

OLIGO NAME	REF#	CHEMISTRY	BASES	
Neomycin Rev	UN-PR085-005	DNA	20	
SEQUENCE (5' → 3')				
5' AGG-TGA-GAT-GAC-AGG-AGA-TC 3'				
BACKBONE (BASES + LINKAGES)				
A	7 GC (%) ¹	50.0	Quantity (OD@260nm)	1.07
C	2 MW (g.mol ⁻¹) ¹	6255.2	(nmol)	5.0
G	8 Ext. coeff. (l.mol ⁻¹ .cm ⁻¹) ¹	213500	(µg)	31.3
T	3 Tm (°C) ¹	60.0		
Wobble	0			
¹ Theoretical values				
PURIFICATION(S)				
HPLC-RP				
QUALITY CONTROL(S)				
MS - CGE				Passed

DELIVERY CONDITION
Room Temperature
RECONSTITUTION
To make up to 100 µM concentration, add 50 µl of recommended buffer. To make up to 20 µM concentration, add 250 µl of recommended buffer. To make up to 5 µM concentration, add 1000 µl of recommended buffer.
LONG TERM STORAGE
-20 °C

COMMENTS	
General: Tm (°C) calculated with 50 mM NaCl and 0 % formamide (PCR conditions) Specific:	
TECHNICAL SUPPORT	
If you have any questions feel free to call our Oligo Centre	
EUROPE	NORTH AMERICA
☎ 00 800 666 00 123 (European toll free number), ✉ oligo@eurogentec.com	☎ +1 (510) 896-1350 ✉ oligo.na@egt-biotech.com
Kaneka Eurogentec S.A. Liège Science Park Rue Bois Saint-Jean 5 - 4102 SERAING BELGIUM Tel: +32(0)4 372 74 00 - Fax: +32(0)4 264 07 88 E-mail: info@eurogentec.com Web: www.eurogentec.com	EGT North America, Inc. 34801 Campus Drive - Fremont, CA 94555 Tel.: +1 (510) 896 -1350 * Fax: +1 (510) 791-9572 E-mail: ivd.na@egt-biotech.com * Web: www.egt-biotech.com
RPM Liège T.V.A.-(BE)-0427.348.346 - ING Belgique Bank - IBAN: BE86 3400 2118 6050 BIC: BBRUBEBB	