



Hi-Flow™ Plus HF120 Membrane

GENERAL INFORMATION

HF12002 and HF12004 are Hi-Flow™ Plus Cellulose Ester membranes with nominal capillary flow time of 120 seconds/4 cm and nominal membrane thickness of 135 µm on 50 µm (2 mil) or 100 µm (4 mil) polyester backing. Hi-Flow™ Plus membranes are manufactured to create a superior surface morphology that allows for the consistent and even flow distribution of reagents across the membrane surface. This excellent surface quality improves reagent addition and the clarity of the test results.

MATERIALS OF CONSTRUCTION

Membrane is made of nitrocellulose.

DESIGN PROPERTIES

Visual Quality

Membrane is white in color with no visual defects visible under normal lighting.

Capillary Flow Time (sec/4 cm) membrane coupon size = 1.5 cm x 4.0 cm

Master roll average:	120 +/- 12 seconds
Individual values:	120 +/- 30 seconds
Composite membrane thickness with 2 mil backing:	185 µm +/- 19 µm
Composite membrane thickness with 4 mil backing:	235 µm +/- 19 µm

Protein Binding (µg/cm²)

Insulin:	>30
BSA:	>45
Goat IgG:	>95
Length:	100 meters (328 feet)
Format:	Roll

Hi Flow™ Plus membranes are available in the standard slit widths, however custom widths and sheet sizes are also available.

Slit width (cm)	Slit rolls/Master	Linear feet/100 meters	Slit width (cm)	Slit rolls/Master
1.60	18	5904	3.175	9
1.70	17	5576	3.20	9
1.80	16	5248	3.50	8
2.00	14	4592	3.81 (1.5 in)	7
2.20	13	4264	4.00	7
2.25	13	4264	4.5	6
2.50	11	3608	4.80	6
2.54 (1 in)	11	3608	5.00	5
2.80	10	3280	8.00	3
3.00	9	2952	13.20	2

HF12002XSS (2mil backed) and HF12004XSS (4mil backed) are 5 sheet packs of membrane; 30.5 cm x 25 cm

SHF1200225 (2 mil backed) and SHF1200425 (4 mil backed) are slit sample rolls; 25 mm wide x 100 meters long



To place an order or receive technical assistance

In the U.S. and Canada, call toll-free
1-800-645-5476

For other countries across Europe
and the world, please visit:
[EMDMillipore.com/offices](https://www.EMDMillipore.com/offices)

For Technical Service, please visit:
[EMDMillipore.com/techservice](https://www.EMDMillipore.com/techservice)

[EMDMillipore.com](https://www.EMDMillipore.com)

MilliporeSigma
400 Summit Drive
Burlington, MA 01803

