

OLIGO NAME	CATALOG #	CHEMISTRY	BASES		
AP® 5-T10	UN-CT001-005	DNA	10		
SEQUENCE (5' → 3')					
5' AP® 5 - TTT-TTT-TTT-T 3'					
BACKBONE (BASES + LINKAGES)					
A	0	GC (%) ¹	0.0	Quantity (OD@260nm)	0.61
C	0	MW (g.mol ⁻¹) ¹	3841.6	(nmol)	5.0
G	0	Ext. coeff. (l.mol ⁻¹ .cm ⁻¹) ¹	122600	(µg)	19.2
T	10	Tm (°C) ¹	20.0		
Others	1				
¹ Theoretical values					
PURIFICATION(S)					
HPLC-RP					
QUALITY CONTROL(S)					
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

 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 Web: www.eurogentec.com

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