



Your lab,
Your reputation,
Your choice.

Bio-Rad **A1c**

BE THE DIFFERENCE

BIO-RAD



Working together for better patient care

At Bio-Rad, we recognize your need to balance high quality care with the operational demands of your laboratory. That's why we're always looking for ways to go the extra mile with support and services designed to keep your office, laboratory or clinic moving forward—so you can focus on being the difference through quality patient care.

Designed for unmatched performance

Bio-Rad A1c solutions deliver the unmatched performance and quality results you need to deliver the best patient care possible. When you use equipment from Bio-Rad, you can be confident that you're doing your best to help healthcare professionals treat their patients more effectively.

Quality results that fit your needs

Because every patient deserves accurate, reliable results, Bio-Rad has developed solutions for any workload and testing environment.

Today's most demanding laboratories can turn to Bio-Rad A1c assays for a range of fully automated testing solutions, whether they're running a few samples a day or hundreds. We offer analyzers for point of care settings, where immediate results are required, as well as hospitals and reference labs where turnaround times are crucial.

No matter what you require, count on Bio-Rad A1c to help you be the difference for the people who depend on your results.

Depend on Bio-Rad A1c for:

- Quality and innovation backed by 30 years of expertise
- Trusted leadership in the industry
- A broad, comprehensive line of A1c products to serve any environment
- Dedicated customer service that ensures support whenever you need it

Bio-Rad A1c

Be the difference in quality patient care with precise A1c solutions designed for your specific testing needs.

Helping you be the difference

Quality patient care is at the heart of what you do—and you choose your A1c assays accordingly. At Bio-Rad, we share this commitment to excellence and are dedicated to helping you be the difference.

Like you, we believe that what some consider “acceptable” margins of error can make a big difference in the level of care a patient receives. That's why we've never settled for “good enough” since pioneering A1c testing over 30 years ago. We're constantly working to improve our line of A1c testing solutions—so you can always look to Bio-Rad A1c for the laboratory results you need to be the difference for the people who count on you most.



