

Primers and Oligonucleotides

DNA and RNA Oligos

Content

| | |
|---|---|
| Unmodified DNA Oligonucleotides | 2 |
| Modified DNA Oligonucleotides - 5' Fluorescent Labels | 2 |
| Modified DNA Oligonucleotides - 5' Non-Fluorescent Labels | 3 |
| Modified DNA Oligonucleotides - 3' Fluorescent Labels | 3 |
| Modified DNA Oligonucleotides - 3' Non-Fluorescent Labels | 4 |
| Modified DNA Oligonucleotides - Internal Fluorescent Labels | 4 |
| Modified DNA Oligonucleotides - Internal Non-Fluorescent Labels | 5 |
| Antisense DNA Oligos - Phosphorothioates (PTOs) | 5 |
| Additional Documentation | 5 |
| Unmodified RNA Oligonucleotides | 6 |
| Modified RNA Oligonucleotides - 5' Labels | 6 |
| Modified RNA Oligonucleotides - 3' Labels | 6 |
| Shipping Charge | 6 |

Primers and Oligonucleotides

DNA and RNA Oligos

Unmodified DNA Oligonucleotides

| Scale [µmol] | Standard Purification * | | OPC Purification * | | HPLC Purification | |
|---------------------------------|-------------------------|----------------------------|--------------------|----------------------------|-------------------|----------------------------|
| | price per base | Yield [OD ₂₆₀] | price per base | Yield [OD ₂₆₀] | price per base | Yield [OD ₂₆₀] |
| 0.02 ** (<34 bases only) | 0.29 € | 3 | 0.75 € | 1 | 0.85 € | 1 |
| 0.04 | 0.37 € | 5 | 0.80 € | 2.5 | 0.98 € | 2.5 |
| 0.2 | 0.98 € | 16 | 1.20 € | 8 | 1.49 € | 8 |
| 1.0 | 2.40 € | 80 | - | - | 2.98 € | 25 |
| 10 | - | - | - | - | 20.00 € | - |

* Standard and OPC purification can only be ordered for oligos <45 bases.

** The 0.02 µmol scale can only be ordered for oligos <34 bases.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

No extra charge for 5' and internal wobbles (degenerated bases). 3' wobbles require a setup fee of 20 €.

For technical reasons we have to double the price per base for oligos >80 bases.

2

Modified DNA Oligonucleotides - 5' Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|-----------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|---------------------------------|----------------------------------|
| | 0.02 µmol 3 OD ₂₆₀ | 0.04 µmol 5 OD ₂₆₀ | 0.2 µmol 12 OD ₂₆₀ | 1.0 µmol 50 OD ₂₆₀ | 0.02 µmol 1 OD ₂₆₀ | 0.04 µmol 2 OD ₂₆₀ | 0.2 µmol 5 OD ₂₆₀ | 1.0 µmol 12 OD ₂₆₀ |
| 6-FAM, TET | 27 € | 29 € | 45 € | 85 € | 30 € | 33 € | 55 € | 100 € |
| HEX, TAMRA | - | - | - | - | 35 € | 40 € | 70 € | 150 € |
| JOE, ROX | - | - | - | - | 58 € | 65 € | 100 € | 175 € |
| Fluorescein | 27 € | 29 € | 45 € | 85 € | 30 € | 33 € | 55 € | 100 € |
| Cy3, Cy5 | - | - | - | - | 50 € | 56 € | 95 € | 202 € |
| IRD 700/800 | - | - | - | - | 38 € | 42 € | 50 € | 180 € |
| CAL Fluor Gold 540 | - | - | - | - | 75 € | 85 € | 99 € | 180 € |
| CAL Fluor Orange 560 | - | - | - | - | 50 € | 55 € | 75 € | 180 € |
| CAL Fluor Red 610 | - | - | - | - | 80 € | 90 € | 120 € | 150 € |
| Rhodamine ITC | - | - | - | - | 37 € | 42 € | 50 € | 115 € |
| Rhodamine Green | - | - | - | - | 95 € | 110 € | 140 € | 240 € |
| Texas Red | 150 € | 165 € | 190 € | 300 € | 155 € | 170 € | 200 € | 320 € |

Standard purification can only be ordered for oligos <45 bases.

