

June, 2022

Keywords:

UniSart® Membrane, Lateral Flow Assay, Surfactants

Surfactants Selection Guide for Lateral Flow Assays

Dr. Shuo Liang

Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079
Goettingen

Correspondence

E-Mail: shuo.liang@sartorius.com

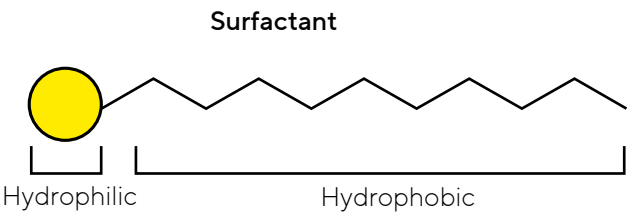


Introduction

Lateral flow assays (LFA) are a type of immunographic assay widely used for the rapid detection of pathogens, toxins, environmental contaminants, etc. Their ease of use and accuracy make these assays well suited for the point-of-care market. For purposes of this study, three lateral flow tests were selected. The effects of surfactants on test performance were demonstrated by adding different surfactants (surface active agent) during the pre-treatment of sample pads.

1. Classification of surfactants and their function in lateral flow assay:

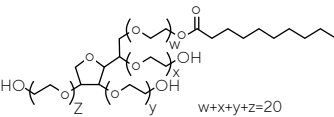
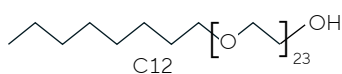
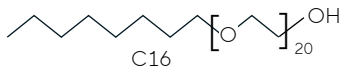
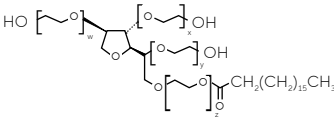
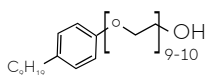
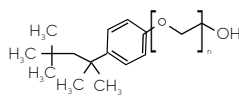
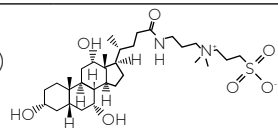
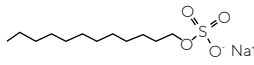
Figure 1: Different forms of surfactants (Surface Active Agent)



Classification		Function in Lateral Flow Assay
Non-ionic		▪ Enhance or decrease the test line intensity
Cationic		▪ Accelerate or decelerate the background clearance
Anionic		▪ Reduce the unspecific binding
Zwitterionic		▪ Increase wettability of the pad material ▪ Support the blocking

As listed in Table 1 the following surfactants were selected. All surfactants were used as additives in pre-treatment of samples pads of the lateral flow tests.

Table 1. Overview of surfactants used for this study

Commercial Name	Structure	Type	GHS	HLB	Application
Tween 20 (Polysorbate 20)		Non-ionic		16.7	Gentle surfactant and commonly used in LFA
Brij 35		Non-ionic		16.9	Performance optimization ¹
Brij 58		Non-ionic		15.7	Performance optimization
Tween 60 (Polysorbate 60)		Non-ionic		14.9	Gentle surfactant
Tergitol™ type NP-40		Non-ionic	  	17.8	Background clearance and conjugate pad release
IGEPAL®CA-630		Non-ionic	  	13.4	Background clearance and conjugate pad release
CHAPS (3-[(3-cholamidopropyl) dimethylammonio]-1- propanesulfonate)		Zwitterionic		/	Protein solubilization
SDS (sodium dodecyl sulfate)		Anionic	  	40.0	Commonly used in protein chemistry

HLB: hydrophilic-lipophilic balance, high value hydrophilic (water soluble), low value hydrophobic (oil soluble)

2. Effect of different surfactants on the performance of UniSart® diagnostic membranes

2.1 hCG pregnancy test: an example for impacts of surfactants on signal intensity and background clearance

The hCG test is known to be robust and highly reproducible. Thus, it was selected to demonstrate the effects of surfactants on signal intensity and background clearance. The surfactants listed in Table 1 were applied in three categories of concentration: Low concentration: $\leq 0.1\%$ (w/w) or (v/w), moderate concentration: 0.2-0.5 % (w/w) or (v/w), and high concentration: $\geq 0.5\%$ (w/w) or (v/w). The half-strips assay was used to simplify the test process. This method is commonly used for the initial screening of buffer conditions (Table 2).

