



## 신제품 연구용 TFF 필터, Vivaflow SU 지금 특별 할인으로 만나보세요 !



기간 : 2025. 04. 21 ~ 2025. 08. 31

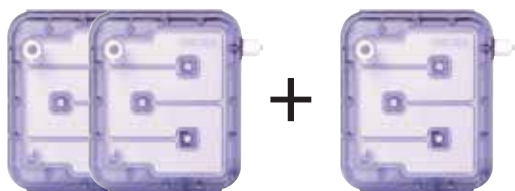
대상 : Vivaflow SU 구매 고객 \*EVENT1, 2 중 택1



프로모션 신청하기 ▲

### EVENT 1

첫 구매 시,  
Vivaflow SU  
2+1 증정



2box 구입 시, 1box 추가 증정 !  
(box당 2개입)

### EVENT 2

패키지 구매 시,  
펌프와 필터  
모두 30% 할인



패키지 구성 (펌프+필터)

## Ordering Information

## Vivaflow® SU Cassettes

| Membrane Material                     | MWCO      | Recommended Feed Volume |                 |
|---------------------------------------|-----------|-------------------------|-----------------|
|                                       |           | 100 – 500 mL            | 500 – 1,000 mL  |
| Hydrosart® regenerated cellulose (RC) | 2 kDa     | VF-S050H0002-IV         | VF-S050H0002-SV |
|                                       | 5 kDa     | VF-S050H0005-IV         | VF-S050H0005-SV |
|                                       | 10 kDa    | VF-S050H0010-IV         | VF-S050H0010-SV |
|                                       | 30 kDa    | VF-S050H0030-IV         | VF-S050H0030-SV |
|                                       | 100 kDa   | VF-S050H0100-IV         | VF-S050H0100-SV |
|                                       | 300 kDa   | VF-S050H0300-IV         | VF-S050H0300-SV |
| Polyethersulfone (PES)                | 5 kDa     | VF-S050P0005-IV         | VF-S050P0005-SV |
|                                       | 10 kDa    | VF-S050P0010-IV         | VF-S050P0010-SV |
|                                       | 30 kDa    | VF-S050P0030-IV         | VF-S050P0030-SV |
|                                       | 50 kDa    | VF-S050P0050-IV         | VF-S050P0050-SV |
|                                       | 100 kDa   | VF-S050P0100-IV         | VF-S050P0100-SV |
|                                       | 300 kDa   | VF-S050P0300-IV         | VF-S050P0300-SV |
|                                       | 1,000 kDa | VF-S050P1000-IV         | VF-S050P1000-SV |
|                                       | 0.2 µm    | VF-S050P2000-IV         | VF-S050P2000-SV |

## Peristaltic Pump

| Description                                  | Package Contents                    | Order No.    |
|----------------------------------------------|-------------------------------------|--------------|
| Peristaltic Pump, 100-120/200-240V           | 1 unit <sup>5</sup><br>1 user guide | VF-APD0001-1 |
| Peristaltic Pump Head for 1.6 mm W.T. tubing | 1 unit<br>1 user guide              | VF-APH0001-1 |
| Peristaltic Pump Head for 2.4 mm W.T. tubing | 1 unit<br>1 user guide              | VF-APH0002-1 |

<sup>5</sup> Includes power cable with region-specific plug.

