>®</sup> antibody working solution for 1 h.
5. Wash the membrane twice with TBST for 30 min.
6. Incubate with the secondary antibodies of choice (e.g. goat anti-rabbit HRP-conjugate) according to recommendation of the respective manufacturer.
7. Wash the membrane twice with TBST for 30 min.
8. Wash the membrane twice with TBST for 5 min.
9. Perform the appropriate detection reaction depending on the used secondary antibody type. For detailed instructions refer to the manual of the secondary antibody manual.

\* **TBST:** Tris-buffered saline with Tween

**Note:** This protocol is just an example and can be regarded as supporting guideline. Protocols for other applications may vary.