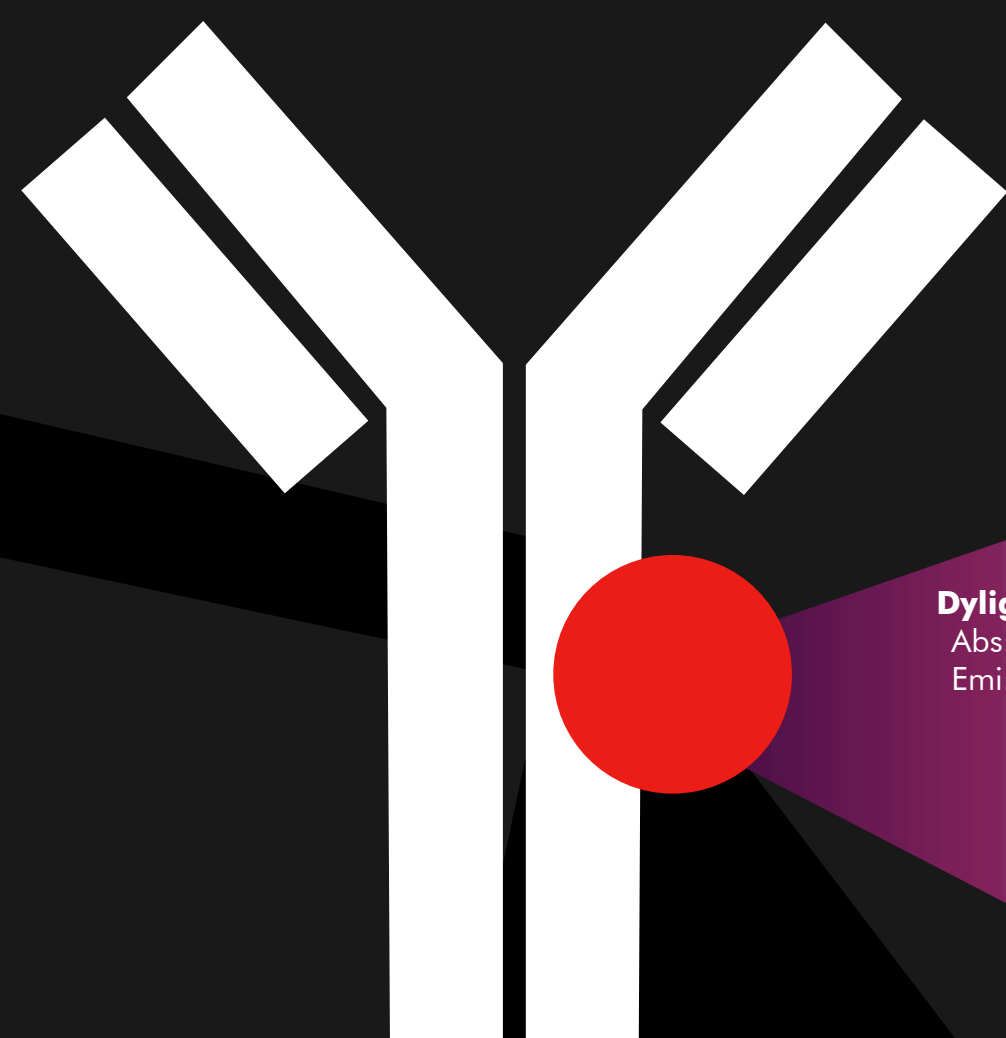


FLUORESCENT LABELS

Y Browse pre-labeled antibodies and Lightning-Link™ Antibody Labeling Kits at www.novusbio.com/Fast-Antibody-Labeling.html Y



Dylight™ 405
Abs: 400 nm
Emi: 420 nm

Dylight™ 350
Abs: 353 nm
Emi: 432

AMCA 405
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

Atto465
Abs: 453 nm
Emi: 508 nm

FITC
Abs: 494 nm
Emi: 518 nm

Fluorescein
Abs: 494 nm
Emi: 521 nm

Atto488
Abs: 501 nm
Emi: 523 nm

HiLyte Fluor™ 488
Abs: 497 nm
Emi: 525 nm

Atto532
Abs: 533 nm
Emi: 553 nm

HiLyte Fluor™ 555
Abs: 550 nm
Emi: 566 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

Dyomics™ 547
Abs: 557 nm
Emi: 574 nm

DyLight™ 550
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 3.5 (Cy3.5™)
Abs: 581 nm
Emi: 596 nm

Atto 565
Abs: 563 nm
Emi: 592 nm

DyLight™ 594
Abs: 583 nm
Emi: 618 nm

Atto594
Abs: 601 nm
Emi: 627 nm

Atto 590
Abs: 600 nm
Emi: 627 nm

Texas Red®
Abs: 596 nm
Emi: 615 nm

DyLight™ 633
Abs: 638 nm
Emi: 658 nm

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

Atto637
Abs: 630 nm
Emi: 659 nm

Dyomics™ 647
Abs: 653 nm
Emi: 672 nm

HiLyte Fluor™ 647
Abs: 649 nm
Emi: 672 nm

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

DyLight™ 650
Abs: 654 nm
Emi: 673 nm

Atto680
Abs: 670 nm
Emi: 710 nm

DyLight™ 680
Abs: 692 nm
Emi: 712 nm

PerCP
Abs: 490 nm
Emi: 675 nm

Atto700
Abs: 700 nm
Emi: 725 nm

FluoProbes™ 674H
Abs: 653 nm
Emi: 675 nm

FluoProbes® 682
Abs: 690 nm
Emi: 709 nm

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

Atto655
Abs: 663 nm
Emi: 684 nm

FluoProbes® 752
Abs: 748 nm
Emi: 772 nm

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

DyLight™ 755
Abs: 754 nm
Emi: 776 nm

HiLyte Fluor™ 750
Abs: 750 nm
Emi: 782 nm

DyLight™ 800
Abs: 777 nm
Emi: 794 nm