



CELEBRATING 20 YEARS OF
PARTNERSHIP • INNOVATION • EXCELLENCE

A Letter From Our CEO

Dear Team, Customers, and Partners,

Twenty years ago, we started this company with a clear goal in mind: to advance assay and test technologies (AAT). Two decades later, those three letters have grown into something far beyond my expectations. That is not to say the journey has been without challenges; quite the contrary. But time and again, we have risen to the occasion together, guided by our passion, ingenuity, and mutual respect.

To my esteemed colleagues and fearless team members, thank you for making these past twenty years such a meaningful and rewarding journey. Your dedication, creativity, and teamwork have shaped AAT into what it is today. It has been a true pleasure working alongside all of you, and I look forward with great excitement to the discoveries still to come.

To our valued customers and partners, thank you for your trust, support, and collaboration over the years. We are deeply grateful for the role you have played in our journey, and we remain committed to

serving you with the same respect, innovation, and transparency that have guided us from the beginning.

The past twenty years have been filled with bright moments, meaningful achievements, and lasting partnerships. Yet I believe our brightest days still lie ahead. With gratitude for all we have accomplished together, and excitement for all that is still to come, let us continue to move forward with purpose, curiosity, and shared commitment. Together We Shine!

Sincerely,



Zhenjun Diwu, Ph.D.
CEO, AAT Bioquest, Inc.



Our Story

At AAT Bioquest, we develop the tools that help researchers solve the most challenging problems through innovation.

Started in 2006 by three passionate scientists, the company has since grown exponentially. Today, we distribute to over 60 countries across the globe, while fielding a catalog of over 3500 products. The secret to our success? The dedicated and talented individuals who have come to form the AAT family.

Our team brings together deep scientific expertise, creativity, and a shared vision of pushing the frontiers of life science research. Through continuous investment in research and development, and more importantly, our people, we have been able to stay at the forefront of bioreagent technologies, designing high-quality assay kits, fluorescent probes, biochemicals, and custom solutions that help scientists work more efficiently and achieve more reliable results. From academic laboratories to biotechnology and pharmaceutical companies, we are proud to support discoveries that move science forward.



MISSION

To empower scientists with bright tools through sustainable innovation

VISION

To unlock the frontiers of life science research, drug discovery, and in vitro diagnostics

PHILOSOPHY

Mutual respect fosters a community where passion and ingenuity can thrive

Core Values

01 | INNOVATION

We advance life science research by leveraging our expertise to develop reliable, high-quality tools that help scientists solve complex problems with confidence.

02 | PARTNERSHIP

We seek to build lasting relationships with our colleagues, customers, and community through respect, trust, responsiveness, and a shared commitment to helping each other succeed.

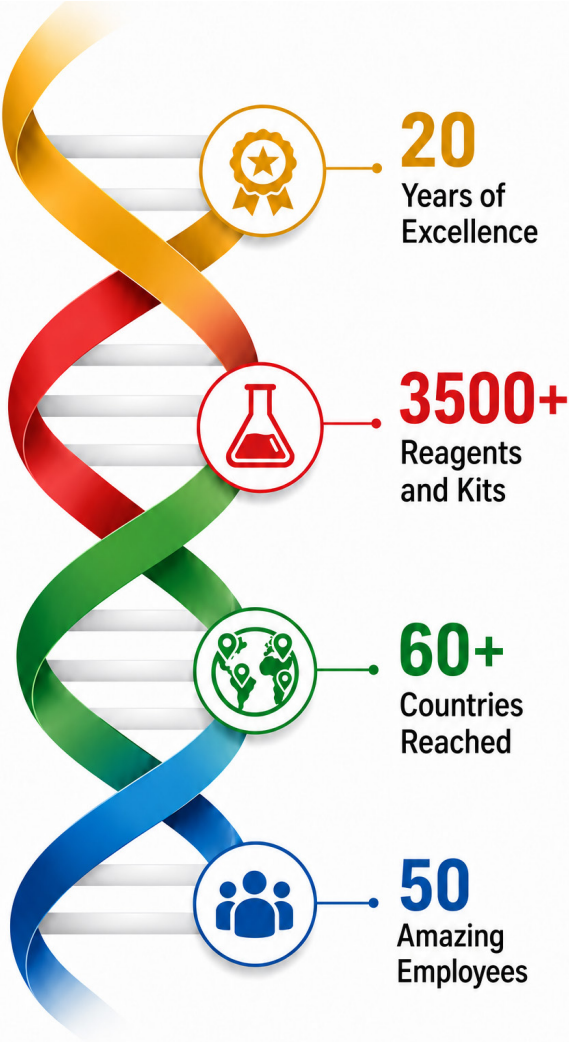
03 | EXCELLENCE

We hold ourselves to the highest standards in product development, service, and support so researchers can achieve dependable results.

