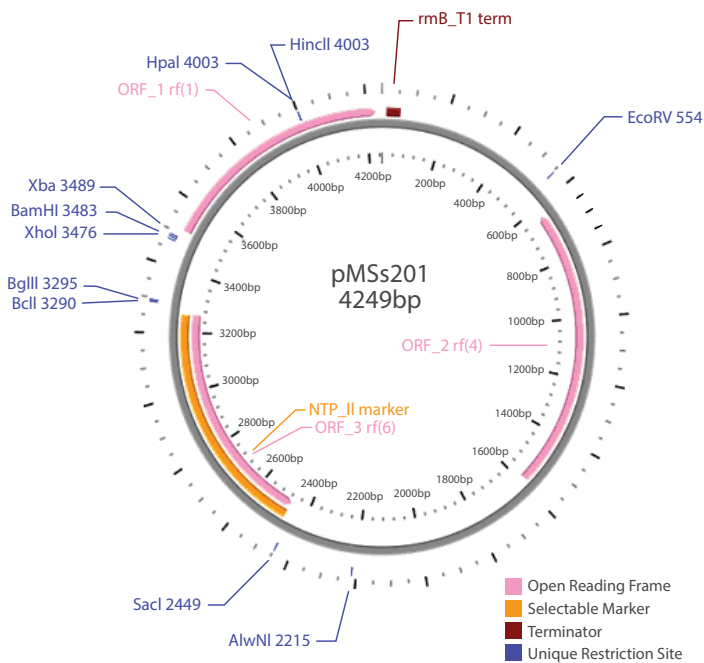


# *E. coli* promoter collection



**Figure 1. pMS201 vector map.**

**Table 1. Antibiotic resistances conveyed by pMS201.**

Antibiotic	Concentration	Utility
Kanamycin	25 µg/mL	Bacterial selection marker

## Product description

Researchers at the Weizmann Institute of Science have produced a collection of *E. coli* strains that enables monitoring of gene expression at high temporal resolution in living cells. Each of the reporter strains has a bright, fast-folding green fluorescent protein (GFP) fused to a full-length copy of an *E. coli* promoter in a low-copy plasmid. This collection includes more than 1900 promoters (out of 2500 in the entire genome) for *E. coli* K12 strain MG1655 and enables measurement of gene expression at a resolution of minutes with high accuracy and reproducibility. Performing experiments in a multi-well fluorimeter using FACS or time-lapse fluorescence microscopy has the necessary sensitivity to measure gene expression in individual cells.

## Protocol I—replication

### Materials for replication

2x LB broth (low salt)  
Peptone, granulated, 2 kg—Difco  
Yeast Extract, 500 g, granulated  
Glycerol  
Kanamycin

### 2x LB broth (low salt) medium preparation

1 x LB broth	10 g/L
Peptone	10 g/L
Yeast Extract	5 g/L
Appropriate antibiotic(s)	at recommended concentration(s)

For archive replication, grow all clones at 37 °C in 2x LB broth (low salt) medium plus 25 µg/mL kanamycin. Prepare medium with 8% glycerol\* and the appropriate antibiotics.

\*Glycerol should be omitted from the medium if you are culturing for plasmid preparation. If making copies of the constructs for long-term storage at -80 °C, 8% glycerol is required.

Freeze at -80 °C for long term storage. Avoid long periods of storage at room temperature or higher in order to control background recombination products.

## Protocol II—plasmid preparation

### Materials for replication

For plasmid preparation, grow all clones at 37 °C in 2x LB broth (low salt) medium plus 25 µg/mL kanamycin.

Most plasmid mini-prep kits recommend a culture volume of 1–10 mL for good yield. For these constructs, 5 mL of culture can be used for one plasmid mini-prep generally producing 5–10 µg of plasmid DNA.

1. Upon receiving your glycerol stock(s), store at –80 °C until ready to begin.
2. To prepare plasmid DNA, first thaw your glycerol stock culture and pulse vortex to resuspend any *E. coli* that may have settled to the bottom of the tube.
3. Take a 10 µL inoculum from the glycerol stock into 3–5 mL of 2x LB broth (low salt) with 25 µg/mL kanamycin. Return the glycerol stock(s) to –80 °C.



If a larger culture volume is desired, incubate the 3–5 mL culture for 8 hours at 37 °C with shaking and use as a starter inoculum. Dilute the starter culture 1:500–1:1000 into the volume.

4. Incubate at 37 °C for 18–19 hours with vigorous shaking.
5. Pellet the 3–5 mL culture and begin preparation of plasmid DNA.
6. Run 3–5 µL of the plasmid DNA on a 1% agarose gel. pMS201 without ORF is 4260 bps.

### What clones are part of my collection?

A USB drive containing the data for this collection will be shipped with each collection. This data file can be downloaded from the *E. coli* Promoter product page on our [website](#).

### What antibiotic should I use?

You should grow all *E. coli* Promoter clones in 2x LB broth (low salt) with 25 µg/mL kanamycin for archive replication.

## References

1. Kalir S *et al.*, (2001). Ordering genes in a flagella pathway by analysis of expression kinetics from living bacteria. *Science* **292**, 2080–2083.
2. Ronen M *et al.*, (2002). Assigning numbers to the arrows: parameterizing a gene regulation network by using accurate expression kinetics. *Proc. Natl. Acad. Sci. USA* **99**, 10555–10560.
3. Setty *et al.*, (2003). *Proc. Natl. Acad. Sci. USA* **100**, 7702–7707.
4. Zaslaver A *et al.*, (2004). Just-in-time transcription program in metabolic pathways. *Nature Genetics* **36**, 486–491.
5. Zaslaver A, Anat Bren, Michal Ronen, Shalev Itzkovitz, Ilya Kikoin, Seagull Shavit, Wolfram Liebermeister, Michael G Surette & Uri Alon (2006). A comprehensive library of fluorescent transcriptional reporters for *Escherichia coli*. *Nature Methods* **3**, 623–628.

### If you have any questions, contact

**t** +44 (0) 1223 976 000 (UK) **or** +1 800 235 9880 (USA); +1 303 604 9499 (USA)

**f** +44 (0)1223 655 581

**w** [horizondiscovery.com/contact-us](http://horizondiscovery.com/contact-us) **or** [dharmacon.horizondiscovery.com/service-and-support](http://dharmacon.horizondiscovery.com/service-and-support)

**Horizon Discovery**, 8100 Cambridge Research Park, Waterbeach, Cambridge, CB25 9TL, United Kingdom

**horizon**<sup>TM</sup>  
INSPIRED CELL SOLUTIONS

V2-0418