



Random Primers

Cat. No. 48190-011

Size: 9 A₂₆₀ units

Conc.: 3 µg/µl

Store at -20°C

Description:

The solution contains Random Primer oligonucleotides (mostly hexamers), at a concentration of 0.09 OD₂₆₀ units/µl, in 3 mM Tris-HCl (pH 7.0), 0.2 mM EDTA.

Quality Control:

Incorporation of radioactively labeled nucleotides is verified using control DNA in a random primers labeling reaction.

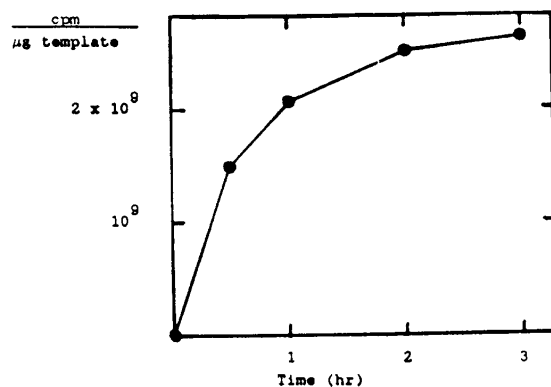
Notes:

Random Primers can be used to prepare labeled DNA probes from DNA templates (1,2,3,) or cDNA from RNA (4,5).

Doc. Rev.: 062001

This product is distributed for laboratory research only. **CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.**

For technical questions about this product, call the Invitrogen Tech-Line™ U.S.A. 800 955 6288



References:

1. Feinberg, A. P. and Vogelstein, B. (1983) *Anal. Biochem.* 132, 6.
2. Feinberg, A. P. and Vogelstein, B. (1984) *Anal. Biochem.* 137, 266.
3. Thomas-Cavallin, M. and Ait-Ahmed, O. (1988) *J. Histochem. Cytochem.* 36, 1335.
4. Haymerle, H., Herz, J., Bressan, G. M., Frank R. and Stanley, K. K. (1986) *Nucl. Acids Res.* 14, 8615.
5. Persons, D.A. and Finn, O. J. (1986) *BioTechniques* 4, 398.