04 | SUSTAINABILITY

We pursue growth with responsibility, supporting scientific discovery while respecting our people, our communities, and the world we share.

Key Achievements

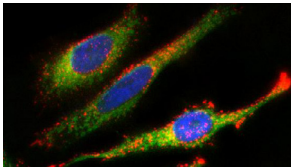


A History of Innovation

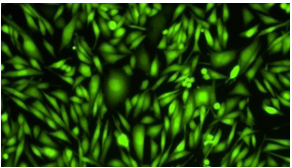


2006
AAT Bioquest is founded by a trio of passionate scientists looking to advance assay and test technologies.

2008
AAT Bioquest files its first patent. It is for **Fluo-8®**, a novel calcium indicator that unlocks high-throughput screening of GPCRs for drug discovery by enabling room temperature cell loading. It is sold as kits under the **Screen Quest™** brand.



2009
AAT Bioquest enters the fluorescence microscopy space with the first of several innovative dye families - **iFluor®** - which offers brighter and more photostable signals. **Cell Navigator®** kits are also released, simplifying cell and subcellular imaging.



2012
Continued R&D efforts lead to the patented **Cal-520®**, a next-generation calcium indicator that enables probenecid-free experimentation while exceeding previously achievable S/R ratios.

2007
AAT Bioquest trademarks its first series of assays for biochemical detection in microplate readers, sold under the brand, **Amplite®**.



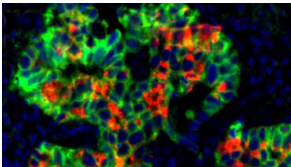
2010
AAT Bioquest enters the flow cytometry space with a family of long Stokes shift dyes for multiplexing, sold under the **mFluor™** brand name.

2011
The addition of **ReadiLink™** kits demonstrates AAT Bioquest's position as not only a leader in fluorescent dyes but also in conjugation technology.

2013
Dissatisfied with the poor reporting of CFSE, AAT Bioquest develops the **CytoTell®** dyes to properly "tell" the story of cell proliferation on flow cytometry and fluorescence microscopy platforms.

2017
With the release of **Calbryte™**, the brightest calcium indicator on the market, AAT Bioquest cements its position as *the* leader in calcium detection.

2018
AAT Bioquest adds a new linking technology to its toolkit. Sold under the **Buccutite™** brand, this technology makes bioconjugation of challenging proteins and enzymes as easy as "buckling your seatbelt."

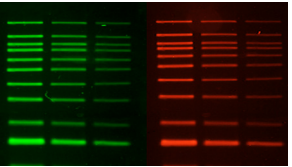


2019
Expanding into the multiplex immunofluorescence space, AAT Bioquest releases the **Styramide™** dyes, enabling ultra-sensitive detection of tissue targets. Simultaneously, **Transfectamine™** creates synergies in the lipofection and genomics space.

2020
AAT Bioquest mobilizes all resources to support the COVID-19 response. Viral protease substrates (**Covidyte™**) aid in the vaccine development effort.



2021
With an eye on safety and sustainability, AAT Bioquest commits to the development and release of **Gelite™ Safe**, a low-toxicity, high-sensitivity DNA gel electrophoresis dye.



2022
AAT Bioquest moves to a newly renovated, state-of-the-art facility in Pleasanton, California, greatly increasing its research capacity.

2024
Strengthening its core conjugation chemistry, AAT Bioquest releases **FastClick™**, greatly expanding the availability of click chemistry technology in its portfolio.

