

## LabSorb Polyethylene-Coated Paper for Surface Protection



### Filter Papers

The LabSorb grade is a highly absorptive paper coated on one side with polyethylene. Used with the cellulose side up, the paper absorbs liquids, which are stopped by their polyethylene layer and thus prevented from soaking through. Used with the polyethylene side up, the paper is highly useful for recover valuable or toxic liquids.

# Applications

- Prevent radioactive contamination of work surfaces (in radiochemical laboratories)
- Recover spilled solutions containing expensive reagents
- Protect laboratory bench surfaces from spillage or splashes of liquids by preventing absorption and seepage of these liquids through work surfaces
- Line animal cages for protection and improvement of hygiene
- Protect objects from falling on hard surfaces to reduce impact

# Properties

- Weight: 140 g/m<sup>2</sup>
- Water capacity: 150 %

# Standard Sizes

LabSorb

Part number	Dimensions	Quantity per pack
FT-1-601-400050	400 mm × 50 m	1 roll
FT-1-601-400100	400 mm × 100 m	1 roll
FT-1-601-460050	460 mm × 50 m	1 roll
FT-1-601-600050	600 mm × 50 m	1 roll
FT-1-601-600100	600 mm × 100 m	1 roll
FT-2-601-460570K	460 mm × 570 mm	50 sheets
FT-2-601-480600K	480 mm × 600 mm	50 sheets



## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Straße 20  
37079 Göttingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 631 254 4249  
Toll-free +1 800 635 2906

For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)