

The pETBlue™ vectors are designed to identify recombinants by traditional blue/white screening while also allowing T7 *lac* promoter based expression of target genes. Screening is independent of expression because the T7 *lac* expression promoter is in an opposed orientation relative to the *E. coli* promoter that mediates blue/white screening. pETBlue-1 is a vector that facilitates the expression of native unfused proteins and allows convenient subcloning of target genes already fused to existing detection and purification tags. The *EcoR* V cloning site is appropriately spaced downstream of an *E. coli* ribosome binding site. Inserts must encode an ATG start codon at their 5' end if expression is desired. The cloning/expression region of the coding strand transcribed by T7 RNA polymerase is shown below. The sequence is numbered from the first base of the T7 promoter sequence. Unique sites are shown on the circle map. The f1 origin in pETBlue-1 is oriented so that infection with helper phage will produce virions containing single stranded DNA that corresponds to the T7 RNA polymerase coding strand. Therefore, single stranded sequencing should be performed using the pETBlueDOWN primer (Cat. No. 70603-3).

pETBlue-1 sequence landmarks

<i>lac</i> operator	3429–3448
T7 promoter	1–17
<i>lac</i> operator	22–42
T7 transcription start	18
multiple cloning region (<i>EcoR</i> V– <i>Srf</i> I)	276–297
<i>lacZ</i> start codon	314
<i>lacZ</i> α -peptide ORF	57–314
<i>E. coli</i> promoter	364–392
f1 origin	919–1374
<i>bla</i> coding sequence	1492–2349
pUC origin	3029



