

OLIGO NAME	CATALOG #	CHEMISTRY	BASES		
ROX-T10	UN-CT035-005	DNA	10		
<b>SEQUENCE (5' → 3')</b>					
5' ROX™ - TTT-TTT-TTT-T 3'					
<b>BACKBONE (BASES + LINKAGES)</b>					
A	0	GC (%) <sup>1</sup>	0.0	Quantity (OD@260nm)	0.52
C	0	MW (g.mol <sup>-1</sup> ) <sup>1</sup>	3675.8	(nmol)	5.0
G	0	Ext. coeff. (l.mol <sup>-1</sup> .cm <sup>-1</sup> ) <sup>1</sup>	104200	(µg)	18.4
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 <span style="float: right;">Passed</span>					

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
✉ <a href="mailto:oligo@eurogentec.com">oligo@eurogentec.com</a>
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: <a href="mailto:info@eurogentec.com">info@eurogentec.com</a> Web: <a href="http://www.eurogentec.com">www.eurogentec.com</a>
RPM Liège T.V.A.-(BE)-0427.348.346 - ING Belgique Bank - IBAN: BE86 3400 2118 6050 BIC: BBRUBEBB