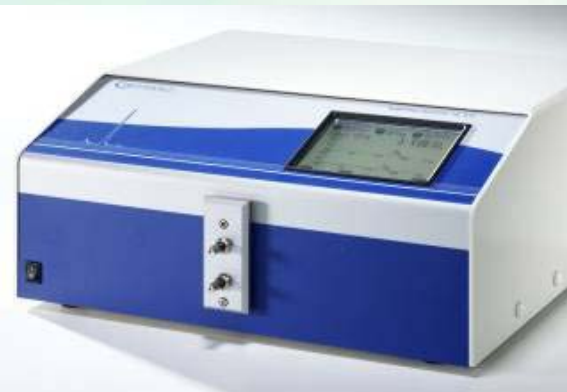


## FlowStar LB 513

Radioactivity flow-through detector for HPLC



### Touch screen for easy operation

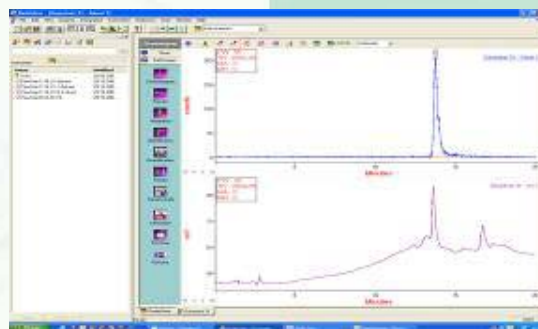


The FlowStar represents the latest enhancements in flow-through detectors for Radio HPLC. It combines the latest technology such as touch screen for easy operation and the highest sensitivity.

[Learn more about FlowStar](#)

### RadioStar software

FlowStar can be controlled by the RadioStar software. A control library to implement FlowStar into other software packages is available on request. When operated in stand alone mode the instrument can be interfaced with all available HPLC systems.



### Measuring cells with ID chip



A wide range of highly sensitive measuring cells for all types of applications has been developed. All cells are equipped with an internal cell ID chip to store cell related data.

[Get in touch with Berthold Technologies](#)

[www.Berthold.com/Bio](http://www.Berthold.com/Bio)