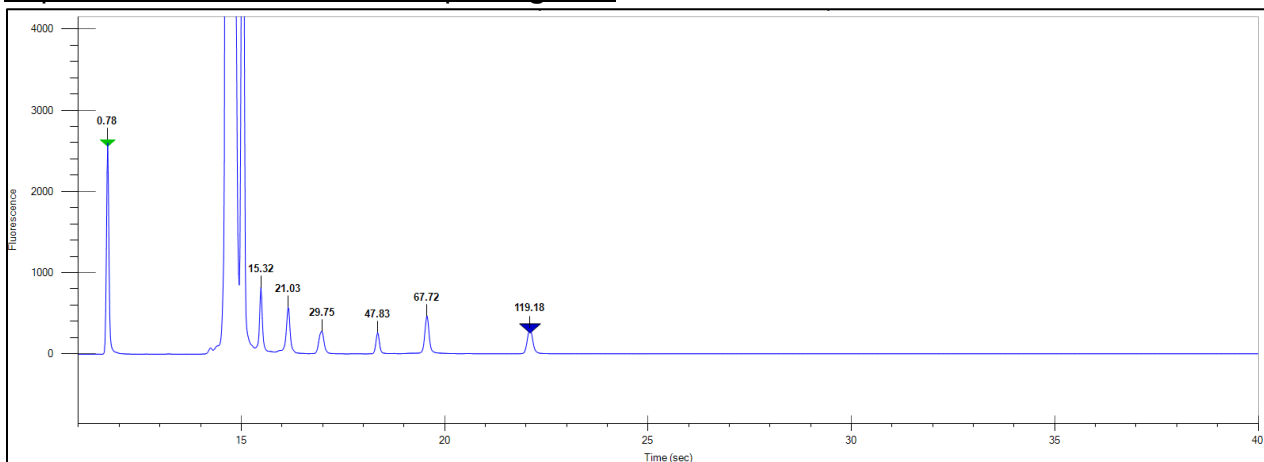


HT Pico Protein Express Reagent Kit

Product Number: 760498
 Lot Number: CJ03RK01
 Expiration Date: 03 JUN 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Resolution	Ladder	Minimum resolution value > 0.83	Pass
Purity	Antibody	Heavy chain/Light chain > 1.6	Pass
Sensitivity	CA	Signal/Noise ≥ 3	Pass
Sizing Accuracy	Ladder	Each peak +/- 20%	Pass
Impurities	Ladder	Ladder peak height after the 1 st ladder peak, $\leq 50\%$ before the 1 st ladder peak	Pass
Impurities	Ladder	Total impurities < 10% of total ladder peak heights	Pass
Peak to Valley ratio	NIST antibody	Ratio ≥ 1.01	Pass

Representative Ladder Electropherogram*



* Peaks shown are typical Protein Express ladder results. Peak height and width may vary.

Reagents	Vial	Quantity
Pico 5x Labeling Buffer	Clear ○	1 vial, 1.0mL each
Lyophilized Labeling Dye	Blue ●	4 vials
Sample Buffer	White ○	5 vials, 1.8mL each
Protein Gel Matrix	Red ●	2 vials, 1.8mL each
Protein Ladder	Yellow ●	1 vial, 0.08mL
Lower Marker	Green ●	1 vial, 0.5mL
Wash Buffer	Purple ●	4 vials, 1.8mL each
Stop Buffer	Orange ●	1 vial, 1mL
DMSO	Brown ●	1 vial, 1.0mL each
Consumable Items	PN 701553	1

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

760498-R Rev01

revvity