

## **m<sup>7</sup>G(5')ppp(5')G RNA Capping Analog**

**Cat. No. 15619-018**

**Size: 25 OD<sub>260</sub> units**

**Store at -20°C.**

### Description:

The 5' terminal m<sup>7</sup>G cap present on most eukaryotic mRNAs promotes translation *in vitro* at the initiation level (1,2,3). For most RNAs, elimination of the cap structure causes a loss of stability, especially against exonuclease degradation (4), and a decrease in the formation of the initiation complex of mRNAs for protein synthesis (4,5). Certain prokaryotic mRNAs containing a 5'-terminal cap structure are translated as efficiently as or more efficiently than eukaryotic mRNAs in a eukaryotic cell-free protein synthesizing system (5). Also a cap requirement has been observed for splicing eukaryotic substrate RNAs (6).

A method using *E.coli* RNA polymerase primed with m<sup>7</sup>G(5')ppp(5')G or m<sup>7</sup>G(5')ppp(5')A for an efficient *in vitro* synthesis of capped RNAs has been developed by Contreas (7). Larger amounts of capped RNAs are produced by transcription systems using SP6 RNA polymerase primed with m<sup>7</sup>G(5')ppp(5')G (6).

### Unit Definition:

Addition of 137 µl of water gives approximately a 10 mM solution.

MW = 846 (2 Na, water not determined)  
λ<sub>max</sub> = 254 nm

E = 18,200  
21.5 OD/mg

25 OD units = 1.16 mg = 1.37 × 10<sup>-6</sup> moles  
25 OD units dissolved in 137 µl of water = 10 mM solution

Doc. Rev.: 05/18/01

Quality Control:

This product has passed the following quality control assays: Determination of optical density and performance evaluation in an *in vitro* translation assay using CAT mRNA.

References:

1. Shatlin, A.J. (1978) *Cell* 9, 645.
2. Filipowicz, W. (1978) *FEBS Lett* 96, 1.
3. Banerjee, A.K. (1980) *Microbiol. Rev.* 44, 175.
4. Miura, K. (1981) *Adv. Biophys.* 14, 205.
5. Shatkin, A.J., *et al* (1977) *Nucleic Acids Res.* 4, 3065.
6. Konarska, M.M., *et al* (1984) *Cell* 38, 731.
7. Contreas, R., *et al* (1982) *Nucleic Acids Res.* 10, 6353.
8. Paterson, B.M. and Rosenberg, M. (1979) *Nature* 279, 696.