

## HT Protein QC Reagent Kit

Product Number: CLS960014  
 Lot Number: CC20RK01  
 Expiration Date: 19 DEC 2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Resolution	All Ladders	5% resolution $\geq 0.83$	Pass
Peak height	Lower Marker	LM > Ladder peaks (Ratio $\geq 1$ )	Pass
Sensitivity	5 ng/ul CA samples	All detected (Min height $\geq 5$ rfu)	Pass
Sizing	All ladder	All peaks within 15%	Pass
Largest Impurities	New Ladder	$\leq 20\%$ of smallest peak	Pass
Total Impurities	New Ladder	$\leq 10\%$ of total % purity	Pass
Standard Calibration	VeriMAb™ Standard	Final calibration range: 9-11%	Pass
Standard Precision	VeriMAb™ Standard	CV of minor peak: <5%	Pass

Reagent	Vial	Quantity
Protein QC Dye Solution	Blue ●	1 vial of 0.13 mL
Sample Buffer	White ○	7 vials, 1.5 mL each
Protein Gel Matrix	Red ●	3 vials, 1.7 mL each
Protein QC Ladder	Yellow ●	1 vial, 0.11 mL
Lower Marker	Green ●	2 vials, 0.5 mL each
Wash Buffer	Purple ●	3 vials, 1.8 mL each
VeriMAb™ Standard	Orange ●	1 vial, 0.04mL
Consumable Items	PN 960008	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.

CLS960014-R Rev01