Construction of hCG test

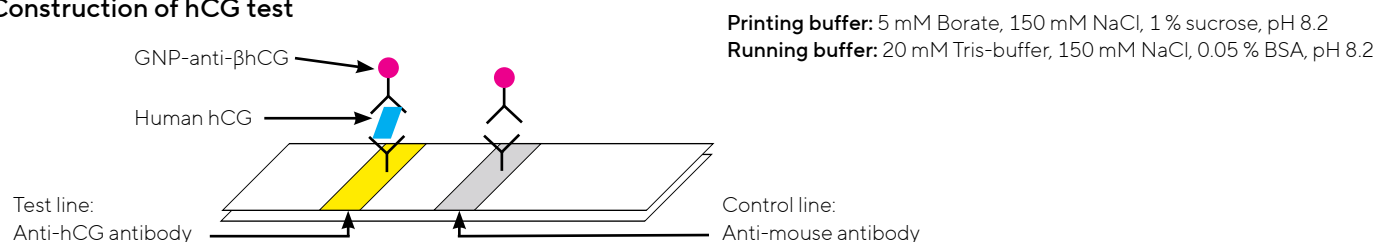


Table 2 . Summary of test results of hCG test with different surfactants

Criteria	Color	Description
Bad		No signal/strong background
Neutral		Low signal
Good		Strong signal/clear background

UniSart® Membrane	Surfactant Concentration	Test line and control line signal (naked eye detection)							
		Tween 20	Brij 35	Brij 58	Tween 60	Tergitol™ type NP-40	IGEPAL® CA-630	CHAPS	SDS
CN95 backed	Low								
	Moderate								
	High								
CN110 backed	Low								
	Moderate								
	High								
CN140 unbacked	Low								
	Moderate								
	High								
CN140 backed	Low								
	Moderate								
	High								
CN150 backed white	Low								
	Moderate								
	High								
CN150 backed clear	Low								
	Moderate								
	High								
CN180 backed	Low								
	Moderate								
	High								

Low concentration: $\leq 0.1\%$; Moderate concentration: 0.2-0.5 %; High concentration: $\geq 0.5\%$

2.2 The flu A test was chosen to demonstrate the effect of surfactants on non-specific binding

Figure 2. *Effect of surfactants on flu A test*

Printing buffer: 5 mM Borate, 150 mM, NaCl, 1 % sucrose, pH 8.2
Sample pad pre-treatment buffer: 100 mM Tris-buffer, 0.5 % BSA, 0.25 % surfactants, pH 8.0
Running buffer: 20 mM Tris-buffer, 150 mM NaCl, 0.05 % BSA, pH 8.2

CN95 backed membrane, 10000x

The analytes bind to nitrocellulose membrane non-specifically. Then, the nanoparticle conjugates can bind to these analytes³

The aggregates of nanoparticle conjugates are easily trapped in the pores of nitrocellulose membrane³

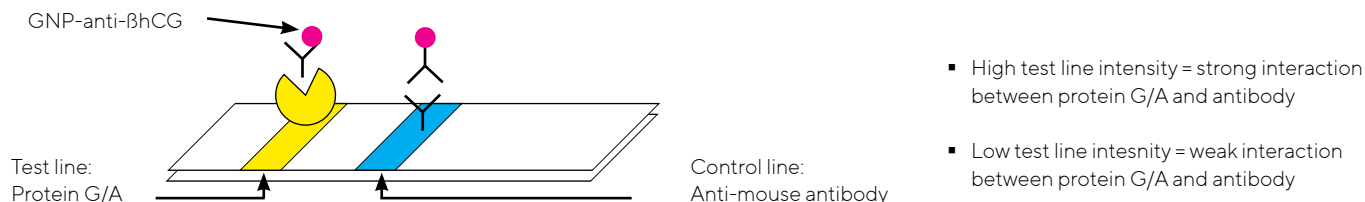
The nanoparticle conjugates bind non-specifically to nitrocellulose membrane through hydrophobic and electrostatic force³

