

HT Protein Charge Variant Kit

Product Number: CLS760670
 Lot Number: BM15RK01
 Expiration Date: 14SEPT2024

PARAMETER	SAMPLE	SPECIFICATION	RESULT
Histidine Migration	Average value of histidine sample wells	Meets spec	Pass
Lysine Migration	Average value of lysine sample wells	Meets spec	Pass
Dye Migration Time	Average value of all sample wells	Meets spec	Pass
His/Dye Peak Height Ratio	Average value of histidine sample wells	≥ 4.7	Pass
Lys/Dye Peak Height Ratio	Average value of lysine sample wells	≥ 36.9	Pass

Items at -20 °C (HT Protein Charge Variant Labeling Kit)	Vial	Quantity
Labeling Buffer	Orange ●	1 vial, 1 mL
Dye Concentrate	Blue ●	5 vials, 5 µL (24 samples / vial)

Items at 4 °C (HT Protein Charge Variant Buffer Kit)	Bottle	Quantity
Running Buffer pH 5.6	Red ●	1 bottle, 25 mL
Running Buffer pH 7.2	Green ●	1 bottle, 25 mL
Storage Buffer	Purple ●	1 bottle, 25 mL
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.

CLS760670-R Rev01