

OLIGO NAME	REF#	CHEMISTRY	BASES		
Anchored Oligo dT (22)	UN-PR020-005	DNA	22		
<b>SEQUENCE (5' → 3')</b>					
5' TTT-TTT-TTT-TTT-TTT-TTT-TTV-N 3'					
<b>BACKBONE (BASES + LINKAGES)</b>					
A	0	GC (%) <sup>1</sup>	5.3	Quantity (OD@260nm)	0.91
C	0	MW (g.mol <sup>-1</sup> ) <sup>1</sup>	6641.52	(nmol)	5.00
G	0	Ext. coeff. (l.mol <sup>-1</sup> .cm <sup>-1</sup> ) <sup>1</sup>	181800	(µg)	33.2
T	20	Tm (°C) <sup>1</sup>	31.1		
Wobbles	2				
<sup>1</sup> Theoretical values					
<b>PURIFICATION(S)</b>					
HPLC-RP					
<b>QUALITY CONTROL(S)</b>					
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**

☎ 00 800 666 00 123 (European toll free number),

 ✉ [oligo@eurogentec.com](mailto:oligo@eurogentec.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](mailto:info@eurogentec.com) Web: [www.eurogentec.com](http://www.eurogentec.com)

 RPM Liège T.V.A.-(BE)-0427.348.346 - ING Belgique Bank - IBAN: BE86 3400 2118 6050  
 BIC: BBRUBEBB

**NORTH AMERICA**

☎ +1 (510) 896-1350

 ✉ [oligo.na@egt-biotech.com](mailto:oligo.na@egt-biotech.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](mailto:ivd.na@egt-biotech.com) \* Web: [www.egt-biotech.com](http://www.egt-biotech.com)