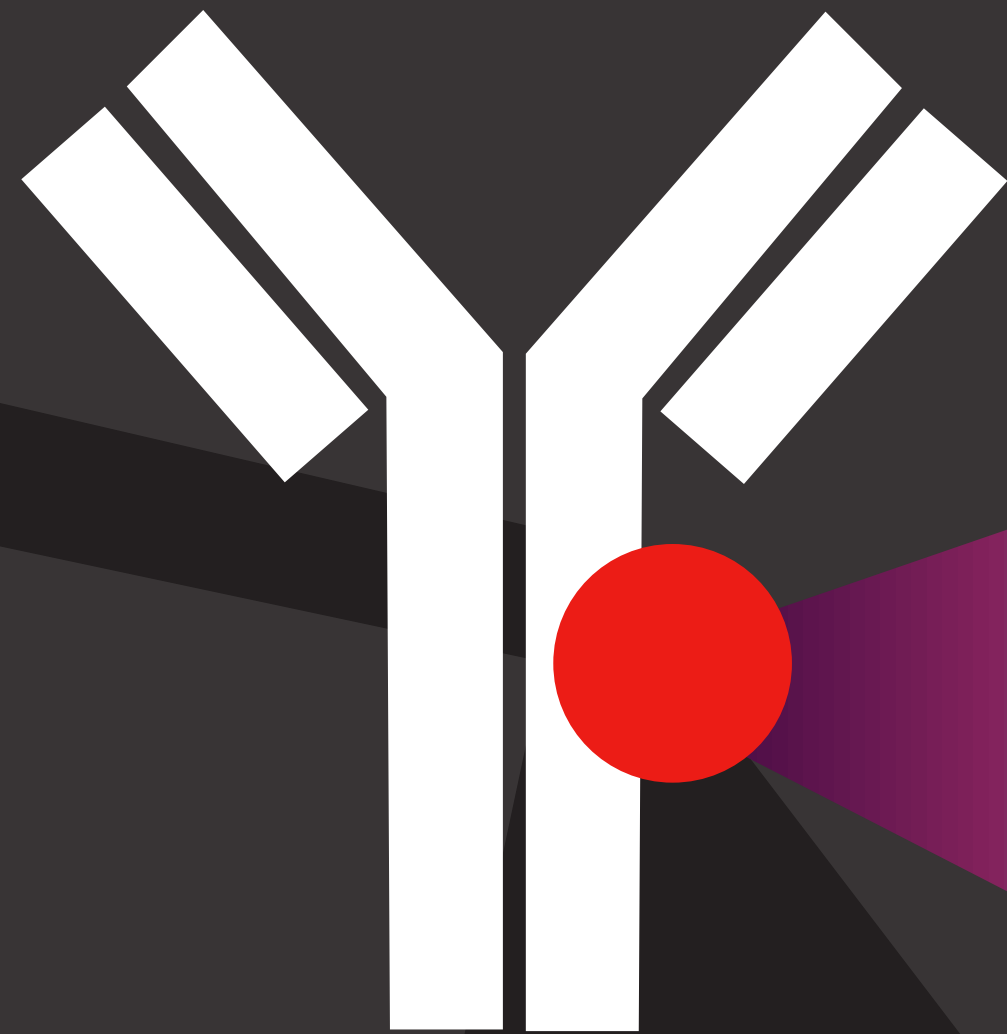


# FLUORESCENT LABELS

Browse pre-labeled antibodies and Lightning-Link™ Antibody Labeling Kits at [www.novusbio.com](http://www.novusbio.com)



**AMCA**  
Abs: 346 nm  
Emi: 442 nm

**Atto390**  
Abs: 390 nm  
Emi: 479 nm

**Atto425®**  
Abs: 436 nm  
Emi: 484 nm

**Cyanine Dye 2 (Cy2)**  
Abs: 489 nm  
Emi: 506 nm

**DyLight™ 488**  
Abs: 493 nm  
Emi: 518 nm

**Atto488**  
Abs: 501 nm  
Emi: 523 nm

**FITC**  
Abs: 494 nm  
Emi: 518 nm

**Atto465**  
Abs: 453 nm  
Emi: 508 nm

**Alexa Fluor® 488**  
Abs: 495 nm  
Emi: 519 nm

**Atto532**  
Abs: 532 nm  
Emi: 553 nm

**DyLight™ 549**  
Abs: 562 nm  
Emi: 576 nm

**R-Phycoerythrin (R-PE)**  
Abs: 490; 565 nm  
Emi: 578 nm

**Rhodamine**  
Abs: 552 nm  
Emi: 575 nm

**Cyanine Dye 3 (Cy3)**  
Abs: 550 nm  
Emi: 570 nm

**TRITC**  
Abs: 544 nm  
Emi: 572 nm

**FluoProbes®547H**  
Abs: 557 nm  
Emi: 572 nm

**Cyanine Dye 3.5 (Cy3.5)**  
Abs: 581 nm  
Emi: 596 nm

**Texas Red**  
Abs: 596 nm  
Emi: 615 nm

**Atto590**  
Abs: 600 nm  
Emi: 627 nm

**Atto594**  
Abs: 601 nm  
Emi: 627 nm

**Atto565**  
Abs: 563 nm  
Emi: 592 nm

**Atto610**  
Abs: 575 nm  
Emi: 640 nm

**Atto620**  
Abs: 619 nm  
Emi: 643 nm

**Atto633**  
Abs: 629 nm  
Emi: 657 nm

**Atto637**  
Abs: 630 nm  
Emi: 659 nm

**Atto655**  
Abs: 663 nm  
Emi: 684 nm

**DyLight™ 649**  
Abs: 654 nm  
Emi: 673 nm

**Cyanine Dye 5 (Cy5)**  
Abs: 643 nm  
Emi: 667 nm

**Allophycocyanin (APC)**  
Abs: 650 nm  
Emi: 661 nm

**FluoProbes®647H**  
Abs: 653 nm  
Emi: 675 nm

**PerCP**  
Abs: 490 nm  
Emi: 675 nm

**Atto611**  
Abs: 611 nm  
Emi: 681 nm

**Cyanine Dye 5.5 (Cy5.5)**  
Abs: 675 nm  
Emi: 694 nm

**FluoProbes®682**  
Abs: 690 nm  
Emi: 709 nm

**FluoProbes®752**  
Abs: 748 nm  
Emi: 772 nm

**Cyanine Dye 7 (Cy7)**  
Abs: 747 nm  
Emi: 776 nm

**Atto680**  
Abs: 670nm  
Emi: 710 nm

**Atto700**  
Abs: 700nm  
Emi: 725 nm



Fluorescent dyes are environmentally sensitive and may produce varying peak excitation and emission wavelengths due to buffer type, instrument settings, etc. Please refer to product datasheets and your instrument handbooks for guidance prior to designing and conducting your experiments. Human vision is in the range from 380-790 nm, thus special equipment is required to visualize the far-red fluorescent dyes. Alexa Fluor® is a Registered Trademark of Molecular Probes Inc, USA. ATTO Dyes are manufactured by ATTO-TEC GmbH. CyDye® is a Registered Trademark of GE Healthcare. DyLight™ is a Trademark of Thermo Fisher Scientific and its subsidiaries. Fluorescent Dyes Poster © Novus Biologicals, LLC 2010. All Rights Reserved.

[www.novusbio.com](http://www.novusbio.com) • [novus@novusbio.com](mailto:novus@novusbio.com)