






HT Glycan Reagent Kit

Product Number: 760525
Lot Number: DD02RK01
Expiration Date: 01 Jan 2026

| PARAMETER | DESCRIPTION | SPECIFICATION | RESULT |
|---------------------------------|--|---|--------|
| Marker/Ladder Peak Height Ratio | Marker/tallest ladder peak height in new ladder | 3 to 5 | Pass |
| Marker Purity | No extra peak on new marker | (<2s from LM) or $\geq 50\%$ LM height (>2s from LM) and $\geq 10\text{rfu}$ (after LM) | Pass |
| Marker Height | Average marker peak height | Average $\leq 6000\text{rfu}$ | Pass |
| New Ladder Peak Height | Last ladder peak height is called | $\geq 100\text{rfu}$ | Pass |
| Ladder Purity | Smallest ladder peak/largest impurity peak of new ladder | Ratio > 5 | Pass |
| LM Migration Time | New ladder | Meets time spec | Pass |
| Last Ladder Peak Migration Time | New ladder | Meets time spec | Pass |
| Size % CV | All ladder peaks | $\leq 2.5\%$ | Pass |

| Reagents | Label | Quantity |
|--------------------------|--|----------------------|
| HT Glycan Ladder Diluent | Purple  | 1 vial, 0.75 mL |
| HT Glycan Marker Diluent | White  | 1 vials, 0.65 mL |
| HT Glycan Gel Matrix | Red  | 2 vials, 1.8 mL each |
| HT Glycan Ladder | Yellow  | 4 vials, 0.3 uG each |
| HT Glycan Marker | Green  | 4 vials, .6 uG 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.

760525-R Rev01