T20: Tween 20; NP-40: Tergitol™ type NP-40

2.3 Protein G and protein A test: an example for surfactants' effect on the interaction of carrier proteins and antibodies

Protein G and protein A are commonly used in lateral flow assays to enhance the sensitivity. Both proteins interact with the Fc fragment of the antibody. It allows for better binding of the detection protein to the membrane surface. In such test, protein G/A-antibody complex is immobilized on the nitrocellulose membrane via protein G/A. Therefore, analysis of protein G/A on membrane immobilization and their interaction with antibodies are important for the optimization of the assay. Here, the effects of surfactants on protein G/A and their antibody interaction were evaluated (Table 3). The concentration of surfactants used 0.25 % (w/w) or (v/w) was adapted from the hCG test due to the good performance.

Construction of protein G and protein A test



Printing buffer for protein G: 100 mM Citrate, pH 5

Printing buffer for protein A: 5 mM Borate, 150 mM NaCl, 1 % sucrose, pH 8.2

Sample pad pre-treatment buffer: 100 mM Tris-buffer, 0.5 % BSA, 0.25 % surfactants, pH 8.0

Running buffer: 20 mM Tris-buffer, 150 mM NaCl, 0.05 % BSA, pH 8.2

Table 3. Summary of test results for protein G/A and antibody-interaction with different surfactants

Unisart® Membrane	Quantitative Signal Intensity (a.u.)											
	Tween 20		Brij 35		Brij 58		Tween 60		Tergitol™ type NP-40		IGEPAL® CA-630	
	Protein G	Protein A	Protein G	Protein A	Protein G	Protein A	Protein G	Protein A	Protein G	Protein A	Protein G	Protein A
CN95 backed	+++++	+++++	++	+++++	++++	+++	+++++	++++	+	+	+++	++
CN110 backed	+++++	+++++	+++	++++	++++	++	+++++	+++++	++	+	+	+++
CN140 unbacked	+++++	+++++	++++	+++++	+++	++++	+++++	+++	+	+	++	++
CN140 backed	+++++	+++++	+++	+++++	+++++	++++	++++	+++	+	++	++	+
CN150 backed white	+++++	+++	+++	+++	++++	++++	+++++	+++++	+	++	++	+
CN150 backed clear	+++++	++++	++++	++++	+++	+++	+++++	+++++	+	+	++	++
CN180 backed	+++++	+++++	+++++	+++++	+++	++	++++	+++	+	+	++	++++

+++++ : strongest interaction, + : weakest interaction

Conclusion

The effect of surfactants on test performance varied depending on the assay and the membrane, but also the type and concentration of surfactant used. Production costs must be considered if using a surfactant at high concentrations. It is therefore recommended to screen a variety of surfactants and concentrations to ensure the best performance.

Additional reading:

1. L. Zeng et al., Development of ic-ELISA and lateral-flow immunochromatographic strip for detection of vitamin B2 in an energy drink and vitamin tablets, *Food and Agricultural Immunology*, (2018) 29:1, 121-132.
2. Y. Zhao et al., Adsorbed Tween 80 is unique in its ability to improve the stability of gold nanoparticles in solutions of biomolecules, *Nanoscale* (2010) 2, 2114-2119.
3. L. Zhan et al., Development and optimization of thermal contrast amplification lateral flow immunoassays for ultrasensitive HIV p24 protein detection, *Microsystems & Nanoengineering* (2020) 6:543.
4. Global Lateral Flow Assays Market By type, By Technique, By Sample Type, By Application, By End User, By Regional Outlook, Industry Analysis Report and Forecast, 2021-2027.
5. G.A. Posthuma-Trumpie et al., Lateral flow (immune)assay: its strengths, weaknesses, opportunities, and threats. A literature survey, *Anal Bioanal Chem* (2009) 393:569-582
6. Y. Liu et al., Ultrasensitive and Highly Specific Lateral Flow Assays for Point-of-Care Diagnosis, *ACS Nano* (2021) 15 (3), 3593-3611
7. F. Di Nardo et al., Ten Years of Lateral Flow Immunoassay Technique Applications: Trends, Challenges and Future Perspectives. *Sensor* (2021) 2921, 21, 5185

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906



For additional information
visit www.sartorius.com