The 0.02 µmol scale can only be ordered for oligos <34 bases.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Primers and Oligonucleotides

DNA and RNA Oligos

Modified DNA Oligonucleotides - 5' Non-Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|-----------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|---------------------------------|----------------------------------|
| | 0.02 µmol 3 OD ₂₆₀ | 0.04 µmol 5 OD ₂₆₀ | 0.2 µmol 12 OD ₂₆₀ | 1.0 µmol 50 OD ₂₆₀ | 0.02 µmol 1 OD ₂₆₀ | 0.04 µmol 2 OD ₂₆₀ | 0.2 µmol 5 OD ₂₆₀ | 1.0 µmol 12 OD ₂₆₀ |
| Biotin | 30 € | 37 € | 47 € | 120 € | 33 € | 40 € | 55 € | 135 € |
| C8-Alkyne-dC | 25 € | 28 € | 33 € | 60 € | 26 € | 29 € | 37 € | 65 € |
| C8-Alkyne-dU | 15 € | 17 € | 27 € | 57 € | 16 € | 18 € | 30 € | 62 € |
| C6-Amino | 10 € | 12.50 € | 15 € | 20 € | 15 € | 20 € | 25 € | 30 € |
| C12-Amino | 23 € | 32 € | 40 € | 80 € | 26 € | 35 € | 50 € | 95 € |
| 2-Deoxyuridine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Digoxigenin | 65 € | 70 € | 85 € | 140 € | 69 € | 75 € | 95 € | 155 € |
| Hexynyl | 8 € | 10 € | 13 € | 27 € | 9 € | 11 € | 17 € | 32 € |
| Inosine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Methylcytosin | 15 € | 17 € | 20 € | 40 € | 17.50 € | 20 € | 25 € | 50 € |
| Phosphate | 15 € | 22 € | 30 € | 35 € | 18 € | 25 € | 35 € | 40 € |
| Thiol * | 49 € | 55 € | 65 € | 140 € | 53 € | 60 € | 75 € | 155 € |

Standard purification can only be ordered for oligos <45 bases.

The 0.02 µmol scale can only be ordered for oligos <34 bases.

* A set up fee of 110 € per order will be charged.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Modified DNA Oligonucleotides - 3' Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|--------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|---------------------------------|----------------------------------|
| | 0.02 µmol 3 OD ₂₆₀ | 0.04 µmol 5 OD ₂₆₀ | 0.2 µmol 12 OD ₂₆₀ | 1.0 µmol 50 OD ₂₆₀ | 0.02 µmol 1 OD ₂₆₀ | 0.04 µmol 2 OD ₂₆₀ | 0.2 µmol 5 OD ₂₆₀ | 1.0 µmol 12 OD ₂₆₀ |
| 6-FAM | - | - | - | - | 30 € | 33 € | 55 € | 200 € |
| HEX | - | - | - | - | 50 € | 60 € | 115 € | 400 € |
| TAMRA | - | - | - | - | 50 € | 60 € | 75 € | 220 € |
| Fluorescein | - | - | - | - | 30 € | 33 € | 45 € | 150 € |
| Cy3, Cy5 | - | - | - | - | 50 € | 61 € | 95 € | 213 € |

Standard purification can only be ordered for oligos <45 bases. The 0.02 µmol scale can only be ordered for oligos <34 bases.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Prices for 3' - fluorescent modifications are valid up to 40 bases, for longer oligonucleotides or other modifications please inquire.

