

# Micro-Flow Imaging View System Suite (V5.0)

## Title 21 Code of Federal Regulations (CFR) Part 11

### Compliance Checklist

This document supports the use of Micro-Flow Imaging® systems (MFI 5100, MFI 5200) in a QC and GMP environment. All functionality is integrated in MVSS (V5.0) for compliance with 21 CFR Part 11. The information presented assumes that the appropriate system configuration parameters have been set.

MVSS is a Closed System where access is controlled by the person(s) responsible for electronic records content. MVSS uses an ID/password combination and doesn't use any tokens or biometrics.

## Part 11, Subpart B — Electronic records

If you use the MVSS to create, modify, maintain, or transmit electronic records, you'll need procedures and controls to ensure authenticity, integrity, and, when appropriate, the confidentiality of the electronic record. Procedures and controls should also ensure that signers can't easily deny the validity of signed records.

### Section 11.10 — Controls for Closed Systems

SECTION 11.10(a)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is MFI View System Software (MVSS) validated?	✓			MVSS is validated and comes with a certificate of validation with software installation.
Does the validation documentation confirm Part 11 requirements have been met and are functioning correctly?	✓			MVSS software lets you be compliant with 21 CFR Part 11 guidelines, but complete compliance can only occur within a validated electronic records environment. Validation documentation is available upon request.
Are invalid records detected where applicable, like when there are invalid field entries, fields left blank that should contain data, or values outside of limits?	✓			Software utilities are in place to detect input errors/invalid data types, range checking for numeric entries, and marking of mandatory fields.

SECTION 11.10(b)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Can I view the entire contents of the records?	✓			
Can I print the entire contents of the records?	✓			
Can I generate all the records electronically in a format that can be put on a portable medium (e.g., diskette or CD) or transferred electronically?	✓			Methods, batches and reports as well as instrument and software data can be exported in an electronic format.

# Micro-Flow Imaging View System Suite (V5.0) Title 21 Code of Federal Regulations (CFR) Part 11 Compliance Checklist

SECTION 11.10(c)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are my records protected against intentional or accidental modification or deletion?	✓			The ability to modify data is a specifically assigned privilege. When you create or modify data, the action is logged to the audit trail and requires user confirmation before changes are saved. File management and keeping users from deleting data files is a customer responsibility.
Is my data archived off MVSS? If so, is the meta data (including the audit trail) archived as well? Can all the archived data be accurately retrieved after system upgrades?	✓			MVSS offers export functions for the data files, which include the metadata, as well as the audit trail for the system.

SECTION 11.10(d)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are there different levels of access based on user responsibilities (e.g., user, administrator)? Is this documented and controlled?	✓			
Are user access levels approved by management or the system owner before assignment to a user?	✓			
Is there is a controlled, documented process for granting access to a new user, for changing privileges for an existing user and for deleting user accounts?	✓			
Are there physical security and procedures to protect the server, database and system components from unauthorized access?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.

SECTION 11.10(e)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Can I automatically generate an electronic audit trail function all operator entries?	✓			
Is the audit trail completely outside the control and access of users (except for read-only access of the audit trail file)?	✓			
Is it impossible to disable the audit trail function?	✓			The audit trail is always active in MVSS.
Is MVSS's date and time protected from unauthorized change?			✓	Date and time are taken from the local date and time for the system's computer. The ability to change the system date and time is controlled through the computer operating system.
When my data is changed, are all previous values still electronically available?	✓			
Is the audit trail data protected from accidental or intentional modification or deletion?	✓			The audit trail can't be modified or deleted.
Are electronic audit trails maintained and retrievable for at least as long as its respective electronic records?	✓			
Are electronic audit trails readily available for inspections and audits?	✓			Audit trail entries are available for inspection by authorized users in the MVSS interface.

