

Random Octamers

Random Primers

Primers and Oligonucleotides

	Cat.-No.	Size	Conc.
	PM-302S	200 µl	100 µM
	PM-302L	1 ml	100 µM

For *in vitro* use only

Quality guaranteed for 12 months

Store at -20 °C, avoid frequent thawing and freezing

Description

Random Octamers are a mixture of oligonucleotides representing all possible sequences for a hexamer. Random Octamers are used in DNA labeling by PCR (DOP-PCR) or cDNA synthesis by RT-PCR.

Sequence

5'-NNN NNN NN-3'

Molecular weight

2410

Package size and conversion

Random Hexamers			
PM-302S	200 µl / 100 µM	20 nmol	50 µg
PM-302L	1 ml / 100 µM	100 nmol	250 µg

Random Octamers (white cap)

100 µM in aqueous solution