



## Random Octamers

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

**For general laboratory use.**

**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'