

## Product Information

**M-MLV Reverse Transcriptase,**  
recombinant,  
expressed in *E. coli*

Product Code **M 1302**

Storage Temperature -20 °C

## TECHNICAL BULLETIN

### Product Description

M-MLV (Moloney Murine Leukemia Virus) Reverse Transcriptase is a DNA polymerase that uses DNA, single-stranded RNA, or an RNA:DNA hybrid (using a primer) to synthesize a complementary DNA strand. This enzyme is isolated from *E. coli* expressing a portion of the *pol* gene of M-MLV on a plasmid.<sup>1</sup> M-MLV is used for the preparation of cDNA libraries or for first strand cDNA synthesis for use in RT-PCR<sup>†</sup> reactions.

Unit Definition: One unit incorporates 1 nmole of TTP into acid precipitable material in 10 minutes at 37 °C.<sup>2</sup>

### Components

- M-MLV Reverse Transcriptase, Product Code M1427  
200 units/μl in 20mM Tris-HCl (pH 7.5), 200mM NaCl, 0.1mM EDTA, 1mM DTT, 0.01% Nonidet® P-40 and 50% glycerol.  
Provided as 40,000 or 200,000 units
- 10X M-MLV Reverse Transcriptase Buffer, Product Code B8559  
500 mM Tris-HCl, pH 8.3, with 500 mM KCl, 30 mM MgCl<sub>2</sub>, and 50 mM DTT  
Provided as 0.5 mL vials, 1 vial/40,000 units and 4 vials/200,000 units

### Reagents Required but Not Provided

(Product Codes have been given where appropriate.)

- Deoxynucleotide Mix, Product Code D7295, 10 mM dATP, 10 mM dCTP, 10 mM dGTP, 10 mM dTTP
- Water, Product Code W1754
- Choice of specific primer (user defined), anchored oligo d(T)<sub>23</sub> primers (Product Code O4387) or random nonamers (Product Code R7647).
- RNA template
- Ribonuclease Inhibitor, Product Code R2520

### Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Procedure

The following procedure uses 1-5 μg of total RNA or approximately 50 ng of mRNA.

1. In a clean, nuclease-free 0.5 ml or 0.2 ml thin-walled microcentrifuge tube, add the following:
  - 1 μl 10 mM dNTP mix
  - 1 μl Specific primer, random nonamers or oligo(dT)<sub>23</sub> (final concentration 1-5 μM)
  - 1 μl RNA template
  - q.s. Nuclease-free water
  - 10 μl Total volume
2. Mix gently and briefly centrifuge to collect all components to the bottom of the tube.

3. Incubate at 70 °C for 10 minutes.
4. Remove the tubes and place on ice.
5. Add the remaining components to the microcentrifuge tube:
  - 2 µl 10x M-MLV Reverse Transcriptase Buffer
  - 1 µl M-MLV Reverse Transcriptase
  - 0.5 µl RNase Inhibitor (40 units/µl).
  - 6.5 µl Nuclease-free water
  - 20 µl Final Volume

6. Incubate at 37 °C for 50 minutes.  
Optional step: Incubate reaction at room temperature for 10 minutes prior to the 37 °C incubation for 50 minutes. This will ensure elongation of random primers or anchored oligo(dT)<sub>23</sub> primers before the higher reverse transcriptase temperature.
7. The cDNA strand has now been produced. Heat the reaction tube between 80 °C and 94 °C for 10 minutes to denature the M-MLV reverse transcriptase.

## References

1. Gerard, G.F., *et al.*, Influence on stability in *Escherichia coli* of the carboxy-terminal structure of cloned Moloney murine leukemia virus reverse transcriptase. *DNA*, **5**:271-9 (1986).
2. Houts, G.E., *et al.*. Reverse transcriptase from avian myeloblastosis virus. *J. Virol.* **29**:517-22 (1979).
3. Howland, P., *et al.* Positive- and negative-acting promoter sequences regulate cell type-specific expression of the rat synapsin I gene. *Mol. Brain Res.* **11**:345-53.
4. Gerard, G.F. and D'Alessio, J.M. *Methods in Molecular Biology* Vol. 16: Enzymes of Molecular Biology, (Buwell, M. Ed.) Humana Press, Totowa, N.J., p. 73 (1993)

†The PCR process is covered by patents owned by Hoffman-LaRoche, Inc.

TRITON is a registered trademark of Union Carbide Corp.

VS,JLL,MAM,SM 08/20-1