

AACR

SAN DIEGO, CA

APRIL 17-22 2026

FIND US AT
#1819

Decoding the Tumor Microenvironment

Fluorescence reagents for multiplexed tissue profiling, immunophenotyping and oncology assay workflows

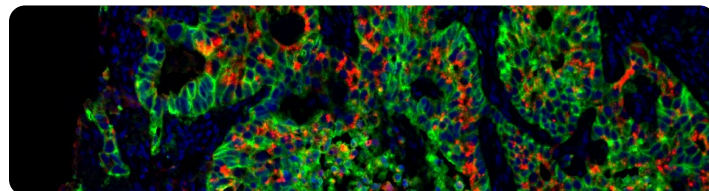
AAT Bioquest provides fluorescence reagents for multiplex immunofluorescence and spatial immuno-oncology workflows, including cyclic immunofluorescence and antibody conjugation. These tools support biomarker characterization and downstream therapeutic response studies, with applications in antibody-drug conjugate (ADC) research.

Map the Tumor Immune Landscape

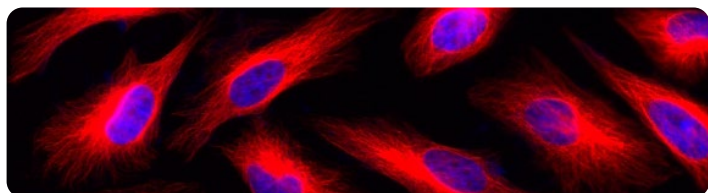
Power Styramide™ (PSA™) Platform

- Enhanced signal amplification for detection of low-abundance targets
- Compatible with FFPE tissue and cyclic IF workflows
- Enables detection of immune checkpoint markers such as PD-L1, LAG-3, and TIM-3

01



PSA™ iFluor® 488 multiplex IF



ReadiLink™ Rapid iFluor® 633 Antibody Labeling Kit

02

Profile Infiltrating Immune Populations

iFluor® Secondaries and Antibody Labeling Tools

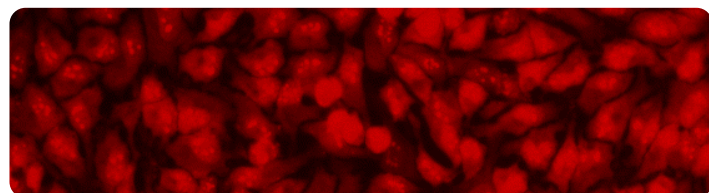
- Over 100 pre-conjugated iFluor® secondary antibodies spanning UV to NIR spectral range
- ReadiLink™ antibody labeling for flexible generation of custom conjugates
- Buccutite™ PE, APC, and tandem dye kits for high-parameter flow cytometry

Assess Cell Death and Stress Responses

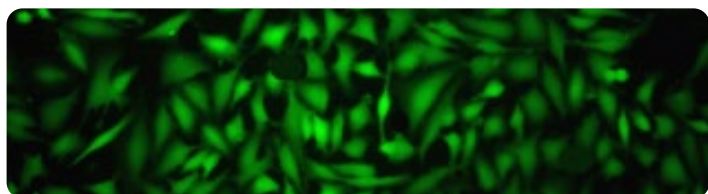
Apoptosis and Oxidative Stress Assays

- Multiplex caspase 3/7, 8, and 9 detections in a single assay for apoptosis signaling
- MitoROS™ and total ROS assays for measuring oxidative stress
- Annexin V-iFluor® conjugates with multiplex compatibility

03



Superoxide measurement using MitoROS™ 580



HeLa cells stained with Calcein UltraGreen™ AM

04

Evaluate Cell Viability and Proliferation

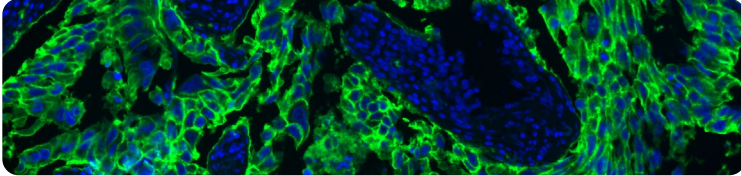
CCK-8, Calcein UltraGreen™ and CytoTell® Assays

