

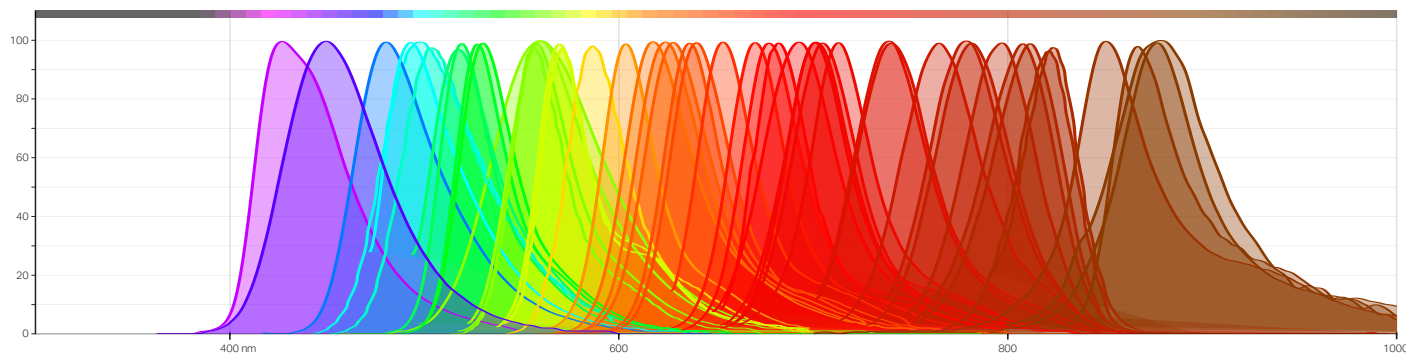
iFluor® Dyes

Brighter. More Stable. The Next Generation of Fluorescent Dyes.

AAT Bioquest's iFluor® dyes are a comprehensive family of fluorescent labeling dyes designed as improved alternatives to classic fluorophores like FITC, Alexa Fluor®, Cy®, Texas Red®, and IRDye®. Compatible with standard laser lines, the family was built to give fluorescence microscopy, flow cytometry, immunohistochemistry, Western blotting, and in vivo imaging workflows a single dye platform with broad reactivity coverage across the visible and infrared spectrum. Whether you are labeling your own antibodies, constructing a high-plex spectral flow panel, or running sensitive tissue imaging, iFluor® provides well-characterized options in every spectral window.

KEY FEATURES

- **45+ spectrally distinct colors** spanning 345–860 nm
- **Extinction coefficients up to 275,000 M⁻¹cm⁻¹** for high-sensitivity detection
- **Improved photostability** for extended imaging sessions
- **Water-soluble formulations** for aqueous labeling buffers
- **10+ reactive forms** — NHS esters, maleimides, azides, alkynes, tetrazines, BCN, etc.
- **Pre-validated conjugates** — antibodies, phalloidin, streptavidin, annexin V, WGA, CD markers
- **Spectral channels matched** for multicolor flow and imaging panels



Spectra image: Emissions of iFluor® dyes 350–860

SELECT PRODUCT CATEGORIES

IMMUNOFLUORESCENCE

iFluor® Secondary Antibodies

Pre-validated conjugates with optimized fluorophore-to-protein ratios for maximum brightness and minimal background in indirect immunofluorescence, Western blotting, and tissue imaging.

20+ Colors Available

CYTOSKELETAL IMAGING

Phalloidin-iFluor®

High-affinity F-actin probes for visualizing actin filaments in fixed cells and tissue sections. Effective at nanomolar concentrations for high-contrast cytoskeletal imaging with minimal non-specific binding.

10+ Colors Available

RAPID ANTIBODY LABELING

ReadiLink™ Kits

Label your own antibodies or proteins with iFluor® dyes in under 60 minutes—no purification required. Quenching technology eliminates free dye background without column cleanup.

