

FFA120 Reagent Kit

Product Number: CLS158081
 Lot Number: CA29RK01
 Expiration Date: 28NOV2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Lower marker migration Time	FFA120 ladder sample	Mean is within [19, 24] seconds	Pass
120 nt ladder migration time	FFA120 ladder sample	Mean is within [32, 38] seconds	Pass
Peak quality	FFA120 ladder sample	Mean of FWHM (nt) <=5 nt for all ladder samples	Pass
LM peak height	FFA120 ladder sample	Mean of LM peak height >= 50 for all ladder samples	Pass
Detection sensitivity	40 nt fragment, 20nM	Mean of peak height >= 5 RFU	Pass
Ladder quality	FFA120 ladder	All 5 peaks detected	Pass
Ladder purity	FFA120 ladder sample	No extra peak (with peak height repeatedly more than 5 RFU), and only five visible peaks	Pass

Reagent	Vial	Quantity
FFA120 Ladder	● Yellow	1 vial, 0.12 mL
FFA Chip Storage Buffer	○ White	5 vials, 1.8mL
FFA Gel Matrix	● Red	6 vials, 0.510mL
FFA Marker	● Green	1 vial, 0.8 mL
Consumable Items	PN 701672	1

Additional FFA120 Ladder can be ordered separately using Part Number CLS158127.

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.

CLS158081-R Rev01