chemgineering

ERES Certificate

Software: SoftMax Pro GxP

Version:



Lagerhausstr. 45 | 5071 Wals | Austria www.MolecularDevices.com

Assessment Date:

11-Apr-2019

Auditor:

Sieghard Wagner, mech. Engineer (grad.), Chemgineering Business Design GmbH

Assessment objective: Compliance check against the Electronic Records and Electronic Signatures (ERES) requirements of FDA 21 CFR Part 11 and EudraLex, Volume 4, Annex 11 for

SoftMax Pro GxP 7.1

Object description:

Software to process preconfigured protocols and custom assay workflows for microplate data

acquisition and analysis.

Approach:

The functions and properties of the above software system are audited against the ERES requirements of 21 CFR Part 11 and EU GMP Annex/11 with its current interpretations. The assessment is guided by a requirements checklist, which is based on the respective 21 CFR Part 11 and Annex 11 paragraphs.

Assessment results:

Operator responsibility

ERES compliance requires an appropriate operational environment.

It is the operations responsibility to provide a compliant environment regarding:

- Technical environment, data management
- Training
- Administration
- Standard Operating Procedures (SOP).

Summary

The software is compliant with the following ERES requirements:

21 CFR Part 11:

11.10 (b), (d), (f), (g), (h); 11.50, 11.70; 11.300 (a)

EU GMP Annex 11:

Paragraphs: 4.8, 6, 8, 9, 12, 14b, 14c

The software is compliant with the following ERES requirements with support of the operator:

21 CFR Part 11:

11.10 (a), (c), (e), (i), (j), (k); 11.100 (a), (b);

11.200 (a); 11.300 (b), (c), (d)

EU GMP Annex 11:

Paragraphs: Principles, 1 to 4.7, 7, 9, 10, 14a, 17

The software supports electronic signature.

Sieghard Wagner