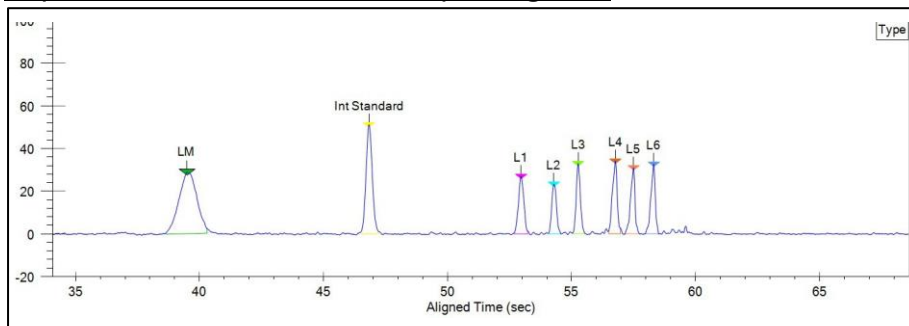


## HT pDNA Assay Reagent Kit

Product Number: CLS160450  
 Lot Number: DB19RK01  
 Expiration Date: 03 DEC 25

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Ladder Peak Height	Ladder	Height > 15 rfu	Pass
Ladder Peak Amount	Ladder	Max and Min Height < 30% from Median	Pass
Carry-over	TE Buffer	Impurities < 2.5 rfu	Pass
LOD	5 kb Linear DNA	Signal/Noise $\geq 3$	Pass
Resolution	Ladder	Resolution Scores > 1.5	Pass
SC Sizing Accuracy	SC 5.2kb	< 15%	Pass
Linear Sizing Accuracy	Revvity Standard 5.3kb	< 15%	Pass
Lower Marker Arrival Time	Ladder	Lower Marker (45-51 s),	Pass
Upper Marker Arrival Time	Ladder	Upper Marker (56-63 s)	Pass

### Representative Ladder Electropherogram\*



\* Peaks shown are typical pDNA ladder results. Peak height and width may vary.

Reagents	Vial	Quantity
Gel	Red ●	2 vial, 1.1 mL each
Dye	Blue ●	4 vials, 0.09 mL each
Sample Buffer	Bottle	1 Bottle, 30 mL each
Gel Dilution Buffer	Bottle	1 Bottle, 25 mL each
Ladder	Yellow ●	1 vial, 0.26 mL
Lower Marker	Green ●	1 vial, 1.5mL
Storage Buffer	White ○	9 vials, 1.8mL each

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

CLS160450-R Rev02

revvity