

Glycan Release and Labeling Kit

Product Number: CLS155434
 Lot Number: CA24RK01
 Expiration Date: 23OCT2024

PARAMETER	DESCRIPTION	SPECIFICATION	RESULT
Dye Purity	No extra peak on new dye blank	Meets Spec	Pass
Dye Reactivity	Signal-to-Noise of G0f samples	Average S/N > 6	Pass

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Digestion Limits	% Impurity of NGHC in 9.5 mg/ml antiTNF + enzyme sample	>95%	Pass
Complete Digestion	% Impurity of NGHC in 7.5 mg/ml antiTNF + enzyme sample	= 100%	Pass

Glycan Release and Labeling Kit Contents	Quantity
Denaturing Reagent (4 x 0.15 mL) for denaturing glycoprotein	1
PNGase F Reagent (4 x 0.15 mL) for releasing the oligosaccharides.	1
Dye Reagent (4 x 0.15 mL) for labeling the oligosaccharides	1
Store at -20°C	

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.

CLS155434-R Rev01