



Random Hexamers

100 μ M

| Cat. No. | Amount |
|----------|-------------|
| PM-301S | 200 μ l |
| PM-301L | 1 ml |

For *in vitro* use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 1793 g/mol

Purity: by HPLC

Form: clear aqueous solution

Concentration: 100 μ M

Description:

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

Sequence:

5'- NNN NNN -3'

Package size and conversion:

| | | | |
|---------|---------------------------|----------|-------------|
| PM-301S | 200 μ l / 100 μ M | 20 nmol | 37 μ g |
| PM-301L | 1 ml / 100 μ M | 100 nmol | 185 μ g |