



APC-Antibody Conjugation Kit

P-9903

Product Images

Additional Information

Conjugation Targets	Antibody
Label/Modifier Type	APC
Reactivity	Amine
Components	S-HyNic Crosslinker (DMF Soluble), Modification Buffer (10X), Anhydrous DMF, Zeba Desalting Columns, 1.5 mL Collection Tubes, 10X MES Buffer, 2-Sulfobenzaldehyde, Conjugation Buffer (10X), 4FB-Modified APC
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)

Products in this set



APC-Antibody Conjugation Kit, 1.00 Kit

