

OLIGO NAME	CATALOG #	CHEMISTRY	BASES		
TET-T10	UN-CT020-005	DNA	10		
<b>SEQUENCE (5' → 3')</b>					
5' TET - TTT-TTT-TTT-T 3'					
<b>BACKBONE (BASES + LINKAGES)</b>					
A	0	GC (%) <sup>1</sup>	0.0	Quantity (OD@260nm)	0.49
C	0	MW (g.mol <sup>-1</sup> ) <sup>1</sup>	3655.3	(nmol)	5.0
G	0	Ext. coeff. (l.mol <sup>-1</sup> .cm <sup>-1</sup> ) <sup>1</sup>	97900	(µg)	18.3
T	10	Tm (°C) <sup>1</sup>	20.0		
Others	1				
<sup>1</sup> Theoretical values					
<b>PURIFICATION(S)</b>					
HPLC-RP					
<b>QUALITY CONTROL(S)</b>					
MS					
					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**

Store at -20 °C

**COMMENTS**

General: Tm % GC calculated with 50 mM NaCl and 0 % formamide (PCR conditions)

**TECHNICAL SUPPORT**

If you have any questions feel free to call our Oligo Centre

 ✉ [oligo@eurogentec.com](mailto:oligo@eurogentec.com)

 Kaneka Eurogentec S.A. Liège Science Park  
 Rue Bois Saint-Jean 5 - 4102 SERAINING 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