Primers and Oligonucleotides

DNA and RNA Oligos

Modified DNA Oligonucleotides - 3' Non-Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|-----------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|---------------------------------|----------------------------------|
| | 0.02 µmol 3 OD ₂₆₀ | 0.04 µmol 5 OD ₂₆₀ | 0.2 µmol 12 OD ₂₆₀ | 1.0 µmol 50 OD ₂₆₀ | 0.02 µmol 1 OD ₂₆₀ | 0.04 µmol 2 OD ₂₆₀ | 0.2 µmol 5 OD ₂₆₀ | 1.0 µmol 12 OD ₂₆₀ |
| Alkyne | 9 € | 11 € | 17 € | 33 € | 10 € | 12 € | 20 € | 38 € |
| C8-Alkyne-dC | 25 € | 28 € | 33 € | 60 € | 26 € | 29 € | 37 € | 65 € |
| C8-Alkyne-dU | 15 € | 17 € | 27 € | 57 € | 16 € | 18 € | 30 € | 62 € |
| C7-Amino | 12 € | 15 € | 30 € | 90 € | 15 € | 20 € | 40 € | 105 € |
| BHQ-1 | - | - | - | - | 55 € | 75 € | 115 € | 355 € |
| BHQ-2 | - | - | - | - | 65 € | 85 € | 135 € | 395 € |
| BHQ-3 | - | - | - | - | 70 € | 95 € | 140 € | 400 € |
| Biotin | 30 € | 37 € | 47 € | 120 € | 33 € | 40 € | 55 € | 135 € |
| Dabcyl | - | - | - | - | 50 € | 70 € | 95 € | 195 € |
| ddC | 28 € | 48 € | 68 € | 200 € | 30 € | 50 € | 70 € | 205 € |
| 2-Deoxyuridine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Inosine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Methylcytosin | 15 € | 17 € | 20 € | 40 € | 17.50 € | 20 € | 25 € | 50 € |
| Phosphate | 15 € | 22 € | 30 € | 35 € | 18 € | 25 € | 35 € | 40 € |
| Thiol | 30 € | 40 € | 50 € | 95 € | 35 € | 45 € | 60 € | 110 € |

Standard purification can only be ordered for oligos <45 bases.

The 0.02 µmol scale can only be ordered for oligos <34 bases.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Prices for 3' - non-fluorescent modifications are valid up to 40 bases, for longer oligonucleotides or other modifications please inquire.

Modified DNA Oligonucleotides - Internal Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|----------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|----------------------------------|---------------------------------|----------------------------------|
| | 0.02 µmol 3 OD ₂₆₀ | 0.04 µmol 5 OD ₂₆₀ | 0.2 µmol 12 OD ₂₆₀ | 1.0 µmol 50 OD ₂₆₀ | 0.02 µmol 1 OD ₂₆₀ | 0.04 µmol 2 OD ₂₆₀ | 0.2 µmol 5 OD ₂₆₀ | 1.0 µmol 12 OD ₂₆₀ |
| FITC-dT | - | - | - | - | 150 € | 200 € | 240 € | 400 € |
| Fluorecein-dT | - | - | - | - | 150 € | 200 € | 240 € | 400 € |
| TAMRA-dT | - | - | - | - | 150 € | 200 € | 240 € | 400 € |
| Cy5-dT | - | - | - | - | 168 € | 225 € | 270 € | 449 € |

Standard purification can only be ordered for oligos <45 bases. The 0.02 µmol scale can only be ordered for oligos <34 bases.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Prices for internal fluorescent modifications are valid up to 40 bases. All prices for internal fluorescent modifications include HPLC purification.

Primers and Oligonucleotides

DNA and RNA Oligos

Modified DNA Oligonucleotides - Internal Non-Fluorescent Labels

Prices per modification are added to the price of the oligo synthesized in standard purification.

