

ambr® 250 high throughput