30+ Colors Available

iFluor® Dyes – Technical Specifications & Ordering Guide

UV: 355 nm Violet: 405 nm Blue: 488 nm Green: 532 nm Yellow: 561 nm Orange: 594 nm Red: 640 nm Far-Red: 730 nm NIR: 785 nm							
iFluor® Dye	Ex (nm)	Em (nm)	QY	EC (M ⁻¹ cm ⁻¹)	Laser	Channel	Replaces
iFluor® 350	345	450	0.95	20,000	355 nm	DAPI / Hoechst	AMCA, Alexa Fluor® 350
iFluor® 405	403	427	0.91	37,000	405 nm	BV421 / Pacific Blue	No direct equivalent
iFluor® 430	433	497	0.78	20,000	405 nm	BV510 / AmCyan	Alexa Fluor® 430
iFluor® 440	434	480	0.67	40,000	405 nm	BV510 / AmCyan	DEAC, SpectrumAqua
iFluor® 445	446	558	N/D	N/D	405 nm	FITC / GFP	No direct equivalent
iFluor® 450	451	502	0.82	40,000	405 nm	BV510 / AmCyan	Alexa Fluor® 450
iFluor® 488	491	516	0.90	75,000	488 nm	FITC / GFP	FITC, Alexa Fluor® 488, Cy2
iFluor® 500	501	520	N/D	80,000	488 nm	FITC / GFP	No direct equivalent
iFluor® 510	511	530	N/D	N/D	488 nm	FITC / GFP	No direct equivalent
iFluor® 514	511	527	0.83	75,000	488 nm	FITC / GFP	Alexa Fluor® 514
iFluor® 532	537	560	0.68	90,000	532 nm	Cy3 / TRITC	Alexa Fluor® 532
iFluor® 540	540	557	N/D	N/D	532 nm	Cy3 / TRITC	No direct equivalent
iFluor® 546	541	557	0.67	100,000	532 nm	Cy3 / TRITC	Alexa Fluor® 546
iFluor® 555	557	570	0.64	100,000	532 nm	PE	Cy3, TRITC, Alexa Fluor® 555
iFluor® 560	560	571	0.57	120,000	561 nm	PE	No direct equivalent
iFluor® 568	568	587	0.57	100,000	561 nm	PE	Alexa Fluor® 568, Rhodamine Red
iFluor® 570	557	570	0.71	120,000	561 nm	PE	No direct equivalent
iFluor® 594	587	603	0.53	200,000	594 nm	Texas Red	Texas Red, Alexa Fluor® 594
iFluor® 597	598	618	0.70	100,000	594 nm	Texas Red	No direct equivalent
iFluor® 605	603	623	N/D	N/D	594 nm	Texas Red	No direct equivalent
iFluor® 610	610	628	0.85	110,000	594 nm	Texas Red	No direct equivalent
iFluor® 620	621	636	N/D	N/D	640 nm	Texas Red	No direct equivalent
iFluor® 625	624	640	N/D	N/D	640 nm	Texas Red	No direct equivalent
iFluor® 633	640	654	0.29	250,000	640 nm	APC / Cy5	Alexa Fluor® 633
iFluor® 647	656	670	0.25	250,000	640 nm	APC / Cy5	Cy5, Alexa Fluor® 647
iFluor® 660	663	678	0.26	250,000	640 nm	APC / Cy5	Alexa Fluor® 660
iFluor® 665	667	692	0.22	110,000	640 nm	Cy5.5	Alexa Fluor® 660
iFluor® 670	671	682	0.55	200,000	640 nm	Cy5.5	Alexa Fluor® 670, Cy5B
iFluor® 675	683	700	N/D	N/D	640 nm	Cy5.5	No direct equivalent
iFluor® 680	684	701	0.23	220,000	640 nm	Cy5.5	Alexa Fluor® 680, IRDye® 680, Cy5.5
iFluor® 690	685	704	0.30	220,000	640 nm	Cy5.5	No direct equivalent
iFluor® 700	690	713	0.23	220,000	640 nm	NIR (700+)	Alexa Fluor® 700
iFluor® 720	716	740	0.14	240,000	730 nm	NIR (700+)	No direct equivalent
iFluor® 740	740	764	0.20	225,000	730 nm	NIR (700+)	No direct equivalent
iFluor® 750	757	779	0.12	275,000	730 nm	NIR (700+)	Cy7, Alexa Fluor® 750, HiLyte™ 750
iFluor® A7	764	782	0.10	275,000	785 nm	NIR (700+)	APC-H7 (as tandem acceptor)
iFluor® 770	777	797	0.16	250,000	785 nm	NIR (700+)	IRDye® 800 RS
iFluor® 780	784	808	0.16	250,000	785 nm	NIR (700+)	IRDye® 800 CW
iFluor® 790	787	812	0.13	250,000	785 nm	NIR (700+)	IRDye® 800 CW, Alexa Fluor® 790, ICG
iFluor® 800	801	820	0.11	250,000	785 nm	NIR (700+)	ICG, IRDye® 800 CW
iFluor® 810	811	822	0.05	250,000	785 nm	NIR (700+)	No direct equivalent
iFluor® 820	822	850	N/D	250,000	785 nm	IR (830+)	No direct equivalent
iFluor® 830	830	867	N/D	N/D	785 nm	IR (830+)	No direct equivalent
iFluor® 840	836	879	N/D	200,000	785 nm	IR (830+)	No direct equivalent
iFluor® 860	853	878	N/D	250,000	785 nm	IR (830+)	No direct equivalent