



ORDER FORM HEADER

MANDATORY
YOUR ORDER REFERENCE:

MENTION OBLIGATOIRE : VOTRE N° DE BON DE COMMANDE

DATE:

Page of

Our offer Nr.

Standing order - Bon Provisionnel - Guthaben: YES

Contact person:

Phone:

INVOICE ADDRESS

Institute: _____

Department/Lab.: _____

NAME: _____

Address: _____

Phone: _____ Fax: _____

E-mail: _____ VAT Nr. _____

VAT Exemption Nr. (UK only) _____

DELIVERY ADDRESS

Institute: _____

Department/Lab.: _____

NAME: _____

Address: _____

Phone: _____ Fax: _____

E-mail: _____

**For any question about ordering procedure, do not hesitate to contact us at:
+ 32 4 372 76 65 or info@eurogentec.com**

ORDERING FAX:

Benelux and Switzerland (french): + 32 (0)4 264 07 88
Germany, Austria and Switzerland (german): + 49 221 258 94 54

France: 02 41 73 10 26
UK: + 44 (0)1794 522 417



Oligonucleotide Order form

Order ref.:

Name:

Date:

Page of

OLIGO NAME _____ (Max 15 characters)

DNA RNA

5' _____ 3'

_____ 3	_____ 6	_____ 9	_____ 12	_____ 15	_____ 18	_____ 21	_____ 24	_____ 27	_____ 30	_____ 33	_____ 36
_____ 39	_____ 42	_____ 45	_____ 48	_____ 51	_____ 54	_____ 57	_____ 60	_____ 63	_____ 66	_____ 69	_____ 72
_____ 75	_____ 78	_____ 81	_____ 84	_____ 87	_____ 90	_____ 93	_____ 96	_____ 99	_____ 102	_____ 105	_____ 108

Purification : HPLC IE - HPLC PAGE UltraPureGold®

Comments:

5' modification :

3' modification :

Internal modification 1 (indicate with 1: AA1AA):

Internal modification 2 (indicate with 2: AA2AA):

OLIGO NAME _____ (Max 15 characters)

DNA RNA

5' _____ 3'

_____ 3	_____ 6	_____ 9	_____ 12	_____ 15	_____ 18	_____ 21	_____ 24	_____ 27	_____ 30	_____ 33	_____ 36
_____ 39	_____ 42	_____ 45	_____ 48	_____ 51	_____ 54	_____ 57	_____ 60	_____ 63	_____ 66	_____ 69	_____ 72
_____ 75	_____ 78	_____ 81	_____ 84	_____ 87	_____ 90	_____ 93	_____ 96	_____ 99	_____ 102	_____ 105	_____ 108

Purification : HPLC IE - HPLC PAGE UltraPureGold®

Comments:

5' modification :

3' modification :

Internal modification 1 (indicate with 1: AA1AA):

Internal modification 2 (indicate with 2: AA2AA):

ABBREVIATIONS : R = [A,g] - Y = [C,T] - M = [A,C] - K = [g,T] - S = [g,C] - H = [A,T,C] - B = [g,T,C] - D = [g,A,T] - N = [A,g,C,T] - V = [g,A,C] - X = [C,] - W = [A,T] - I = Inosine - U = Uracil - O = 5Me-dC - Q = 5Br-dU

5' Modifications	3' Modifications	Internal Modifications
None	None	None
5' A-LNA®	3' A-LNA®	5-iodo-dU
5' C-LNA®	3' C-LNA®	A-LNA®
5' dInosine	3' dInosine	Amino C6-dT
5' G-LNA®	3' dUracil	Biotin-dT
5' T-LNA®	3' G-LNA®	Biotin-TEG
6-FAM	3' T-LNA®	BODIPY®-FLX
Acridine	3' Wobble B	BODIPY®-TMR
Amine-C12	3' Wobble D	C2-Psoralen
Amine-C3	3' Wobble H	C6-Psoralen
Amine-C6	3' Wobble K	Cholesterol-TEG
Biotin-TEG	3' Wobble M	C-LNA®
BODIPY®-FLX	3' Wobble N	DABCYL
BODIPY®-TMR	3' Wobble R	DABCYL-dT
C2-Psoralen	3' Wobble S	Digoxigenin
C6-Psoralen	3' Wobble V	Dinitrophenol
Cascade Blue	3' Wobble W	dInosine
Coumarin	3' Wobble X	dR-Cholesterol
Cy®3	3' Wobble Y	dSpacer
Cy®5	3'dA blocked	Fluorescein
Cy®5.5	3'dC blocked	G-LNA
Digoxigenin	3'ddC	JOE
Dinitrophenol	3'dG blocked	Pyrene
dR Biotin	3'dT blocked	Rhodamine
Dragonfly Orange	5-Br-dU	Rhodamine-Green
dSpacer	5-iodo-dU	ROX
EDANS	Acridine	Spacer-C12
Fluorescein	Amine-C6	Spacer-C18
Glyceryl	BHQ-1®	Spacer-C3
HEX	BHQ-2®	Spacer-C8
JOE	BHQ-3®	Spacer-C9
Phosphate	Biotin-TEG	TAMRA
Pyrene	Bodipy-FLX	TET
Rhodamine	C2-Psoralen	Texas Red®
Rhodamine Green	C6-Psoralen	Thiol
Rhodamine-6G	Cascade Blue	T-LNA®
ROX	Cholesterol-TEG	
Staby™ Tag	Cy®3	
TAMRA	Cy®5	
TEG Cholesterol	DABCYL	
TET	Deep Dark Quencher 1	
Texas Red®	Deep Dark Quencher 2	
Thiol-C6	Digoxigenin	
Yakima Yellow®	Dinitrophenol	
	dSpacer	
	Eclipse Dark Quencher	
	Fluorescein	
	Glyceryl	
	HEX	
	JOE	
	Phosphate	
	Pyrene	
	Rhodamine	
	Rhodamine-6G	
	Rhodamine-Green	
	ROX	
	Spacer-C18	
	Spacer-C3	
	Spacer-C9	
	TAMRA	
	TET®	
	Texas Red®	
	Thiol-C3	