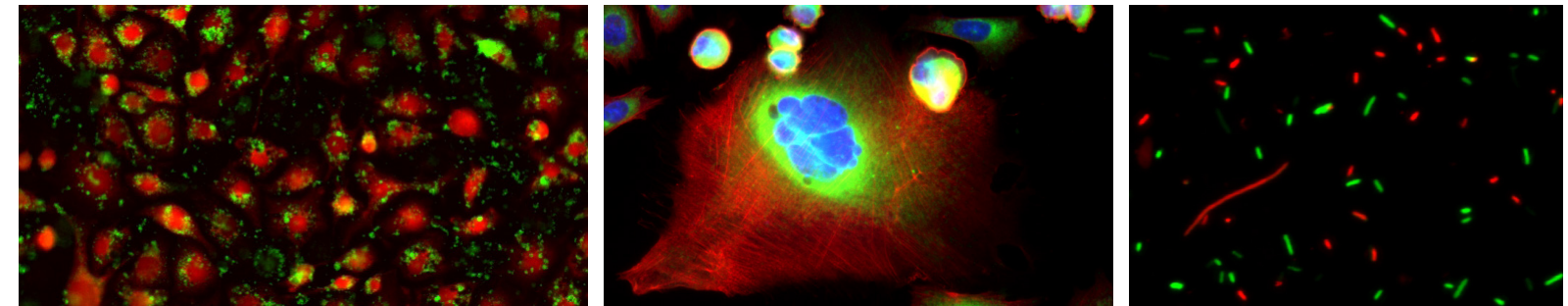
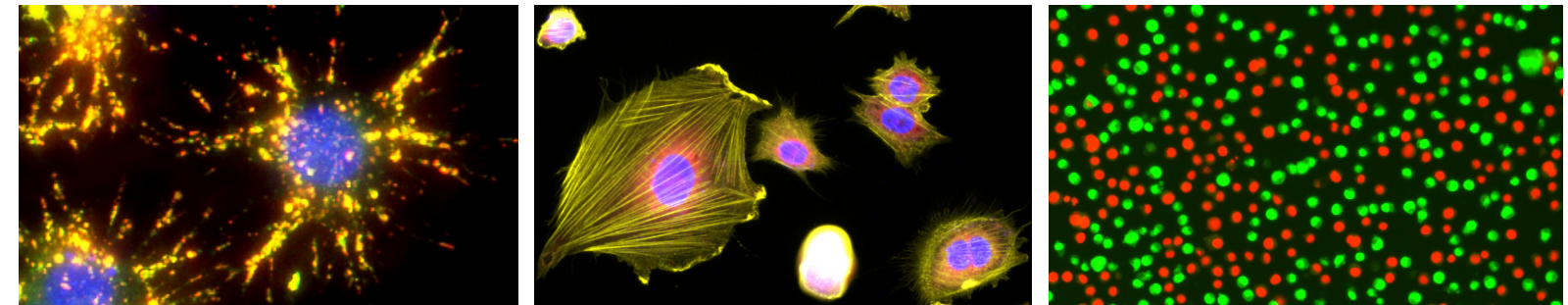
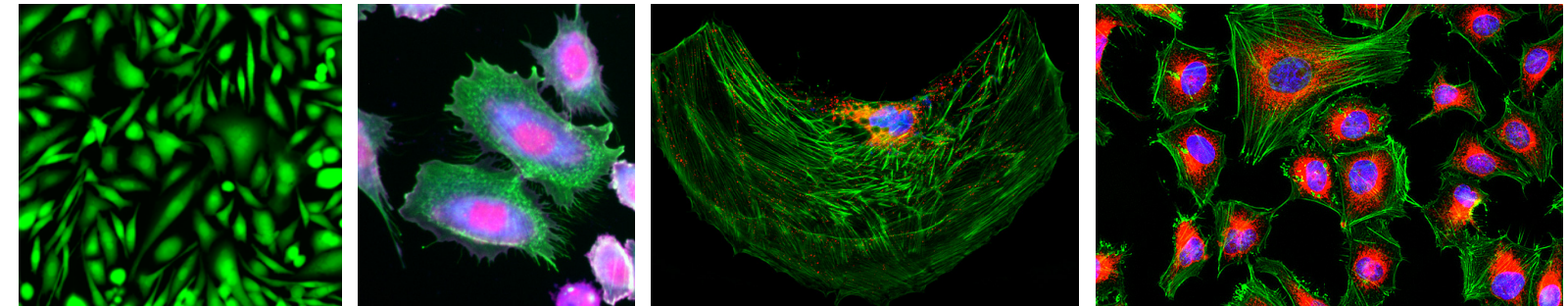
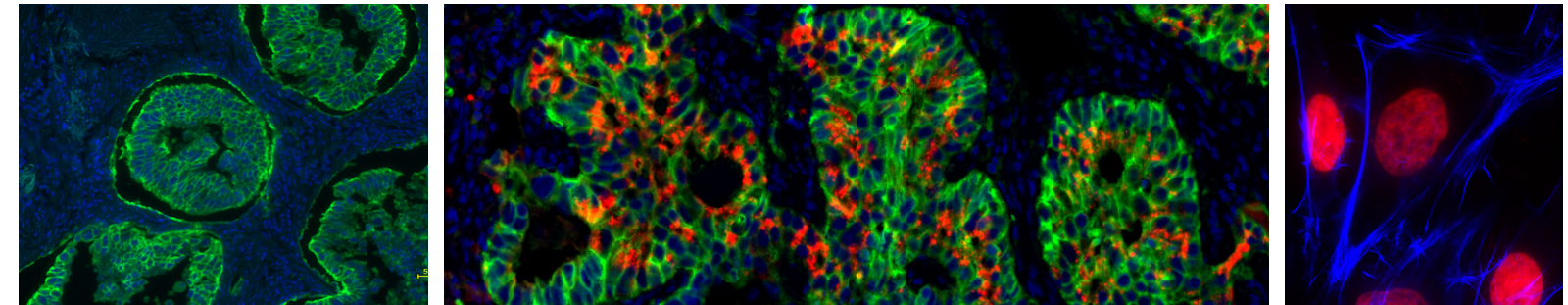
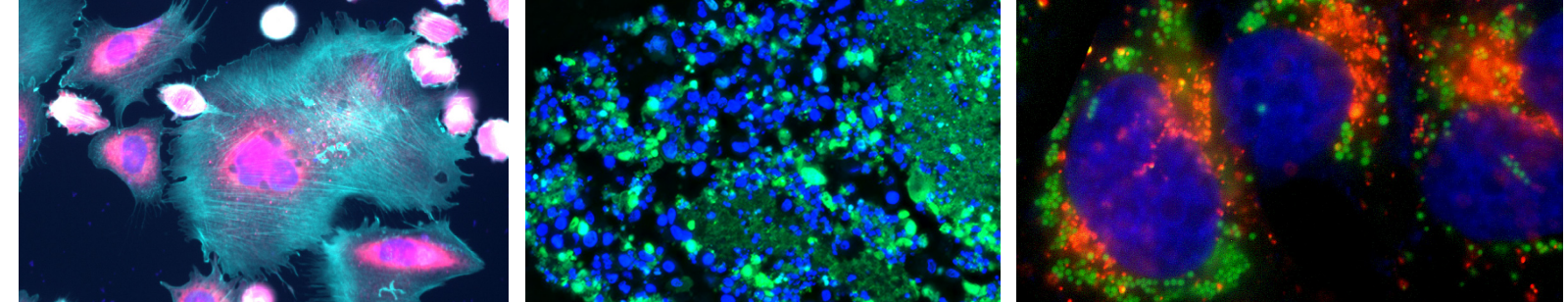
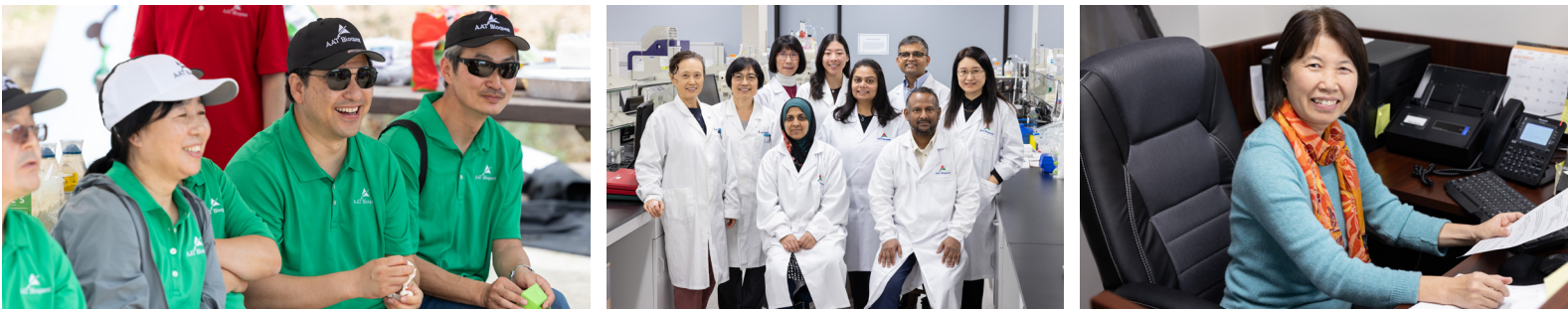


2026
AAT Bioquest expands upon the success of its **Buccutite™** technology by adding Fc-specific labeling capabilities, extending support to biologics development, in particular, antibody-drug conjugates (ADCs).

2025
With the release of **MitoDNA™** dyes and **fluorescently labeled beta-amyloid peptides**, AAT Bioquest looks to expand its disease-specific product portfolio, tackling challenging areas such as mitochondrial diseases and aging.

AAT Bioquest®

Together We Shine®





5775 W Las Positas Blvd.
Pleasanton, CA 94588 USA
Tel: 1-800-990-8053 • 1-408-733-1055
Fax: 1-800-609-2943 • 1-408-733-1304
sales@aatbio.com • www.aatbio.com

