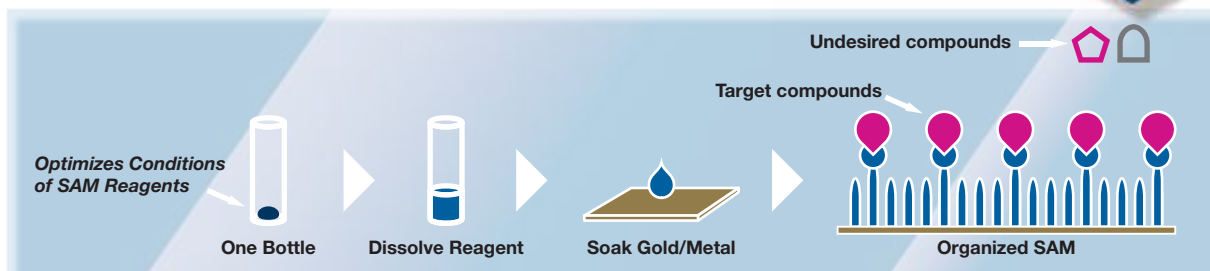


SAM FORMATION REAGENT

SAM (Self-Assembled Monolayer) is a highly organized thin layer of molecules formed on a solid surface by chemisorption. Due to its high applicability, SAM can be used on biosensors. However, unspecific binding is commonly observed while biomaterial samples are used. In the most of these cases, the concentration and combination of SAM reagents are not optimized for preparing an organized layer on a metal surface. Dojindo provides SAM Formation Reagents which are already optimized in one bottle and simple to use. The optimized reagent can reduce unspecific binding significantly.



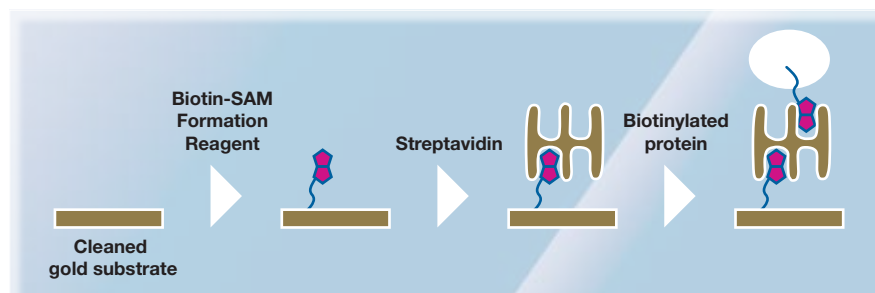
BIOTIN-SAM FORMATION REAGENT

Code: B564-10

Unit: 2 μmol x 3

Avidin-Biotin method has been widely used in the field of enzyme immunoassay (EIA). This method allows the rapid and strong immobilization of proteins such as Biotinylated-antibodies, -peptides, and -DNA on a biosensor.

For more details, please visit our web-site at www.dojindo.eu.com



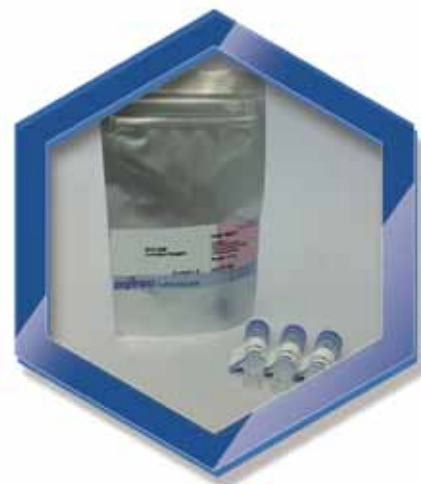
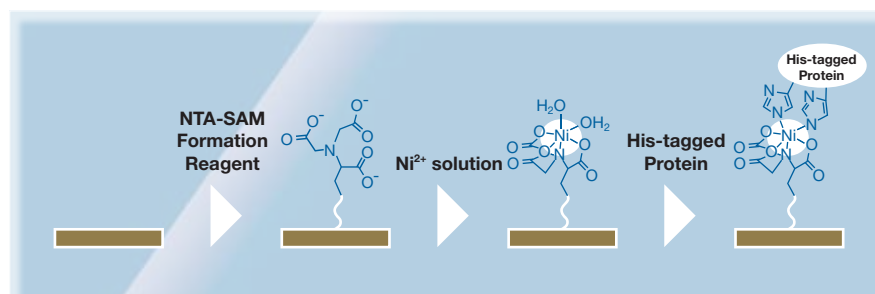
NTA-SAM FORMATION REAGENT

