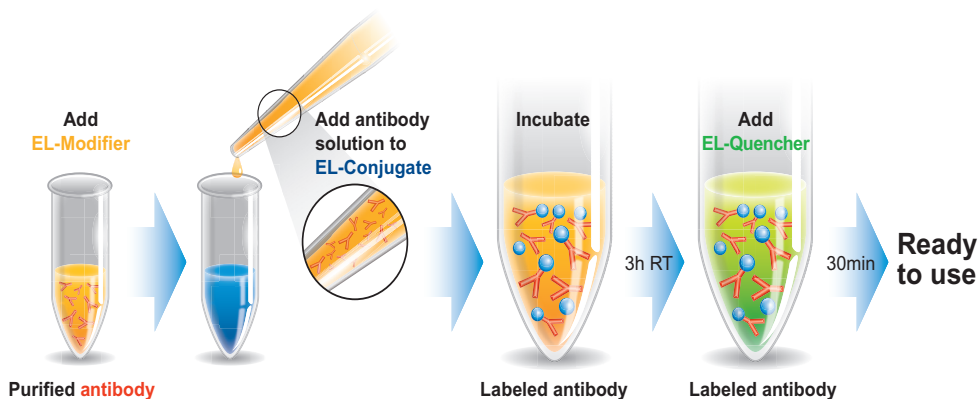


# EasyLink antibody conjugation kits

The easiest way to conjugate your antibodies



The EasyLink process involves free amine groups

- Suitable for **micrograms** to **milligrams** of antibodies
- In the presence of up to 0.5% BSA, 20mM Tris and 0.1% sodium azide
- One-step antibody labeling method
- No separation steps

Concentration and purification kits are also available at:  
[www.abcam.com/EasyLink](http://www.abcam.com/EasyLink)


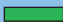
















## Conjugates for enzymatic detection

Conjugation	3x10µg	1x100µg	3x100µg	1x1mg
AP	ab102851	ab102849	ab102850	ab102848
Glucose oxidase (GOx)	ab102888	-	ab102887	ab102886
HRP	ab102892	ab102890	ab102891	ab102889

## Conjugates for biotin/streptavidin detection

Conjugation	3x10µg	1x100µg	3x100µg	1x1mg
Avidin	ab102862	-	ab102861	ab102860
Biotin (type A)	ab102865	-	ab102864	ab102863
Biotin (type B)	ab102868	-	ab102867	ab102866
Streptavidin	ab102922	-	ab102921	ab102920

## Conjugates for fluorescent detection

Conjugation	Colour	3x10µg	1x100µg	3x100µg	1x1mg
AMCA		ab102847	-	ab102846	ab102845
FITC		ab102885	-	ab102884	ab102883
Cy3 <sup>®</sup>		ab102874	-	ab102873	ab102872
Rhodamine		ab102915	-	ab102914	ab102913
Cy3.5 <sup>®</sup>		-	-	ab104332	ab104333
B-PE		-	ab102870	ab102871	ab102869
R-PE		ab102919	ab102917	ab102918	ab102916
Texas Red <sup>®</sup>		ab102925	-	ab102924	ab102923
Cy5 <sup>®</sup>		ab102877	-	ab102876	ab102875
Pe/Cy5 <sup>®</sup>		ab102896	ab102894	ab102895	ab102893
PerCP		ab102908	ab102906	ab102907	ab102905
PerCP/Cy5.5 <sup>®</sup>		ab102912	ab102910	ab102911	ab102909
PE/Cy5.5 <sup>®</sup>		ab102900	ab102898	ab102899	ab102897
APC/Cy5.5 <sup>®</sup>		ab102855	ab102853	ab102854	ab102852
Cy5.5 <sup>®</sup>		-	-	ab102879	ab102878
APC/Cy7 <sup>®</sup>		ab102859	ab102857	ab102858	ab102856
PE/Cy7 <sup>®</sup>		ab102904	ab102902	ab102903	ab102901
Cy7 <sup>®</sup>		ab102882	-	ab102881	ab102880

Learn more about our EasyLink conjugation kits at:

[www.abcam.com/EasyLink](http://www.abcam.com/EasyLink)