| Modification | Standard Purification | | | | HPLC Purification | | | |
|-----------------------|---|---|---|---|---|---|--|---|
| | 0.02 μmol 3 OD ₂₆₀ | 0.04 μmol 5 OD ₂₆₀ | 0.2 μmol 12 OD ₂₆₀ | 1.0 μmol 50 OD ₂₆₀ | 0.02 μmol 1 OD ₂₆₀ | 0.04 μmol 2 OD ₂₆₀ | 0.2 μmol 5 OD ₂₆₀ | 1.0 μmol 12 OD ₂₆₀ |
| C8-Alkyne-dC | 25 € | 28 € | 33 € | 60 € | 26 € | 29 € | 37 € | 65 € |
| C8-Alkyne-dU | 15 € | 17 € | 27 € | 57 € | 16 € | 18 € | 30 € | 62 € |
| C2-Amino-dT | - | - | - | - | 70 € | 80 € | 120 € | 290 € |
| Biotin-dT * | - | - | - | - | 95 € | 105 € | 145 € | 360 € |
| Bromo-dC | - | - | - | - | 78 € | 91 € | 104 € | 208 € |
| Bromo-dG | - | - | - | - | 116 € | 116 € | 149 € | 238 € |
| Bromo-dU | - | - | - | - | 78 € | 91 € | 104 € | 208 € |
| 7-Deaza-dA | - | - | - | - | 76 € | 76 € | 105 € | 228 € |
| 2-Deoxyuridine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Inosine | 7.50 € | 10 € | 12 € | 15 € | 12.50 € | 15 € | 17.50 € | 20 € |
| Iodo-dC | - | - | - | - | 116 € | 116 € | 210 € | 332 € |
| Iodo-dU | - | - | - | - | 83 € | 83 € | 105 € | 166 € |
| Methylcytosin | 15 € | 17 € | 20 € | 40 € | 17.50 € | 20 € | 25 € | 50 € |

Standard purification can only be ordered for oligos <45 bases.

The 0.02 μmol scale can only be ordered for oligos <34 bases.

* A set up fee of 250 € per order will be charged.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Prices for internal non-fluorescent modifications are valid up to 40 bases.

All prices for internal non-fluorescent modifications include HPLC purification.

Antisense DNA Oligos - Phosphorothioates (PTOs)

| Scale [μmol] | € per Base | Yield [OD ₂₆₀] |
|---------------------------|------------|----------------------------|
| 0.2 | 2.90 | 5 |
| 1.0 | 5.80 | 15 |
| 10 | 21.40 | 100 |

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Prices include HPLC- and Cartridge-Purification if the oligo is a "Full-PTO".

For mixed Phosphorothioate/Phosphordiester-Oligos (i. e. "end capping") Cartridge-Purification will be charged at 15 €.

Additional Documentation

Upon special order, MALDI-ToF documentation is available for 20 € per oligo.

Primers and Oligonucleotides

DNA and RNA Oligos

Unmodified RNA Oligonucleotides

| Scale [μmol] | Yield [OD_{260}] | Oligos <12nt | Oligos $\geq 12\text{nt}$ |
|---------------------------|-----------------------------|-----------------|---------------------------|
| 0.2 | 8 | 99 € per oligo | 9.20 € per base |
| 1.0 | 18 | 111 € per oligo | 10.40 € per base |

All RNA Oligonucleotides are HPLC purified.

Guaranteed yields apply for a 20mer +/- 20 %. For oligos >33 bases we cannot give a yield guarantee.

Modified RNA Oligonucleotides - 5' Labels

Prices per modification are added to the price of the oligonucleotide.

| 5' Modification | 0.2 μmol / 3-5 OD_{260} | 1.0 μmol / 15 OD_{260} |
|-----------------|---|--|
| Fluorescein | 90 € | 127 € |
| FAM | 104 € | 139 € |
| TET | 121 € | 174 € |
| HEX | 121 € | 174 € |
| Cy3 | 193 € | 221 € |
| Tamra | 150 € | 204 € |
| Cy5 | 193 € | 221 € |
| Amino | 90 € | 127 € |
| Biotin | 76 € | 104 € |
| Phosphate | 38 € | 52 € |

6

Modified RNA Oligonucleotides - 3' Labels

Prices per modification are added to the price of the oligonucleotide.

| 3' Modification | 0.2 μmol / 3-5 OD_{260} | 1.0 μmol / 15 OD_{260} |
|--------------------------|---|--|
| Tamra | 119 € | 195 € |
| Amino | 64 € | 90 € |
| Biotin | 104 € | 159 € |
| Thiol | 127 € | 162 € |
| Deoxynucleotide / Base * | 7.90 € | 8.60 € |

* Price per deoxynucleotide. For 3' dinucleotide overhangs or 5' RNA-DNA 3' chimera molecules.

Shipping Charge

Shipping cost within Germany is 2 Euro. For order values over 250 Euro shipping is free.

If you wish to order from outside of Germany, please inquire for shipping costs.

For custom-made modifications, larger scales or further information, please contact us at oligos@jenabioscience.com