# Micro-Flow Imaging View System Suite (V5.0) Title 21 Code of Federal Regulations (CFR) Part 11 Compliance Checklist

## SECTION 11.10(e), *continued*

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Can I select portions of the audit trail for inspectors to view and print?	✓			The audit trail is available for review by authorized users and can be printed via MVSS.
Can I select portions of the audit trail to extract in a transportable electronic format for regulatory agencies?	✓			The audit trail can be downloaded as a PDF file by authorized users via the MVSS Print Audit trail tool.
If no audit trail is available, can MVSS detect that a record was altered since its last approval?	✓			MVSS software uses the SHA1 hash algorithm to generate a 160 bit hash code that is unique for all files. This hash code guarantees the file history is correct and no other edits were made.
Is the operator name, date, time, and indication of record (or file) creation, modification or deletion recorded in audit trail?	✓			
If the predicate regulation requires it, is the reason for a change included in the audit trail?		✓		

## SECTION 11.10(f)

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
If MFI requires sequenced steps, does it ensure that the actions are performed in the correct sequence?	✓			

## SECTION 11.10(g)

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Does MVSS only let authorized individuals use MFI?	✓			Anyone working with MVSS must have a user account. This account will define the capabilities you'll have on MFI. You won't have any access without an account.
Does MVSS verify that an individual has the authority to electronically sign a record before allowing them to do so?		✓		MVSS does not currently support electronic signatures.

## SECTION 11.10(h)

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Does MVSS require that data input or instructions only come from specific instrument/terminal? Does MVSS check for this?	✓			Instructions to an MFI instrument must come from proprietary Method and Batch files.

## SECTION 11.10(i)

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is there documentation to show that everyone who maintains or uses MVSS has the education, training and experience to perform their assigned tasks?	✓			Full documentation is available as part of an audit of ProteinSimple quality and training processes.
Is there documentation to show that everyone who maintains or uses MVSS has the education, training and experience to perform their assigned tasks?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.

Is there documentation to show that everyone who uses the system has the education, training and experience to perform their assigned tasks?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.
--	--	--	---	--

SECTION 11.10(j)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is there a written policy in place and enforced that holds users fully accountable and responsible for actions initiated under their electronic signatures?			✓	Each organization must their own develop controlled, documented procedures for compliance with this requirement.

SECTION 11.10(k)(1)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is the distribution of, access to, and use of MVSS operation and maintenance documentation controlled?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.
Is access to “sensitive” MVSS documentation restricted e.g., network security documentation, system access documentation?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.

SECTION 11.10(k)(2)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is there a Change Control (or equivalent) SOP governing revisions to system documentation?			✓	Each organization must their own develop controlled, documented procedures for compliance with this requirement.

## Section 11.30 — Controls for open systems

SECTION 11.30				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are there controls to ensure record authenticity, integrity, and confidentiality?			✓	MVSS is a closed system.
Is data encrypted?			✓	MVSS is a closed system.
Are digital signatures used?			✓	Digital signatures are not currently available in MVSS.

## Section 11.50 — Signature manifestations

SECTION 11.50(a)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Do all electronically signed records contain the full printed name of the signer?			✓	Digital signatures are not currently available in MVSS.
Do all electronically signed records contain the date and time of signing?			✓	Digital signatures are not currently available in MVSS.
Do all electronically signed records contain the meaning of the signature (e.g., review, approval)?			✓	Digital signatures are not currently available in MVSS.

Are the date and time stamps applied automatically and not keyed in by the user?			✓	
Are the date and time stamps presented in a consistent way that makes it simple to reconstruct the sequence of events?			✓	

SECTION 11.50(b)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is the above information subject to the same controls as electronic records? (audit trail, access control, etc.)	✓		✓	Digital signatures are not currently available in MVSS.
Are changes to signatures included in the audit trail?			✓	
Do the printed name, date, time, and signature meaning appear in printed reports of the electronic record?			✓	

## Section 11.70 — Signature/record linking

SECTION 11.70				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
If handwritten signatures are executed to electronic records, are the handwritten signatures linked to the electronic records?			✓	Handwritten signatures are not executed to electronic records. Handwritten signatures may be executed to a printed report, and such a report may include information identifying (and providing a link to) the original electronic record.
If the electronic record is changed, is the signer prompted to re-sign (via either manual procedures (SOP) or technical means)?			✓	All changes to an electronic record are audit trailed. The audit trail includes information on the user making the change, the date and time of the change, what was changed and the reason for the change, but there is no support for electronic signatures.
Are the E-signatures linked (via technology, not procedures) to their corresponding electronic records to ensure that the signature can't be excised, copied, or otherwise transferred to falsify an electronic record by ordinary means?			✓	

## Part 11, subpart C— Electronic signatures

### Section 11.100 — General requirements

SECTION 11.100(a)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is each E-signature unique to one individual?			✓	
Are E-signatures ever reused by, or reassigned to, anyone other than the original owner?			✓	

SECTION 11.100(b)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is the individual identified adequately verified prior to issuance of an electronic signature?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.
Is there a procedure for reissuing forgotten passwords that verifies the requestor's identity?			✓	Each organization must develop their own controlled, documented procedures for compliance with this requirement.

