

# iTLC-SG

## Product Information Sheet



**Agilent**

Trusted Answers

### Definition

Instant Thin Layer Chromatography Medium (iTLC) is a binderless, glass microfiber chromatography paper or sheet impregnated with a silica gel. It is a porous, slightly acidic sheet that provides excellent resolution.

### Features and use

- More convenient with faster developing times than traditional thin layer chromatography plates.
- Open porous matrix provides fast developing time (often less than 20 minutes).
- Ideal for evaluating radioisotope purity.
- Can be used to separate lipids and other nonpolar compounds, and is comparable to many other silica gel plates.
- Sample elution is easy and without interference from organic binders.
- Can be cut with scissors or a cutter. Use caution as it is a fragile glass-type material. Handle the product with care.

### Instructions for use

1. Place iTLC-SG sheets in a 110 °C oven.
2. Bake for 30 minutes.
3. Remove from oven, and:
  - Use immediately
  - Store in a desiccator over dry silica gel or calcium chloride at 15 to 30 °C (59 to 86 °F)

This removes unbound water and activates the binding sites on the silica gel, which results in well separated, tight spots. The medium should be used as soon as it is removed from the oven or the desiccator. Exposure to a humid atmosphere may result in poor or irreproducible separation due to the medium adsorbing water from the air.

### Typical QC data

Includes statistical data regarding relative separation of several dye marker ( $R_f$ s or retention factor), run times for nonpolar organic developing solution.

### Ordering information

Part number	Description	Packaging
SGI0001	iTLC-SG (4.5 × 12 inch sheets)	50/pk

[www.agilent.com/chem](http://www.agilent.com/chem)

This information is subject to change without notice.