

Carboxylic acid-SAM Formation Reagent Technical Manual

Technical Manual (Japanese version) is available at <http://www.dojindo.co.jp/manual/c488.pdf>

General Information

Carboxylic acid-SAM Formation Reagent is applied for the preparation of carboxylic acid terminated self-assembled monolayers (SAMs) on a gold surface of biosensors such as for QCM (Quartz Crystal Microbalance), SPR (Surface Plasmon Resonance) or electrical analysis. Carboxylic acid-SAM can be used as an interface to immobilize a protein or a peptide on a gold surface coupled with carboxyl group activation method using NHS/WSC. Amine Coupling Kit (Product code: A515) is available for protein immobilization on the gold surface coated with a carboxylic acid-SAM.

Carboxylic acid-SAM Formation Reagent is especially designed to minimize a non-specific protein adsorption on the surface.

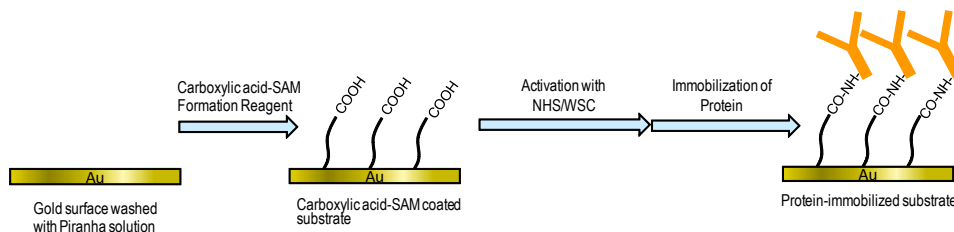


Fig.1 Preparation of a biosensor using Carboxylic acid-SAM Formation Reagent

Contents

1 $\mu\text{mol} \times 3$

Caution

Since a trace amount of colorless or slightly yellowish liquid is in the tube, please centrifuge prior to use.

Storage

0-5 °C

General Protocol for Carboxylic acid-SAM Preparation

Preparation of a Carboxylic acid-SAM surface on a gold substrate

- 1) Add 1 ml ethanol to a tube, and pipette to prepare 1 mmol/l Carboxylic acid-SAM solution. ^{a)} Then, dilute the solution 10-fold with ethanol for Step 2).
- 2) Immerse a gold substrate ^{b)} in the reagent solution prepared at Step 1) at room temperature and leave it overnight.
- 3) Wash the substrate several times with ethanol and purified water sequentially. ^{c)}

a) If the reagent hardly dissolve by pipetting, use an ultrasonic bath. Use freshly prepared reagent solution at Step 2).

b) Clean the gold substrate with Piranha solution, if necessary. Mix 3 volumes of sulfuric acid and one volume of hydrogen peroxide solution to prepare Piranha solution. Since the solution is severely caustic, handle with great care.

c) Store the SAM-coated substrate under nitrogen gas in a tightly sealable glass container at 0-5 °C.

General Protocol for Protein Immobilization

This product does not include necessary reagents and solutions for a protein immobilization. Amine Coupling Kit (Product code: A515) is available for the protein immobilization on the carboxylic acid-SAM surface. If not use the kit, follow this general protocol.

- Required reagents and buffers-

- N-Hydroxysuccinimide (NHS)
- 1-Ethyl-3-(3-dimethylaminopropyl)carbodiimide (WSC)
- pH4-6 buffer solution
 - * MES buffer is generally used.
- pH8-9 buffer solution
 - * Carbonate buffers, phosphate buffers or good's buffers are generally used.

- 1) Prepare 100 mmol/l NHS and 100 mmol/l WSC solutions using pH4-6 buffer solution .
- 2) Mix equal volume of NHS and WSC solutions immediately, and add to the carboxylic acid-SAM coated gold substrate.
- 3) Leave it at room temperature for 1 hr, and wash several times with pH4-6 buffer solution.
- 4) Add a protein solution ^{a)} to the substrate, and leave it at room temperature for 1 hr.
- 5) Wash with PBS, and go to blocking process with a blocking solution ^{b)} if necessary.

a) The solution should be prepared with a buffer solution (pH8-9). Use 10-1,000 $\mu\text{g}/\text{ml}$ of protein solution for immobilization.

b) One mol/l ethanolamine solution (pH8.5) is commonly used for blocking.

Related Products

The following related products are available.

Product Name	Product Code
Amine Coupling Kit	A515
WSC	W001

* Please visit our web site (www.dojindo.co.jp) for the updated prices.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories
2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.
Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>
Dojindo EU GmbH
Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>
Dojindo China Co., Ltd
Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>