



# R-PE Antibody Conjugation Kit

## P-9002-002

Product Images



## Short Description

---

The R-PE Antibody Conjugation Kit is a flexible alternative to the R-PE Antibody All-in-One™ Conjugation Kit. Each kit contains sufficient material for two reactions of up to 1.3 mg antibody each. Based on SoluLINK® bioconjugation technology, the reactions are rapid and high-yielding, converting nearly 100% of antibody to PE conjugate. This kit is ideal for applications where purification to remove excess R-PE is either not required, or purification will be carried out by chromatographic methods (such as size exclusion FPLC).

## Additional Information

---

Conjugation Targets	Antibody
Label/Modifier Type	R-PE
Reactivity	Amine
Recommended Storage	2° - 8° C — Do Not Freeze
Applications	Antibody Labeling

## Products in this set

---



**R-PE Antibody Conjugation Kit, 1.00 Kit**

