

## FFA120 Ladder

Product Number: CLS158127  
 Lot Number: BE02RL01  
 Expiration Date: 03APR2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Ladder peak size	FFA120 ladder sample	Mean of each peak size meets the requirements	Pass
Ladder Height	FFA120 ladder sample	Mean within 30-300 RFU	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

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.

CLS158127-R Rev01