

HT Glycan Reagent Kit

Product Number: 760525
 Lot Number: CF24RK01
 Expiration Date: 23 MAR 2025

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