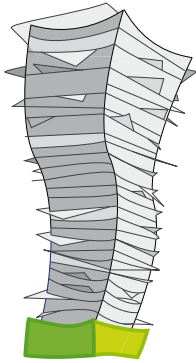
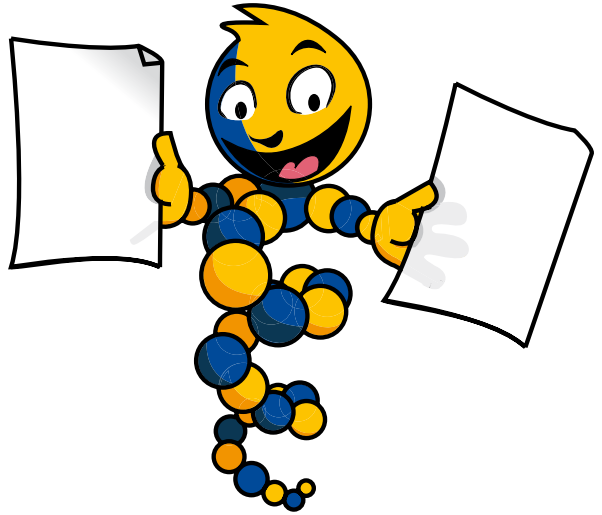


Oli&GO 1•2•3™



WITHOUT Oli&GO 1•2•3™



WITH Oli&GO 1•2•3™

MORE TIME FOR SCIENCE!

**FROM
MAY, 4th
2009**

 **Eurogentec**
Experience true partnership

Oli&GO 1·2·3™

Features

- > 1 Oli&GO integrated counter
- > 2 times more available modifications
- > 3 ranges of Oligonucleotides (Custom Oligonucleotides, Real-Time qPCR or RNAi Oligonucleotides)

Benefits

- > Less administrative burden
- > Bulk order prices
- > Realtime account visualization
- > From PCR primers to highly modified oligonucleotides including Real-Time qPCR probes and siRNA duplexes
- > Online tracking

In 2001, Eurogentec launched an exclusive and revolutionary ordering concept for your Oligonucleotides named Oli&GO Privilege Programme including Genomic and OliGold® Oligonucleotides, extended in 2007 by Oli&GO+ with Double-Dye Probes and siRNA Duplexes.

In 2009, Eurogentec has re-invented Oli&GO to bring you even more!

Oli&GO 1·2·3™ preserves all the specifications and benefits of the original Privilege Programmes, while being simpler than ever.



Oli&GO 1·2·3™ is managed by one and only one Oli&GO counter in Base Units (BU, Eurogentec's currency) according to:

1 BU = 1 DNA Base 10 nmol SePOP Desalted

With only 1 order and 1 invoice, you have access to a new user-friendly interface allowing you to order all your Oligonucleotides (even the most customised ones) whenever you want, even if your needs are spread over several months.

Once your sent order, the online tracking allows you to check the status of your orders at any time.

With Oli&GO 1·2·3™ , you will love ordering your oligonucleotides!

What Oligonucleotides?

Custom Oligonucleotides

- > PCR Primers
- > FISH Probes
- > Pyrosequencing Primers
- > Cloning Linkers
- > Antisense Oligonucleotides

RNAi Oligonucleotides

- > Custom siRNA Duplexes



Real-Time qPCR Oligonucleotides

- > Double-Dye Probes
- > Molecular Beacons
- > Scorpions® Probes
- > LC Hybridization Probes

How much are your Oligonucleotides in Base Units (BU)?

Oligo 1

Custom Oligonucleotide

20 bases 10 nmol $20 \times 1.00 = 20.00$ BU

Purification RP-Cartridge•Gold™

10 nmol 22.00 BU

Custom Oligonucleotides 42.00 BU

Oligo 2

Custom Oligonucleotide

30 bases 200 nmol $30 \times 2.20 = 66.00$ BU

Modification 5' Digoxigenin

200 nmol 490.40 BU

Purification RP-HPLC 200 nmol 180.00 BU

Custom Oligonucleotides 736.40 BU

Examples in Base Units (Catalogue Prices)

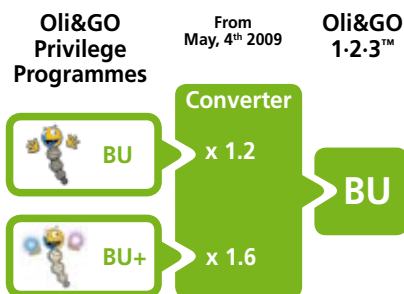
Oligo Type	Bases	Synthesis Scale	Modification(s)	Purification	Base Units (BU)
Custom Oligo	20	10 nmol	-	SePOP	20.00
Custom Oligo	20	40 nmol	-	SePOP	24.00
1 Custom Oligo	20	10 nmol	-	RP-Cartridge	42.00
Custom Oligo	30	200 nmol	5'FAM	RP-HPLC	366.00
2 Custom Oligo	30	200 nmol	Digoxigenin	RP-HPLC	736.40
Double-Dye Probe	25	10 nmol	5'FAM 3'TAMRA	RP-HPLC	460.00
Double-Dye Probe	30	40 nmol	5'FAM 3'BHQ-1™	RP-HPLC	760.00
siRNA Duplex	21 (19+2)	10 nmol	-	SePOP	540.00
siRNA Duplex	21 (19+2)	200 nmol	-	PAGE	1380.00

Those prices in BU may be lower according to dedicated additional discounts

References & Prices

Description	Reference	EUR	CHF	GBP
Oli&GO 1-2-3™ Activation				
444 BU	CD-OG001-0444	80.00	121.60	66.64
1234 BU	CD-OG001-1234	308.50	468.92	256.98
4444 BU	CD-OG001-4444	1111.00	1688.72	925.46
Oli&GO 1-2-3™ Reloads				
1234 BU	CD-OGR01-1234	308.50	468.92	256.98
4444 BU	CD-OGR01-4444	1111.00	1688.72	925.46
8888 BU	CD-OGR01-8888	2133.12	3242.34	1776.89
17776 BU	CD-OGR01-X776	4088.48	6214.49	3405.70

Already an oli&GO Privilege Programme Customer?



Your transition towards **Oli&GO 1-2-3™** will be automatic and totally transparent according to the conversion rules just beside.

From May, 4th 2009 and with the same login (8XXXXX), you will have access to the new **Oli&GO 1-2-3™** interface with only 1 integrated counter in Base Units (BU).

Welcome to **Oli&GO 1-2-3™**!

Eurogentec S.A.

Rue Bois St. Jean 5
4102 Seraing, Belgium
Tel: +32 (0)4 372 76 65
Fax: +49 (0)4 372 75 00
info@eurogentec.com
www.eurogentec.com