S/RICS

Arium® Pro

Application-orientated and flexible to meet the highest demands

- Modular System selection specifically for your application
- Flexible Perfect integration into any laboratory
- Easy to use Display with touch function and intuitive menu
- Fast Favorites function with direct access for recurring volumes
- Compatible with Arium[®] Smart Station

Simplifying Progress