SECTION 11.100(c)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Has certification of the intent to use electronic signatures been submitted to the agency in paper form with a traditional handwritten signature?			✓	
Can additional certification or testimony be supplied to show that an electronic signature is the legally binding equivalent of the signer's handwritten signature?			✓	

## Section 11.200 — Electronic signature components and controls

SECTION 11.200(a)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Does the signature consist of at least two distinct identification components?			✓	
If continuous signing sessions are used, are two (or more) E-signature components required for the initial signing?			✓	
If only one E-signature component is required for subsequent signings: <ul style="list-style-type: none"> <li>• Is the private component, known to and only useable by its owner, used for each subsequent signing?</li> <li>• Is the user required to stay in close proximity to the workstation for the entire session?</li> <li>• Is there an automatic logoff, or password-protected screen saver that launches after a short period of inactivity (with the password known only by one user)?</li> </ul>			✓	
If I leave the workstation, do procedures and/or automatic controls ensure that it's treated as a non-continuous session?		✓		
Are two (or more) E-signature components required for each signing during a non- continuous signing session?			✓	
Are non-biometric signatures only used by their genuine owners (e.g., by procedures or training reinforcing that non-biometric E-signatures are not "loaned" to co-workers or supervisors for overrides)?			✓	

**SECTION 11.200(a)**, *continued*

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are non-biometric signatures administered and executed so that unauthorized use requires the collaboration of two or more individuals?			✓	

**SECTION 11.200(b)**

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are biometric E-signatures designed to ensure that they can be used only by their genuine owners?			✓	MVSS doesn't use biometric E-signatures.

**Section 11.300 — Controls for identification codes/passwords**

**SECTION 11.300(a)**

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are controls in place to maintain the uniqueness of each combined identification code and password, such that no two individuals can have the same combination of identification code and password?	✓			MVSS ensures that there must be unique combination of user names and passwords used on the system.
Are controls (procedural or technical) in place to prevent the re-use of identification codes?	✓			MVSS policies can be used to ensure that passwords may not be reused for individual user accounts, and that user names can't be reused for multiple users.

**SECTION 11.300(b)**

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is the issuance of identification codes and passwords periodically checked, recalled, or revised (e.g., to cover such events as password aging)?	✓			MVSS policies can be used to set password aging based on corporate policies.
Do passwords periodically expire and need to be revised?	✓			MVSS policies can be used to set password aging based on corporate policies.
Is there a procedure for recalling identification codes and passwords if a person leaves or is transferred?	✓			MVSS allows a user account to be removed from active use. Each organization must develop controlled, documented procedures to ensure proper notification of user status changes.

**SECTION 11.300(c)**

REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Is an SOP in place directing action to be taken to electronically deauthorize lost, stolen, missing, or otherwise potentially compromised tokens, cards, and other devices used to carry or generate E-signature components?			✓	MVSS doesn't use tokens, cards or other devices to carry E-signature components.
Does this SOP contain procedures for managing and controlling temporary or permanent token/ card replacements?			✓	MVSS doesn't use tokens, cards or other devices to carry E-signature components.

SECTION 11.300(d)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are any attempts to unauthorized use detected and reported immediately to the system "security unit" (e.g., a system administrator is notified automatically by console message or paper) and, as appropriate, to organizational management?	✓			System records invalid Login attempts to Audit Trail. Five (5) login attempts are allowed upon MVSS startup, after which the program closes. Automatic notification is not handled by MVSS.

SECTION 11.300(e)				
REGULATION	YES	NO	N/A	IMPLEMENTATION AND APPLICATION
Are there procedures covering the initial and periodic testing of devices, such as tokens or cards that bear or generate identification code or password information?			✓	MVSS doesn't use tokens, cards or other devices to carry E-signature components.
Does the testing include checks for proper functioning, performance degradation, and possible unauthorized alteration?			✓	MVSS doesn't use tokens, cards or other devices to carry E-signature components.

## Reference

1. FDA Regulation 21 CFR Part 11 — Electronic Records; Electronic Signatures (FDA, 1997)