



Hi-Flow™ Plus HF135 Membrane

GENERAL INFORMATION

HF13502 and HF13504 are Hi-Flow™ Plus Cellulose Ester membranes with nominal capillary flow time of 135 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)

Master roll average: 135 +/- 15 seconds

Individual values: 135 +/- 34 seconds

Thickness 2 mil backing: 185 µm +/- 19 µm

Thickness 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 standard slit widths, however custom widths and sheet sizes are also available.

Slit width (cm)	Slit rolls/ Master	Linear feet/ 100 meters
1.60	18	5904
1.70	17	5576
1.80	16	5248
2.00	14	4592
2.20	13	4264
2.25	13	4264
2.50	11	3608
2.54 (1 in)	11	3608
2.80	10	3280
3.00	9	2952

Slit width (cm)	Slit rolls/ Master	Linear feet/ 100 meter
3.175	9	2952
3.20	9	2952
3.50	8	2624
3.81 (1.5 in)	7	2296
4.00	7	2296
4.5	6	1968
4.80	6	1968
5.00	5	1640
8.00	3	984
13.20	2	656



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

