

Product Information

N-CMV-30 Sequencing Primer

Catalog Number **P5350**

Store Temperature $-20\text{ }^{\circ}\text{C}$

Product Description

Nucleotide sequence:

5'-AAT-GTC-GTA-ATA-ACC-CCG-CCC-CGT-TGA-
-CGC-3'

N-CMV-30 sequencing primer is a chemically synthesized, 30 base, single strand oligodeoxy-ribonucleotide designed for double-stranded or single-stranded DNA sequencing of FLAG® fusion proteins at the N-terminus of the pFLAG-CMV expression vectors.

The primer is complementary to map position nucleotides 818-847 of positive, coding DNA strand of the pFLAG-CMV vectors.

1

The N-CMV-30 sequencing primer is supplied at an initial concentration of 5 μM in 1 \times TE buffer (10 mM Tris, pH 8.0, 1.0 mM EDTA).

Reagents Required but Not Provided

- 1 \times TE buffer
- 5 M NaOH
- 3 M potassium/5 M acetate: To 60 ml 5 M potassium acetate add 11.5 ml glacial acetic acid and 28.5 ml H₂O. The resulting solution is 3 M with respect to potassium and 5 M with respect to acetate.
- Isopropanol

Procedure

A. Preparation of Denatured DNA Template

(For sequencing single-stranded DNA, proceed to Section B.)

The following protocol is designed to allow sufficient template for two sets of 4 sequencing reactions (2G, 2A, 2T and 2 C reactions).

1. Denature 7 micrograms of highly purified pFLAG-CMV supercoiled DNA in 70 μl of 1 \times TE by adding 3 μl (1/25 volume) of 5 M NaOH and incubating at 37 $^{\circ}\text{C}$ for 5 minutes.

2. Precipitate the denatured pFLAG-CMV supercoiled DNA for 30 minutes at room temperature by adding 150 μl (2 volumes) of 3 M potassium/5 M acetate: isopropanol (1:3 ratio mixture).
3. Collect the precipitated and denatured pFLAG-CMV DNA by centrifugation at 10,000 $\times g$ for 5 minutes.
4. Wash the pFLAG-CMV DNA pellet with 1 ml ethanol and dry.
5. Resuspend the pFLAG CMV DNA in 20 μl of 1 \times TE. The final concentration should be 0.35 $\mu\text{g}/\mu\text{l}$.

The irreversibly denatured pFLAG-CMV DNA template can be stored at $-20\text{ }^{\circ}\text{C}$ at this point if desired.

B. Priming pFLAG-CMV DNA Template with N-CMV-30 Sequencing Primer:

The following protocol uses 10 μl or half of the preceding preparation of irreversibly denatured pFLAG-CMV DNA template. This is sufficient for one set of 4 sequencing reactions (G, A, T and C).

1. Dilute a 3 μl aliquot of the N-CMV-30 sequencing primer with 6 μl of 1 \times TE to make a final concentration of 1.67 pmol/ μl .
2. Add 2-3 μl (3.5-5 pmoles) of N-CMV-30 sequencing primer to 10 μl (1 pmole) of denatured pFLAG-CMV DNA template.
3. To 12-13 μl of N-CMV-30 primer/pFLAG-CMV DNA template add an appropriate volume of sequencing buffer to make the buffer concentration 1 \times . Heat at 70 $^{\circ}\text{C}$ in a wet temperature block for 2 minutes.
4. Slowly cool the mixture to 45 $^{\circ}\text{C}$ by placing the temperature block at room temperature for about 20 minutes.

5. Distribute the N-CMV-30 primer/pFLAG-CMV DNA template to four tubes to be used in the G, A, T, and C DNA sequencing reactions.

The N-CMV-30 primer/pFLAG-CMV DNA template is now ready for supercoil sequencing of the DNA sequence corresponding to the N-terminal FLAG fusion junction.

The DNA sequence corresponding to the fusion junction will be >143 bases away from the 3' end of the N-CMV-30 sequencing primer using pFLAG-CMV-1 template and >98 bases away using pFLAG-CMV-2 template reading from the bottom of an autoradiogram.

FLAG is a registered trademark of Sigma-Aldrich® Biotechnology LP and Sigma-Aldrich Co.

MAM 09/10-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.