

## HT Low MW Protein Express Reagent Kit

Product Number: 760573  
 Lot Number: DC04RK01  
 Expiration Date: 03 DEC 2025

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Resolution	All Ladder	Min Resolution (20% < 7kDa, 10% > 7kDa) $\geq$ 83%	Pass
Peak Height	Lower Marker	LM > Ladder peaks (Ratio $\geq$ 1)	Pass
Migration Time	All Ladder	Migration Time of all Ladder samples within spec	Pass
Sensitivity	30ng/uL CA Samples	All detected (Min Height $\geq$ 5rfu)	Pass
Sizing	All Ladder	All peaks within 20%	Pass
Concentration	125 ng/ul CA Samples	Within 30%	Pass
Largest Impurities	New Ladder	$\leq$ 20% of smallest pk	Pass
Total Impurities	New Ladder	$\leq$ 10% of total % purity	Pass
Stability	New Ladder	Total % purity of 6.5kDa Peak of new ladder > 12%	Pass

Reagents	Vial	Quantity
Protein Dye Solution	Blue ●	1 vial of 0.090 mL
Sample Buffer	White ○	3 vials, 1.4 mL each
Low MW Protein Gel Matrix	Red ●	2 vials, 1.8 mL each
Low MW Protein Ladder	Yellow ●	1 vial, 0.08 mL
Lower Marker	Green ●	1 vial, 0.5 mL
Wash Buffer	Purple ●	4 vials, 1.8 mL 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.

760573-R Rev01