

HT Low MW Protein Express Reagent Kit

Product Number: 760573
 Lot Number: BK16RK01
 Expiration Date: 15Jul2024

| PARAMETER | SAMPLE | SPECIFICATION | RESULT |
|--------------------|----------------------|---|--------|
| Resolution | All Ladder | Min Resolution (20% < 7kDa, 10% > 7kDa) ≥ 83% | Pass |
| Peak Height | Lower Marker | LM > Ladder peaks (Ratio ≥ 1) | Pass |
| Migration Time | All Ladder | Migration Time of all Ladder samples within spec | Pass |
| Sensitivity | 30ng/uL CA Samples | All detected (Min Height ≥ 5rfu) | Pass |
| Sizing | All Ladder | All peaks within 20% | Pass |
| Concentration | 125 ng/uL CA Samples | Within 30% | Pass |
| Largest Impurities | New Ladder | ≤ 20% of smallest pk | Pass |
| Total Impurities | New Ladder | ≤ 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