

Version: 03

Update: 03/12/2024

eStain L1C Protein Staining Kit

Cat.No. L00753

Contents

I. Product Description	1
II. Instructions for use	1
III. Storage Conditions	3
IV. Ordering Information	4

I Product description

eStain L1C Staining Kit (Cat No: L00753) contains reagents that are used with eStain L1 protein staining device (L00657). Each kit contains 1 x 2 L of staining solution, 1 x 2 L of destaining solution, and a pack of filter paper (50 pK).

Table 1. Kit components

Component name	Size	Quantity
Concentrated Staining Solution	2 L	1
Concentrated Destaining Solution	2 L	1
eStain Filter Paper	50 pK	1

II Instructions for use

For shorter time program

1、 Staining Solution

Please dilute the concentrated staining solution as below. Mix well before use.

Table 2. Staining Buffer Preparation for Shorter Time Program

Step	Volume	Total Volume
1	Prepare a clean 5 L empty bottle.	0
2	Add 2 L Concentrated Staining Solution into 5L bottle.	2 L
3	Add 1 L Isopropanol* into step 2 bottle and mix well.	3 L
4	Add 2 L ddH ₂ O into step 3 bottle and mix well	5 L

* Isopropanol must be ACS reagent, ≥99.5%.

2、Destaining Solution

Please dilute the concentrated destaining solution as below. Mix well before use.

Table 3. Destaining Buffer Preparation for Shorter Time Program

Step	Volume	Total Volume
1	Prepare a clean 5 L empty bottle	0
2	Add 1 L Concentrated Destaining Solution into 5 L bottle.	1 L
3	Add 4 L ddH ₂ O into step 2 bottle and mix well.	5 L

3、Program Setting

Refer to eStain L1 Protein Staining System manual for setting new program

1. Press UP or DOWN to select the channel to highlight
2. Press and hold down SETTING until it beeps to enter the program interface
3. Select the method to use or switch by pressing UP or DOWN
4. When finished, press SETTING and return to the Working Interface
(To modify, press and hold SETTING until beeps to enter the Parameters Interface)
5. Modify the program as following parameter

Stain	Equilibrium cycles	0	times
	Staining cycles	1	times
	cycle 1	3.0	min
	Destaining cycles	2	times
	cycle 1	2.0	min
	cycle 2	3.0	min
UP & DOWN to modify cycle or time		Press setting to apply method	
Press and hold setting to save and quit			

For lower background program

1、Staining Solution

Please dilute the concentrated staining solution as below. Mix well before use.

Table 4. Staining Buffer Preparation for Shorter Time Program

Step	Volume	Total Volume
1	Prepare a clean 5 L empty bottle.	0
2	Add 2 L Concentrated Staining Solution into 5 L bottle.	2 L
3	Add 1.5 L Isopropanol* into step 2 bottle and mix well.	3.5 L
4	Add 1.5 L ddH ₂ O into step 3 bottle and mix well	5 L

* Isopropanol must be ACS reagent, $\geq 99.5\%$.

2、Destaining Solution

Please dilute the concentrated destaining solution as below. Mix well before use.

Table 5. Destaining Buffer Preparation for Shorter Time Program

Step	Volume	Total Volume
1	Prepare a clean 5 L empty bottle	0
2	Add 1 L Concentrated Destaining Solution into 5 L bottle.	1 L
3	Add 3 L ddH ₂ O into step 2 bottle and mix well.	4 L

3、Program Setting

Refer to eStain L1 Protein Staining System manual for setting new program

1. Press UP or DOWN to select the channel to highlight
2. Press and hold down SETTING until it beeps to enter the program interface
3. Select the method to use or switch by pressing UP or DOWN
4. When finished, press SETTING and return to the Working Interface
(To modify, press and hold SETTING until beeps to enter the Parameters Interface)
5. Modify the program as following parameter

Stain	Parameter	Value	Unit
	Equilibrium cycles	0	times
	Staining cycles	1	times
	cycle 1	5.0	min
	Destaining cycles	2	times
	cycle 1	3.0	min
	cycle 2	3.0	min

UP & DOWN to modify cycle or time Press setting to apply method
 Press and hold setting to save and quit

III Storage condition

Keep the staining kit in well-ventilated area and away from direct light. The staining kit is stable for up to 12 months if stored at room temperature.

IV Ordering Information

Product Name	Cat No.
eStain L1C Staining Kit	L00753
eStain L1 Protein Staining Device	L00657

For research use only. Not intended for human or animal clinical trials, therapeutic or diagnostic use.

生产商：南京金斯瑞生物科技有限公司 江苏省南京市江宁区科学园雍熙路 28 号

Manufacturer: Nanjing GenScript Biotech Co., Ltd. No. 28 Yongxi Road, Jiangning District, Nanjing, Jiangsu, China