- CCK-8 viability assay for quantitative measurement of metabolic activity
- Calcein UltraGreen™ AM for live cell staining with improved intracellular retention
- CytoTell® for monitoring cell division across multiple generations

The Immuno-Oncology Research Toolkit

SPATIAL BIOLOGY

Power Styramide™ (PSA™) Platform



iFluor® 488 Styramide™

PSA™ Reagents

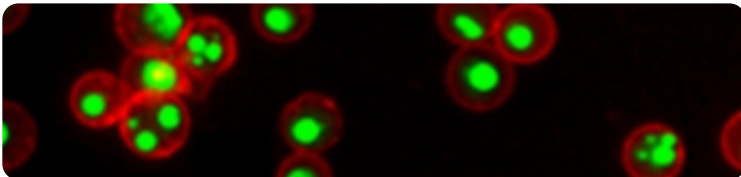
- Available in 19 iFluor® spectral variants (350–860 nm) for multiplex immunofluorescence
- mFluor™ and hapten-based Styramide™ reagents for flexible panel design

PSA™ Multiplex Imaging Kits

- Pre-optimized kits for multicolor staining in FFPE tissue sections
- Compatible with cyclic immunofluorescence workflows

CELL DEATH AND STRESS RESPONSE ASSAYS

Cell Meter™, MitoROS™, ApoSight™ and Annexin V-iFluor®



Jurkat cells stained with Annexin V-iFluor® 555

Apoptosis Assays

- Cell Meter™ Single or Multiplex Caspase Kits
- ApoSight™ Green and other caspase substrates
- Annexin V-iFluor® conjugates available in 10 colors for flow cytometry and imaging

Oxidative Stress Assays

- MitoROS™ reagents for mitochondrial superoxide detection
- Cell Meter™ ROS Assay Kit compatible with plate reader and flow cytometry

EXAMPLE

Multiparametric Functional Response Panel

Apoptosis → Annexin V-iFluor® 647

Mitochondrial superoxide → MitoROS™ 580

Cell viability → Calcein UltraGreen™ AM

IMMUNOFLUORESCENCE REAGENTS

Secondary Antibodies and Antibody Labeling Platforms

Pre-Conjugated Secondary Antibodies

- Over 100 pre-conjugated secondary antibodies- UV to NIR range (350–860 nm)
- Anti-mouse, anti-rabbit, and additional host-specific formats
- Available in iFluor®, Cy dye and more

Antibody Labeling and Conjugation Tools

- ReadiLink™ Kits for antibody labeling using two-step workflow without purification
- Buccutite™ Kits for PE, APC, and tandem dye conjugation (e.g., APC-Cy7, PE-Cy7) for high-parameter flow cytometry

Channel	Dye
FITC	iFluor® 488
PE/Texas Red	iFluor® 594
Cy5	iFluor® 647
NIR	iFluor® 750

CELL VIABILITY AND PROLIFERATION ASSAYS

CCK-8, Calcein UltraGreen™ and CytoTell®

Cell Viability

- CCK-8 (WST-8) assay for quantitative measurement of metabolic activity without cell lysis
- Calcein AM dyes for live-cell staining with improved intracellular retention
- Cell Explorer™ Live and Fixable Viability Kits compatible with intracellular staining workflows

Cell Proliferation

- CytoTell® series for tracking of cell division across multiple generations
- Bucculite™ FdU Cu-Free Cell Proliferation Kits with copper-free click chemistry for DNA synthesis detection

Calcein Variant	Ex/Em	Channel
Calcein UltraGreen™	491/516	FITC
Calcein Orange™	529/545	PE
Calcein Red™	563/588	Texas Red

EXAMPLE

Multiplex IO Panel (PSA™ Platform)

Checkpoint: PD-L1 PSA™ iFluor® 488

Immune infiltration: CD8 PSA™ iFluor® 555

Proliferation: Ki-67 PSA™ iFluor® 647

Tumor cells: PanCK PSA™ iFluor® 594

Nucleus: Nuclear Fragmentation Nuclear Blue™ DCS1

BOOTH
#1819

AAT BIOQUEST

AACR Annual Meeting 2026 | San Diego Convention Center
info@aatbio.com | 1-800-990-8053



aatbio.com