

**Random Octamers**

Cat. No.	Amount
PM-302S	200 µl (20 nmol, 50 µg)
PM-302L	5 x 200 µl (100 nmol, 250 µg)

For *in vitro* use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 2410 g/mol

Purity: by HPLC

Form: clear aqueous solution

Concentration: 100 µM

Description:

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

Sequence:

5'- NNN NNN NN -3'