Code: N475-10

Unit: 2 μmol x 3

NTA forms a stable complex with heavy metals such as Ni, it is, therefore, applied to introduce metal ions onto NTA-biomolecules or NTA-compounds. These metal chelate compounds have been commonly used to detect specific substances which interact with metal ions such as His-Tagged proteins.

For more details, please visit our web-site at www.dojindo.eu.com



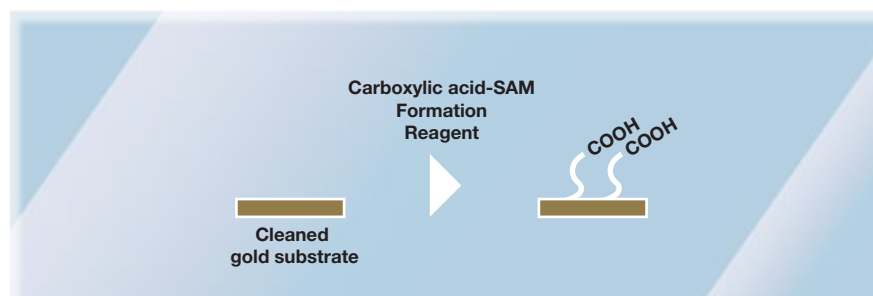
CARBOXYLIC ACID-SAM FORMATION REAGENT

Code: C488-10

Unit: 1 μmol x 3

The carboxylic acid-SAM immobilizes proteins or peptides and other ligands on the metal surfaces by activating a carboxylic acid group. To activate a carboxylic acid group, the Amine Coupling Kit [A515-10] is applicable.

For more details, please visit our web-site at www.dojindo.eu.com



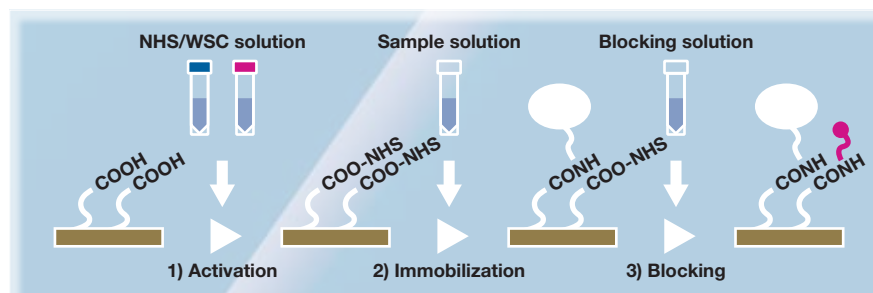
AMINE COUPLING KIT

Code: A515-10

Unit: 2 ml x 4

Amine and carboxylic acid coupling is one of the most common methods to immobilize a protein or a peptide through a covalent bond onto a biosensor surface. Dojindo's Amine Coupling Kit contains all the reagents and buffers necessary for activation of carboxylic acid, protein immobilization, and blocking. A blocking reagent minimizes non-specific protein absorption of the residual activated esters by capping them. Each kit is adequate for approximately 40 immobilizations.

For more details, please visit our web-site at www.dojindo.eu.com



DOJINDO STATE OF THE ART
LIFE SCIENCE TECHNOLOGIES

EUROPEAN HEADQUARTERS
DOJINDO EU GMBH

Leopoldstr. 254, 80807 Munich, Germany
Phone +49 89 3540-4805
Fax +49 89 3540-4806
e-mail info@dojindo.eu.com

www.dojindo.eu.com

DISTRIBUTED BY