Methods for Detecting m6A in RNA

MAPPING

M6A

m6A

RNA

LOWER COST

MeRIP Steps

Most common method for mapping

RNA extraction **RNA** fragmentation Bead-coupled Ab binding Immunoprecipitation of m6A Separation and elution **Reverse transcription** Library construction Sequencing

meRIP

ELISA Steps

Quantifies m6A levels in RNA

RNA extraction RNA binding to assay wells m6A antibody capture **Detection antibody incubation** Addition of enzyme substrate Signal measurement Standard curve calculation

LESS TIME

meRIP

Mapping m6A RNA methylation sites transcriptome-wide methylation sites using imunoprecipitation

miCLIP

ELISA

High-resolution mapping of m6A in RNA at Individual-nucleotide resolution using crosslinking and immunoprecipitation

HIGHER COST

miCLIP Steps

RNA extraction **RNA** fragmentation Incubation with anti-m6A Ab UV-crosslinking Protein A/G affinity purification Membrane transfer Adaptor ligation Proteinase K treatment **Reverse transcription** Library construction Sequencing

HPLC/MS Steps

mRNA isolation Nuclease P1 and alkaline phosphatase digestion **Digested RNA purification** Separation by reverse phase HPLC Mass spectrometry analysis Standard curve calculation

HPLC/MS

miCLIP

M6A MEASURING

ELISA

Enzyme-linked immunosorbent plate-based assay that measures m6A levels in RNA

MORE TIME



analysis



