



R-PE Antibody Conjugation Kit

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

