

When biotinylating a molecule, is there a minimum linker length?

Longer linkers are typically better. In our experience, 29 angstroms or longer should allow the attached molecule to be presented without masking the active site. For details on the biotinylation protocol, refer to Technical Note 28: Biotinylation of Protein for Immobilization onto Streptavidin Biosensors.

Article ID: 103

Posted: Fri, Nov 11, 2016 at 3:07 PM.

Last Updated: Fri, Nov 11, 2016 at 3:07 PM

Filed Under: Assays
