



Hi-Flow™ Plus HF075 Membrane

GENERAL INFORMATION

HF07502 & HF07504 are Hi-Flow™ Plus Lateral Flow membranes with nominal capillary flow time of 75 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 cast on a polyester backing.

DESIGN PROPERTIES

Visual Quality

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

Capillary Flow Time (sec/4 cm)

Master roll average:	75 +/- 8 seconds
Individual values:	75 +/- 19 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

HF07502XSS (2 mil backed) and HF07504XSS (4 mil backed) are 5 sheet packs of membranes; 30.5 cm x 25 cm

SHF0750225 (2 mil backed) and SHF0750425 (4 mil backed) are slit sample rolls; 25 mm wide x 100 meters long each.



To place an order or receive technical assistance

In Europe, please call Customer Service:

France: 0825 045 645

Spain: 901 516 645 Option 1

Germany: 069 86798021

Switzerland: 0848 645 645

Italy: 848 845 645

United Kingdom: 0870 900 4645

For other countries across Europe, please call: +44 (0) 115 943 0840

Or visit: [MerckMillipore.com/offices](https://www.MerckMillipore.com/offices)

For Technical Service visit: [MerckMillipore.com/techservice](https://www.MerckMillipore.com/techservice)

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

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

