



R-PE Antibody All-in-One Conjugation Kit

Product Images

Additional Information

Conjugation Targets	Antibody
Label/Modifier Type	R-PE
Reactivity	Amine
Components	Collection Tubes, S-HyNic Crosslinker (DMF Soluble), Modification Buffer (10X), Anhydrous DMF
Recommended Storage	4°C
Applications	Antibody Labeling, Antisense/RNAi, Aptamers, Immunotherapeutics, In Situ Proximity Ligation, In vitro Diagnostic (IVD), Laboratory Developed Tests (LDT), Analyte Specific Reagents (ASR)
Unit Size	1 Kit

Products in this set



R-PE Antibody All-in-One Conjugation Kit, 1.00 Kit

